

Company Name	Reporting Group	Country	Address
McCrometer, Inc.	McCrometer	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
Aguasin SpA	ChemTreat	Chile	Carey y Cía. Ltda. Isidora Goyenechea 2800, Piso 43 Las Condes / Santiago / Chile Chile
ChemTreat International, Inc.	ChemTreat	United States	CT Corporation System 4701 Cox Road, Suite 301 Glen Allen VA 23060 United States
ChemTreat Mexico, S. de R.L. de C.V.	ChemTreat	Mexico	Via Gustavo Baz 2160 Edificio 2 PB Col. Parque Industrial la Loma Tlalnepantla de Baz Edo. Mex CP 54060 Mexico
ChemTreat Puerto Rico, Inc.	ChemTreat	United States	CT Corporation System 4701 Cox Road, Suite 301 Glen Allen VA 23060 United States
ChemTreat Trinidad Limited	ChemTreat	Trinidad and Tobago	17-20 Queen's Park West Port of Spain Trinidad and Tobago
ChemTreat Venezuela, C.A.	ChemTreat	Venezuela, Bolivarian Republic of	Avenida FCO Miranda C.C. Multicentro Empresarial Del Este EDF Miranda Nivel Torre A, Piso 4, Oficina A-41 Caracas Capital 1060 Venezuela
ChemTreat, Inc.	ChemTreat	United States	CT Corporation System 4701 Cox Road, Suite 301 Glen Allen VA 23060 United States
Lipeq S.A.	ChemTreat	Ecuador	Ferrere Abogados Mirage Building, Avenida 12 de Octubre N24-68 , esq. Lincoln Edf. Mirage Piso 16 Ecuador
Lipesa Colombia S.A.S.	ChemTreat	Colombia	Carretera central Bogota-Tuja frente a Bavaria KM 21 Tocancipa Cundinamarca Colombia
Lipesa del Perú S.A.C.	ChemTreat	Peru	No Registered Agent

Lipesa do Brasil Especialidades Químicas Ltda.	ChemTreat	Brazil	Caçada dos Mancás, No. 26, Qd. 18, Lt. 6, Centro Comercial de Alphaville, city of Barueri São Paulo, 06.453-036 Brazil
U.S. Peroxide, LLC	ChemTreat	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
USP Technologies Canada ULC	ChemTreat	Canada	250 HOWE STREET SUITE 1400 VANCOUVER BC V6C 3S7 Canada
AB Sciex (Thailand) Limited	Hach	Thailand	Chankasem Subdistrict Chatuchak District Bangkok 10900 Thailand
Analytical Developments Limited	Hach	Ireland	Raffeen House, Ringaskiddy, Co. Cork Ireland
AppliTek NV	Hach	Belgium	Venecoweg 19, B-9810 Nazareth Belgium
Aquatic Informatics Australia Pty Ltd	Hach	Australia	HWL EBSWORTH LAWYERS 'AUSTRALIA SQUARE' LEVEL 14 264-278 GEORGE STREET SYDNEY NSW 2000 Australia
Aquatic Informatics Inc.	Hach	United States	C T Corporation System 711 Capitol Way S, Suite 204 Olympia WA 98501 United States
Aquatic Informatics ULC	Hach	Canada	2400 - 1111 West Georgia Street Vancouver BC V6E 4M3 Canada
Aquatic Informatics USA Inc.	Hach	United States	C T Corporation System 711 Capitol Way S, Suite 204 Olympia WA 98501 United States
ATTC Training Center Sdn Bhd	Hach	Malaysia	19-3, Jalan Tasik Selatan 3, Bandar Tasik Selatan 57000 Kuala Lumpur Malaysia
Biotech Analytical Systems Limited	Hach	Ireland	Unit 34 GB Business Park Little Island Co. Cork T45 H681 Ireland
Danaher (Shanghai) Instrumentation and Industrial Technologies R&D Company Limited	Hach	China	Room 101/201/301, Building 10, 518 518 North Fu Quan Road, Shanghai China

DH Water Quality LLC	Hach	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
Dr. Lange Nederland B.V.	Hach	Netherlands	Laan van Westroijen 2a 4003 AZ Tiel Netherlands
Hach Colombia S.A.S.	Hach	Colombia	Calle 60A # 5-77 Bogotá Cundinamarca 110231 Colombia
Hach Company	Hach	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
Hach Korea Yuhanhoesa (Hach Korea, LLC)	Hach	Korea, Republic of	2th FL, 5, Yeongdong-daero 106-gil, Gangnam-gu Seoul Korea, Republic of
Hach Lange AB	Hach	Sweden	Roentgenvaegen 5-7 171 54 Solna Sweden
Hach Lange ApS	Hach	Denmark	Akandevvej 21 DK-2700 Bronshoj Denmark
Hach Lange Controle e Analise de Aguas Ltda.	Hach	Portugal	Avenida do Forte nº 3 Edifício Suécia III – Lote 035 – R/C 2794-044 Carnaxide-Lisboa Portugal
Hach Lange d.o.o. [Croatia]	Hach	Croatia	Ivana Severa 20 Varazdin 42000 Croatia
Hach Lange d.o.o. [Slovenia]	Hach	Slovenia	Fajfarjeva cesta 15 Domzale 1230 Slovenia
Hach Lange EOOD (Bulgaria)	Hach	Bulgaria	45 Kr. Sarafov Str Sofia 1164 Bulgaria
Hach Lange Finance GmbH	Hach	Germany	Königsweg 10 Zehlendorf 14163 Berlin Germany
Hach Lange France S.A.S.	Hach	France	8 MAIL BARTHELEMY THIMONNIER ROND POINT DU MANDINET LOGNES 77437 France

Hach Lange GmbH	Hach	Germany	Königsweg 10 Zehlendorf Berlin 14163 Berlin Germany
Hach Lange GmbH [Austria]	Hach	Austria	Hütteldorferstraße 299/Top 6 1140 Wien Austria
Hach Lange Holding GmbH	Hach	Germany	Königsweg 10 Zehlendorf 14163 Berlin Germany
Hach Lange Kereskedelmi Kft.	Hach	Hungary	Vöröskereszt utca 8-10. Budapest 1222 Hungary
Hach Lange LTD (Ireland)	Hach	Ireland	Unit 1, Chestnut Road Western Industrial Estate Dublin 12 Ireland
Hach Lange LTD (UK)	Hach	United Kingdom	19 Jessops Riverside 800 Brightside Lane Sheffield S9 2RX United Kingdom
HACH LANGE Monoprosopi EPE	Hach	Greece	3rd Irakleitou 15235 Chalandri Greece Greece
Hach Lange N.V.	Hach	Belgium	Venecoweg 19, B-9810 Nazareth Belgium
Hach Lange NZ	Hach	New Zealand	c/o Russell McVeagh Vero Centre 48 Shortland Street Auckland NZ 1140 New Zealand
Hach Lange OOO, Russia	Hach	Russian Federation	21 Stanislavskogo, Bldg 3 Ibaraki 109004 Moscow Russian Federation
Hach Lange S.r.l. (Italy)	Hach	Italy	Via Rossini 1/a 20045 Lainate (Mi) Italy
Hach Lange s.r.l. [Romania]	Hach	Romania	str. Caminului, nr. 3, etaj 1, apartament 1 sector 2 Bucuresti Romania

Hach Lange s.r.o. (Slovakia)	Hach	Slovakia	Roľnícka 21 Bratislava - Vajnory 831 07 Slovakia
HACH LANGE s.r.o. [Czech]	Hach	Czech Republic	Praha 4, Zastrčená 1278/8 141 00 Czech Republic
Hach Lange Sàrl	Hach	Switzerland	6, route de Compois CH 1222 Vesenaz Switzerland
Hach Lange Sensors SAS	Hach	France	Chez HL France: 8 Bvd Barthélémy Thimonnier 77185 Lognes France
Hach Lange Spain SLU	Hach	Spain	Plaça d'Europa 41-43 Torre Realia 4ª Planta 08908 L'HOSPITALET DE LLOBREGAT (Barcelona) Spain
Hach Lange Spółka z.o.o. (Poland)	Hach	Poland	ul. Krakowska 119 50-428 Wrocław Poland
Hach Lange, Su Analiz Sistemleri Ltd. Sti	Hach	Turkey	Mustafa kemal Mah. 2141. Cadde No: 8 Çankaya/ ANKARA 06530 Turkey
Hach Malaysia Sdn Bhd	Hach	Malaysia	19-3, JALAN TASIK SELATAN 3, BANDAR TASIK SELATAN, 57000 KUALA LUMPUR WILAYAH PERSEKUTUAN Malaysia
Hach Pacific Pty Ltd	Hach	Australia	Level 1, 96 Ricketts Road Mount Waverly VIC 3149 Australia
Hach Sales & Service Canada LP	Hach	Canada	PwC Law LLP 18 York Street, Suite 2600 Toronto ON M5J 0B2 Canada
Hach South Africa Pty. Ltd	Hach	South Africa	PO Box 144 Kya Sand Randburg 2163 South Africa
HACH ULTRA JAPAN KK	Hach	Japan	TD Bldg.5F, 1-29-9 Tokyo Takadanobaba, Shinjyuku-ku 169-0075 Japan
Hach Water Analytical Instruments (Shanghai) Co. Ltd.	Hach	China	3F, No. 6, Lane 280, Linhong Road, Shanghai 200335, PRC China

Hach Water Quality Technology (Hangzhou) Co., Ltd.	Hach	China	RM 313 / 315, Block B, No. 90 Wen San Road, Technology Innovation Base Hangzhou PRC China
Hangzhou Hach Environmental Science and Technology Co., Ltd	Hach	China	2F, Building #8, No. 1418 – 41, Mo Ganshan Road, Hangzhou Zhejiang China
Hexis Científica Ltda.	Hach	Brazil	Av. Antonieta Piva Barranqueiros, 385 Distrito Industrial Jundiá São Paulo 13212-000 Brazil
Reactivos y Equipo S.A. de C.V.	Hach	Mexico	José Silvestre Arramberi Poniente 1442 Monterrey Centro Monterrey Monterrey Nuevo León 64000 Mexico
Sea-Bird Electronics, Inc.	Hach	United States	C T Corporation System 711 Capitol Way S, Suite 204 Olympia WA 98501 United States
Sea-Bird GmbH	Hach	Germany	Ludwigstraße 16 87437 Kempten Germany
Sedaru, Inc.	Hach	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
Shanghai Shilu Instrument Co. Ltd.	Hach	China	No. 139 Lane 2638 Hong Mei South Road Min Hang District Shanghai City 201108 China
Termodinamica Limitada	Hach	Chile	Carey y Cía. Ltda. Isidora Goyenechea 2800, Piso 43 Las Condes / Santiago / Chile Chile
Water Quality GmbH	Hach	Germany	Koenigsweg 10 D-14163 Berlin Germany
WaterTrax USA Inc.	Hach	United States	C T Corporation System 711 Capitol Way S, Suite 204 Olympia WA 98501 United States
WESTERN ENVIRONMENTAL TECHNOLOGY LABORATORIES, INC.	Hach	United States	CT Corporation System 780 Commercial Street SE, Ste. 100 Salem OR 97301 United States

X-Ray Optical Systems, Inc.	Hach	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
Alltec Angewandte Laserlicht Technologie GmbH	Videojet	Germany	An der Trave 27/31 23923 Selmsdorf Germany
CD Codificação LTDA	Videojet	Brazil	Av Raul Mourao Guimaraes, 15 Brazil
DH Netherlands BV	Videojet	Netherlands	Strijkviertel 53 3454 PK De Meern Netherlands
DHR MEA GENERAL TRADING LLC	Videojet	United Arab Emirates	EMIRATE OF DUBAI / WAREHOUSE NO 1 /B-416, PLOT NUMBER 416-5483, NADD AL HAMAR, DUBAI United Arab Emirates
FOBA Italia S.r.l.	Videojet	Italy	Via XI Febbraio 2 21052 Busto Arsizio (VA) Italy
J.S.C. Videojet Technologies	Videojet	Russian Federation	108811, Moscow, Kievskoye shosse 22-y km. (s. Moscovskiy) household 4, building 5, floor 7, block E, office 701E 108811 Moscow Russian Federation
Laetus France S.a.r.l.	Videojet	France	9 avenue du Lac, Bois Briard 91080 COURCOURONNES France
Laetus GmbH	Videojet	Germany	Sandwiesenstrasse 27, 64665 Alsbach-Hähnlein Germany
Linx Asia Limited	Videojet	Hong Kong	Room 14-15, 2nd Floor, Block A2, Yau Tong Industrial City, 17 Ko Fai Road, Yau Tong Hong Kong
Marsh Interex Ltd.	Videojet	Jamaica	c/o 1500 North Mittel Blvd. Wooddale IL 60161-1073 Jamaica
Marsh Label Technologies LLC	Videojet	United States	C T Corporation System 120 South Central Avenue Clayton MO 63105 United States
Videojet X-Rite K.K.	Videojet	Japan	Telecom Center West Bldg.6F,2-5-10 Aomi, Koto-ku Tokyo 135-0064 Japan

Videojet Argentina S.R.L.	Videojet	Argentina	Marval Law Firm Av. Leandro N. Alem 928 Buenos Aires C1001AAR Argentina
Videojet Canada Limited Partnership	Videojet	Canada	PwC - Vancouver 250 Howe Street Suite 1400 Vancouver BC V6C3S7 Canada
Videojet Chile Codificadora Limited	Videojet	Chile	EXEQUIEL FERNANDEZ N° 2831 MACUL SANTIAGO Chile
Videojet Do Brasil Comércio de Equipamentos Para Codificação Industrial Ltda.	Videojet	Brazil	Rua São Paulo 261 and 291 Alphaville Barueri São Paulo 06465-130 Brazil
Videojet Italia Srl	Videojet	Italy	Vicolo San Michele n. 15 21100 Varese Italy
Videojet Korea Co. Ltd.	Videojet	Korea, Republic of	Videojet Korea 13F Pantech R&D tower 179 Seongam-ro Mapo-gu Seoul Korea 133-834 Korea, Republic of
Videojet Technologies (I) Pvt. Ltd.	Videojet	India	S-220 A, 2nd Floor, Eastern Business District Neptune Mall, L B S Marg Bhandup West Mumbai Maharashtra 400078 India
Videojet Technologies (Nottingham) Limited	Videojet	United Kingdom	4 & 5 Ermine Centre Lancaster Way Huntingdon England PE29 6XX United Kingdom
Videojet Technologies (S) PTE. LTD.	Videojet	Singapore	1 Kaki Bukit View #04-15/17 Techview 415941 Singapore
Videojet Technologies (Shanghai) Co., Ltd.	Videojet	China	Part J, Building 50, No. 389 YingLun Road Waigaoqiao Free Shanghai City 200131 China

Videojet Technologies B.V.	Videojet	Netherlands	Gildenstraat 33 4143 HV Leerdam Netherlands
Videojet Technologies Czechia s.r.o.	Videojet	Czech Republic	Na Hlavní 166 182 00 Praha 8, Březiněves Czech Republic
Videojet Technologies Europe B.V.	Videojet	Netherlands	Strikviertel 39 PJ DeMeern 3454 Netherlands
Videojet Technologies GmbH	Videojet	Germany	An der Meil 2 65555 Limburg Hessen Germany
Videojet Technologies Inc.	Videojet	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
Videojet Technologies Limited	Videojet	United Kingdom	4 & 5 Ermine Business Centre Lancaster Way Huntingdon Cambridgeshire PE29 6XX United Kingdom
Videojet Technologies Mexico S. de R.L. de C.V.	Videojet	Mexico	Av. Coyoacán # 1213 Col. Del Valle Mexico D.F 03100 Mexico
Videojet Technologies Mexico Services S. de R.L. de C.V.	Videojet	Mexico	Av. Coyoacán # 1213 Col. Del Valle Mexico D.F 03100 Mexico
Videojet Technologies S.A.S.	Videojet	France	SILIC 703, 16 Avenue du Québec, Villebon sur Yvette 91961 Courtaboeuf Cedex France
Videojet Technologies S.L.	Videojet	Spain	C/Valgrande, 8 Nave B1A 28108 Alcobendas 28108 Madrid Spain
Videojet Technologies Slovakia s.r.o.	Videojet	Slovakia	Einsteinova 25 851 01 Bratislava - mestská časť Petržalka Slovakia
Videojet Technologies SP z.o.o.	Videojet	Poland	ul. Kolejowa 5/7 01-217 Warszawa Poland

Videojet Technologies Suisse GmbH	Videojet	Switzerland	Gummertlistrasse 7 CH-4702 Oensingen Switzerland
Videojet Technologies Urun Kodlama ve Etiketleme Danismanlik ve Ticaret Limited Sirketi	Videojet	Turkey	Cubukcuoglu Is Merkezi, Ruya sokak no 11, Kucukbakkalkoy, Özel İşyeri:4 Atasehir / Istanbul 34750 Turkey
VJ HoldCo GmbH & Co. KG	Videojet	Germany	An der Meil 2 65555 Limburg Germany
VJH GmbH	Videojet	Germany	An der Meil 2 65555 Limburg Germany
VTI Sweden AB	Videojet	Sweden	Johannefredsgaten 4 431 53 Möndal Sweden
Willett America, Inc.	Videojet	United States	C T Corporation System 350 North St. Paul Street, Suite 2900 Dallas TX 75201 United States
Willett GmbH	Videojet	Germany	An der Meil 2 65555 Limburg Germany
Willett Holdings B.V.	Videojet	Netherlands	Strijkviertel 53 3454 PK De Meern Netherlands
Willett International Limited	Videojet	United Kingdom	4 & 5 Ermine Business Centre Lancaster Way Huntingdon Cambridgeshire PE29 6XX United Kingdom
Willett S.A.	Videojet	Uruguay	Ruta 8, KM 17, 500 Local 111 Zona America Montevideo 12300 Uruguay
Zhuhai S.E.Z. Videojet Electronics Ltd.	Videojet	China	No.8, Xinghan Road Sanzao Science and Technology Park Zhuhai Guangdong 519040 China
Beckman Australia ApS	Radiometer	Denmark	c/o Radiometer Medical Aps Akandevvej 21 2700 Bronshoj Denmark

Beckman Coulter France ApS	Radiometer	Denmark	c/o Radiometer Medical ApS Åkandevej 21 2700 Brønshøj Denmark
Beckman Finance ApS	Radiometer	Denmark	c/o Radiometer Medical ApS Åkandevej 21 2700 Bronhoj Denmark
Danaher Finance ApS	Radiometer	Denmark	Åkandevej 21 DK 2700 Bronshoj Denmark
Danaher Medical ApS	Radiometer	Denmark	c/o Radiometer Medical ApS Åkandevej 21 2700 Brønshøj Denmark
Danaher Tax Administration ApS	Radiometer	Denmark	Åkandevej 21 DK-2700 Bronshoj Denmark
DH Denmark Holding ApS	Radiometer	Denmark	Akandevej 21 DK-2700 Bronshoj Denmark
DHCDAN Holding ApS	Radiometer	Denmark	Akandevej 21 DK-2700 Bronshoj Denmark
KVF Finance ApS	Radiometer	Denmark	c/o Radiometer Medical ApS Åkandevej 21 DK-2700 Bronshoj Denmark
Radiometer (Suzhou) Medical Technology Co., Ltd.	Radiometer	China	Room 207, No. 11 Modern Industrial Lane, No. 333 Xingpu Road, Suzhou Industrial Park Suzhou Free Trade Zone China
Radiometer America Inc.	Radiometer	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
Radiometer Arabia Co. Ltd.	Radiometer	Saudi Arabia	RAD Saudia Maliha Street, Building No 7546 Secondary No. 2739 District Qurtubah Dist. Riyadh 13245 Kingdom of Saudi Arabia
Radiometer Austria GmbH	Radiometer	Austria	c/o DLA Piper Weiss- Tessbach Rechtsanwaelte GmbH Schottenring 14 1010 Wien Austria

Radiometer Basel AG	Radiometer	Switzerland	Austrasse 25 4051 Basel Switzerland
Radiometer Benelux B.V.	Radiometer	Netherlands	Chroomstraat 8 2718 RR Zoetermeer Netherlands
Radiometer Canada Limited Partnership	Radiometer	Canada	PwC - Vancouver 250 Howe Street Suite 1400 Vancouver BC V6C3S7 Canada
Radiometer d.o.o.	Radiometer	Croatia	Avenija V. Holjevca 40 , 10000 ZAGREB Croatia
Radiometer GmbH	Radiometer	Germany	Europark Fichtenhain A4 47807 Krefeld Germany
Radiometer HK Limited	Radiometer	Hong Kong	22/F. AXA Tower, Landmark East 100 How Ming Street Kwun Tong Kowloon Hong Kong
Radiometer Iberica S.L.	Radiometer	Spain	Avenida Valgrande numero 8 Planta secunda Alcobendas Madrid Spain Spain
Radiometer Ireland Limited	Radiometer	Ireland	Lismeehan, O' Callaghan Mills, Co. Clare Ireland
Radiometer K.K.	Radiometer	Japan	4-7-35, Kita-Shinagawa Shinagawa-ku Tokyo 140-0001 Japan
Radiometer Kazakhstan LLP	Radiometer	Kazakhstan	98 Zhandosov street, 6th floor Auezovsky District Almaty 050042 Kazakhstan
Radiometer Korea Co., Ltd.	Radiometer	Korea, Republic of	(Cheongdam-dong) 741, Yeongdong-daero, Gangnam- gu, Seoul Korea, Republic of
Radiometer Limited	Radiometer	United Kingdom	Manor Court, Manor Royal Crawley West Sussex RH10 9FY United Kingdom
Radiometer Magyarország Korlátolt Felelősségű Társaság	Radiometer	Hungary	Papírgyár u. 58-5 Budapest 1038 Hungary

Radiometer Malaysia Sdn. Bhd.	Radiometer	Malaysia	Suite 1005 10th Floor Wisma Hamzah-Kwong Hing, No.1 Leboh Ampang 50100 Kuala Lumpur Wilayah Persekutuan Malaysia
Radiometer Medical ApS	Radiometer	Denmark	Akandevvej 21 Bronshoj DK-2700 Denmark
Radiometer Medical Equipment (Shanghai) Co. Ltd.	Radiometer	China	Room 311&312 No.4 Building No. 222 Meiyue Road Shanghai Pilot Free Trade Zone China
Radiometer Mexico S.A. de R.L. de C.V.	Radiometer	Mexico	Av. Popocatépetl 396, Gral Anaya Mexico City 03340 Mexico
Radiometer OOO	Radiometer	Russian Federation	125167, Moscow, 37 Leningradskiy Prospekt, bld. 9. Russian Federation
Radiometer Pacific	Radiometer	New Zealand	3/33 Spartan Road Takanini Takanini Auckland 2105 New Zealand
Radiometer Pacific Pty Limited	Radiometer	Australia	1st floor, 96 Ricketts Rd Mount Waverly VIC 3149 Australia
Radiometer RSCH GmbH	Radiometer	Switzerland	Zürcherstrasse 68 8800 Thalwil Switzerland
Radiometer S.A.S.	Radiometer	France	Z.A. la Fontaine du Vaisseau 8, rue Edmond Michelet Neuill F-93360 France
Radiometer S.E.A. PTE. LTD.	Radiometer	Singapore	1 Science Park Road, #05 - 09 East Wing The Capricorn, Singapore Science Park II Singapore 117528 Singapore
RADIOMETER s.r.o.	Radiometer	Czech Republic	Praha 6 Krenova 3 PSC 162 00 Czech Republic
Radiometer Solutions Sp. z o.o.	Radiometer	Poland	Stalowa 6 73-110 Stargard Poland

Radiometer South Africa Pty Ltd	Radiometer	South Africa	MSI Office Business 68 Rigger Road, Spartan Kempton Park 1619 South Africa
Radiometer Spolka z o.o.	Radiometer	Poland	RAD Spolka Aleje Jerozolimskie 181A 02-222 Warsaw
Radiometer Tibbi Malzemeler Sanayi Ve Ticaret Limited Sirketi	Radiometer	Turkey	FSM Mah. Poligon Cad. No:8C Buyaka Kule-3 K:8 D:54, K:-5/A Tepeüstü, Umraniye Istanbul 34771 Turkey
Radiometer Turku Oy	Radiometer	Finland	Biolinja 12 P.O. Box 343 Turku FI-20750 Finland
Sarbel Acquisition ApS	Radiometer	Denmark	c/o Radiometer Medical Aps Akandevej 21 2700 Bronshoj Denmark
SenDx Medical, Inc.	Radiometer	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
Aquafine Corporation	Trojan	United States	C T Corporation System 818 West 7th Street Los Angeles CA 90017 United States
Salsnes Filter AS	Trojan	Norway	Verftsgata 32 Namsos 7800 Norway
Trojan Technologies Corp.	Trojan	United States	1209 Orange Street Corporation Trust Center Wilmington DE 19801 United States
Trojan Technologies Deutschland GmbH	Trojan	Germany	Aschaffener Str 72 63825 Schöllkrippen Germany
Trojan Technologies Espana S.L.	Trojan	Spain	Calle Orense, planta 1 despacho 32 28002 Madrid Spain
Trojan Technologies Group ULC	Trojan	Canada	250 HOWE STREET SUITE 1400 VANCOUVER BC V6C 3S7 Canada
TROJAN TECHNOLOGIES, INC.	Trojan	United States	C T Corporation System 289 S. Culver St. Lawrenceville GA 30046

Trojan UV Technologies UK Limited	Trojan	United Kingdom	19 Jessops Riverside 800 Brightside Lane Sheffield S9 2RX United Kingdom
Viqua Holdings	Trojan	Canada	PwC Law LLP 18 York Street, Suite 2600 Toronto ON M5J 0B2 Canada
Linx Alltec (Shanghai) Trading Limited	Linx	China	Room 1-118, Floor 1, Block 4, Building 1, No. 2001 North Yanggao Road, Shanghai Free Trade Area Shanghai China
Linx Printing Technologies Limited	Linx	United Kingdom	Linx House 8 Stocks Bridge Way Compass Point Business Park St Ives Cambridgeshire PE27 5JL United Kingdom
Linx S.A.S.	Linx	France	2 rue du Bois Sauvage Cedex 91000 Evry France
Artemis Medical, Inc.	LBS	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
Digital Slidebox Limited	LBS	Ireland	BUILDING 1 SWIFT SQUARE NORTH WOOD SANTRY Dublin 9 Ireland
Kreatech Biotechnology B.V.	LBS	Netherlands	Vlierweg 20, 1032LG Amsterdam Netherlands
Kreatech, Inc.	LBS	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
Leica Biosystems (Dublin) Limited	LBS	Ireland	Lismeehan, O' Callaghan Mills, Co. Clare Ireland
Leica Biosystems Deutschland GmbH	LBS	Germany	Heidelberger Strase 17-19 69226 Nussloch Germany
Leica Biosystems Imaging, Inc.	LBS	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States

Leica Biosystems Inc.	LBS	United States	1209 Orange Street Corporation Trust Center Wilmington DE 19801 United States
Leica Biosystems Melbourne Pty Ltd	LBS	Australia	495 Blackburn Road Mount Waverly VIC 3149 Australia
Leica Biosystems Newcastle Limited	LBS	United Kingdom	19 Jessops Riverside 800 Brightside Lane Sheffield S9 2RX United Kingdom
Leica Biosystems Nussloch GmbH	LBS	Germany	Heidelberger Strasse 17-19 69226 Nussloch Germany
Leica Biosystems Richmond Inc.	LBS	United States	C T Corporation System 208 South LaSalle Street, Suite 814 Chicago IL 60604 United States
Leica Microsystems Ltd. Shanghai	LBS	China	Building 1 No. 258 JIn Zang Road PUdong Shanghai 201206 China
Vision Finance Corporation Pty Ltd	LBS	Australia	495 Blackburn Road Mount Waverly VIC 3149 Australia
Vision Properties Pty Ltd	LBS	Australia	495 Blackburn Road Mount Waverly VIC 3149 Australia
Vision Systems Pty Ltd	LBS	Australia	495 Blackburn Road Mount Waverly VIC 3149 Australia
VSL Group Holdings Pty Ltd	LBS	Australia	495 Blackburn Road Mount Waverly VIC 3149 Australia
Cambridge Instruments LTD	LMS	United Kingdom	Lothbury House Newmarket Road Cambridge CB5 8PB United Kingdom
DH Acquisition Holding GmbH	LMS	Germany	Ernst-Leitz Strasse 17-37 35578 Wetzlar Germany
DH Holdings Germany GmbH	LMS	Germany	Ernst-Leitz-Strasse 17-37 D-35578 Wetzlar Germany
DH Holdings Germany LLC	LMS	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States

DH Verwaltungs GmbH	LMS	Germany	Ernst-Leitz Strasse 17-37 35578 Wetzlar Germany
DHRAT GmbH	LMS	Austria	Hernalser Hauptstrasse 219 1170 Wien Austria
DHRAT2 GmbH	LMS	Austria	Hernalser Hauptstrasse 219 1170 Wien Austria
DM FinCo GmbH	LMS	Germany	Ernst-Leitz Strasse 17-37 35578 Wetzlar Germany
DM HoldCo GmbH	LMS	Germany	Ernst-Leitz Strasse 17-37 35578 Wetzlar Germany
GHG GmbH	LMS	Germany	Ernst-Leitz-Strasse 17-37 D-35578 Wetzlar Germany
Kaltenbach & Voigt GmbH	LMS	Germany	Bismarckring 39 88400 Biberach/Riss Germany
KVHG GmbH	LMS	Germany	Ernst-Leitz Strasse 17-37 35578 Wetzlar Germany
Leica do Brasil Importação e Comercio Ltda	LMS	Brazil	Rua Julio Gonzales, 132 Barra Funda Sao Paulo SP Brazil
Leica Instruments (Singapore) Pte Limited	LMS	Singapore	12 Teban Gardens Crescent 608 924 Singapore
Leica Microsistemas - Instrumentos De Precisão, Sociedade Unipessoal Lda	LMS	Portugal	Av. do Forte, Nr.3 Edifício Suécia III- R/c, Dtº 2790-073 Carnaxide Portugal
Leica Microsistemas S.L.U	LMS	Spain	Nicaragua 46 08029 Barcelona Spain
Leica Microsystemes SAS	LMS	France	1 Rue du 1er Mai Immeuble Axe Sur Seine Hall C2, CS 50169 92752 Nanterre Cedex France

Leica Microsystems (Schweiz) AG	LMS	Switzerland	201 Max Schmidheiny Strasse CH-9435 Heerbrugg Switzerland
Leica Microsystems (SEA) Pte Ltd	LMS	Singapore	12 Teban Gardens Crescent 608924 Singapore
Leica Microsystems (Shanghai) Trading Co., Ltd.	LMS	China	Part C No. 127, North Fute Road Waigaoqiao Free Trade Zone Shanghai 200131 China
Leica Microsystems (UK) Limited	LMS	United Kingdom	19 Jessops Riverside 800 Brightside Lane Sheffield S9 2RX United Kingdom
Leica Microsystems A/S	LMS	Denmark	Akandevej 21 KD-2700 Bronshoj Denmark
Leica Microsystems AB	LMS	Sweden	c/o Cepheid, AB Box 1427, 171 27 Solna, Sweden Sweden
Leica Microsystems B.V.	LMS	Netherlands	Vlierweg 20, 1032LG Amsterdam Netherlands
Leica Microsystems Belgium BV	LMS	Belgium	Pegasuslaan 5 1831 Diegem Belgium
Leica Microsystems Cambridge Limited	LMS	United Kingdom	19 Jessops Riverside 800 Brightside Lane Sheffield S9 2RX United Kingdom
Leica Microsystems Canada	LMS	Canada	PwC Law LLP 18 York Street, Suite 2600 Toronto ON M5J 0B2 Canada
Leica Microsystems CMS GmbH	LMS	Germany	Ernst Letiz Strasse 17-37 35578 Wetzlar Germany
Leica Microsystems GmbH	LMS	Germany	Ernst-Leitz-Strasse 17 - 37 35578 Wetzlar Germany
Leica Microsystems Holdings GmbH	LMS	Germany	Ernst-Leitz Strasse 17-37 35578 Wetzlar Germany
Leica Microsystems Holdings Inc.	LMS	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States

Leica Microsystems Inc.	LMS	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
Leica Microsystems IR GmbH	LMS	Germany	Ernst-Leitz Strasse 17-37 35578 Wetzlar Germany
Leica Microsystems KK	LMS	Japan	1-29-9 Takadanobaba Shinjyuku-ku, Tokyo 169-0075 Japan
Leica Microsystems Limited 徠卡儀器有限公司	LMS	Hong Kong	22/F. AXA Tower, Landmark East 100 How Ming Street Kwun Tong Kowloon Hong Kong
Leica Microsystems NC, Inc.	LMS	United States	CT Corporation System 160 Mine Lake CT Suite 200 Raleigh NC 27615 United States
Leica Microsystems Pty Ltd	LMS	Australia	Suite 2, Level 3, Building A 11 Talavera Road Macquarie Park NSW 2113 Australia
Leica Microsystems S.R.L.	LMS	Italy	Via Emilia, 26 20090 Buccinasco MI Italy
Leica Microsystems Suzhou Technology Co., Ltd.	LMS	China	Bldg. B2, Rm. 503, Phase II, 88 Dongchang Rd. Suzhou Industrial Park, Jiang Su China
Leica Mikrosysteme GmbH	LMS	Austria	Hernalser Hauptstrasse 219 1170 Wien Austria
Leica Mikrosysteme Vertrieb GmbH	LMS	Germany	Ernst-Leitz-Straße 17-37 35578 Wetzlar Germany
Leica Technology Acquisition GmbH	LMS	Switzerland	Max Schmidheiny-Strasse 201 9435 Heerbrugg Switzerland
Leica Teknoloji Çözümleri Dağıtım Limited Şirketi	LMS	Turkey	FSM Mah. Poligon cad. no:8, Buyaka2 Sitesi 3 no'lu Blok, daire No:100 Umraniye- İstanbul Turkey
QG HoldCo GmbH	LMS	Germany	Ernst-Leitz-Strasse 17-37 D-35578 Wetzlar Germany

A/B Sciex AB	Sciex	Sweden	c/o Regus Östermalmstorg 1 114 42 Stockholm Sweden
A/B Sciex OY	Sciex	Finland	Biolinja 12 Turku FI-20750 Finland
AB Sciex (Hong Kong) Limited	Sciex	Hong Kong	Suites 406-409, 4/F., Three Pacific Place, 1 Queen's Road East Hong Kong
AB Sciex (Taiwan) Private Limited	Sciex	Taiwan	3F, No. 89, Sec. 2, Tiding Blvd. Neihu District Taipei City 11493 Taiwan
AB Sciex ApS	Sciex	Denmark	c/o Radiometer ApS Åkandevvej 21 2700 Bronshøj Denmark
AB Sciex AS	Sciex	Norway	c/o Regus Cort Adelersgate 16 Oslo 0254 Norway
AB Sciex Australia Pty Ltd	Sciex	Australia	495 Blackburn Road Mount Waverly VIC 3149 Australia
AB Sciex Austria GmbH	Sciex	Austria	Hernalser Hauptstrasse 219 1170 Wien Austria
AB Sciex Comércio de Instrumentos Laboratoriais Ltda.	Sciex	Brazil	Rua Julio Gonzalez, no. 132, 15o andar conjunto 91 Edificio Memorial Office Building Barra Funda CEP 01156-060 Brazil
AB Sciex EDC B.V.	Sciex	Netherlands	1e Tochtweg 11, 3rd Floor,2913LN, PO Box 330 2910 AH Nieuwerkerk aan den IJssel Netherlands
AB Sciex Finance B.V.	Sciex	Netherlands	1e Tochtweg 11, 3rd Floor,2913LN, PO Box 330 2910 AH Nieuwerkerk aan den IJssel Netherlands
AB Sciex Germany GmbH	Sciex	Germany	Landwehrstraße 54 64293 Darmstadt Germany

AB Sciex Holding LP	Sciex	Canada	PwC Law LLP 18 York Street, Suite 2600 Toronto ON M5J 0B2 Canada
AB Sciex Ireland Limited	Sciex	Ireland	Lismeehan, O' Callaghan Mills, Co. Clare Ireland
AB Sciex Kft	Sciex	Hungary	Vöröskereszt utca 8-10 Budapest 1222 Hungary
AB Sciex Korea Limited Company	Sciex	Korea, Republic of	15 FL. Hibrand Living Bldg. 16, Maeheon-ro, Seocho-gu Seoul, 137-924 Korea, Republic of
AB Sciex LLC	Sciex	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
AB Sciex LP	Sciex	Canada	PwC - Calgary 111 5th Avenue SW Suite 3100, East Tower Calgary AB T2P 5L3 Canada
AB Sciex Netherlands B.V.	Sciex	Netherlands	1e Tochtweg 11, 3rd Floor,2913LN, PO Box 330 2910 AH Nieuwerkerk aan den IJssel Netherlands
AB SCIEX NV	Sciex	Belgium	Suikerkaai 1 1500 Halle Belgium
AB Sciex Portugal – Sociedade Unipessoal Lda	Sciex	Portugal	Avenida do Forte n.º 3, Edifício Suécia III Piso 0, Sala 0.35 2794-041 Camaxide, Oeiras Portugal
AB Sciex Pte. Ltd.	Sciex	Singapore	33, #04-06, Marsiling Industrial Estate Road 3 739256 Singapore
AB Sciex Research GP Corp.	Sciex	Canada	71 Four Valley Dr. Concord ON L4K 4V8 Canada
AB Sciex S.r.l.	Sciex	Italy	Via Montenapoleone 8 20122 Milano Italy Italy
AB Sciex s.r.o	Sciex	Czech Republic	Zastrčená 1278/8, Praha 4 141 00 Czech Republic
AB Sciex Sales LP	Sciex	Canada	PwC Law LLP 18 York Street, Suite 2600 Toronto ON M5J 0B2 Canada

AB Sciex South Africa (Proprietary) Limited	Sciex	South Africa	Stand 1A, Growthpoint Business Park Tonnetti Street, Halfway House 1685 Johannesburg ZA South Africa
AB Sciex Spain, S.L.	Sciex	Spain	Calle Valgrande 8 Planta 1a Puerta 1 28108 Madrid Alcobendas Spain
AB Sciex Switzerland GmbH	Sciex	Switzerland	c/o Vosser Rechtsanwälte KIG Stadtturmstrasse 19 5401 Baden Switzerland
AB Sciex UK Limited	Sciex	United Kingdom	19 Jessops Riverside 800 Brightside Lane Sheffield S9 2RX United Kingdom
AB Sciex, S.A. de C.V.	Sciex	Mexico	Cuernavaca #106 Colonia Condesa Delagacion Cuauhtemoc Mexico City CP06140 Mexico
DCHL Holding BV	Sciex	Netherlands	1e Tochtweg 11, 3rd Floor, 2913LN, PO Box 330 2910 AH LN Nieuwerkerk aan den IJssel Netherlands
DH Operations B.V.	Sciex	Netherlands	1e Tochtweg 11, 3rd Floor, 2913LN, PO Box 330 2910 AH LN Nieuwerkerk aan den IJssel Netherlands
DH Technologies Development Pte Ltd.	Sciex	Singapore	33, #04-06, Marsiling Industrial Estate Road 3 739256 Singapore
Intabio, LLC	Sciex	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
K.K. AB Sciex	Sciex	Japan	7-35, Kita Shinagawa 4-chome Shinagawa-ku Tokyo Japan
Shanghai AB Sciex Analytical Instrument Trading Co., Ltd.	Sciex	China	Office 502, 5/F Bldg. 1 North Fuquan Road IBP Changning District Shanghai P.R. China 200335 China

Blueshift Biotechnologies, LLC	Molecular Devices	United States	C T Corporation System 818 West 7th Street Los Angeles CA 90017 United States
ID Business Solutions (Shanghai) Limited	Molecular Devices	China	Room 103, 1/F, No. 51, Lane 588 Shuping Road, Juyuan New Area Jiading District Shanghai China
ID Business Solutions Inc.	Molecular Devices	United States	1301 Marina Village Parkway Suite 320 Alameda CA 94501 United States
ID Business Solutions Limited	Molecular Devices	United Kingdom	2 Occam Court, Occam Road Surrey Research Park Guildford Surrey GU2 5QB United Kingdom
IDBS (Hong Kong) Limited	Molecular Devices	Hong Kong	21st Floor Edinburgh Tower The Landmark 15 Queen's Road Central Hong Kong Hong Kong
IDBS Asia-Pacific Limited	Molecular Devices	United Kingdom	2 Occam Court, Occam Road Surrey Research Park Guildford, Surrey England GU2 5QB United Kingdom
Inforsense Limited	Molecular Devices	United Kingdom	2 Occam Court, Occam Road Surrey Research Park Guildford, Surrey England GU2 5QB United Kingdom
Molecular Devices (Austria) GmbH	Molecular Devices	Austria	Urstein Süd 17, 5412 Puch bei Hallein Austria
Molecular Devices (France) SAS	Molecular Devices	France	Immeuble Rimbaud 22, avenue des Nations Zone Artisanale Paris Nord II 93420 Villepinte France
Molecular Devices (Germany) GmbH	Molecular Devices	Germany	Sauerbruch Strasse 50 81377 Munich Germany
Molecular Devices (Hong Kong) Limited	Molecular Devices	Hong Kong	21st Floor Edinburgh Tower The Landmark 15 Queen's Road Central Hong Kong Hong Kong

Molecular Devices (Shanghai) Corporation	Molecular Devices	China	Room 505, Building 1, No. 389 Gang Ao Road Wai Gao Qiao Free Trade Zone Shanghai City China
Molecular Devices (Shanghai) Limited	Molecular Devices	China	2-1, No. 555 Yuanxi Road Nanhui Industrial Park Shanghai City China
Molecular Devices (UK) Limited	Molecular Devices	United Kingdom	660-665 Eskdale Road Winnersh Triangle Wokingham Berkshir RG41 5TS United Kingdom
Molecular Devices Instrumentação do Brasil Ltda.	Molecular Devices	Brazil	Rua Júlio González, 132 14th floor, conjunto 81 Memorial Office Building, Barra Funda Sao Paulo 01156-060 Brazil
Molecular Devices Japan KK	Molecular Devices	Japan	NihonbashiBakuro-cho Yuraku Bldg., 7F 2-7-8 Nihonbashi Bakuro-cho Chuo-ku Tokyo 103-0002 Japan
Molecular Devices Korea, LLC	Molecular Devices	Korea, Republic of	3rd floor, Gangnam Bldg., 1321-1 Seocho-dong, Seocho-gu, Seoul Korea
Molecular Devices, LLC	Molecular Devices	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
Skyland Analytics Inc.	Molecular Devices	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
SubX, Inc.	Molecular Devices	United States	3rd Fl., 2 Cotton Street Portland ME 04101 United States
Advanced Vision Technology (A.V.T.) Ltd.	Esko	Israel	6 Hanagar St. Hod Hasharon 45241 Israel
Advanced Vision Technology A.V.T (Germany) GmbH	Esko	Germany	32 Donnersberger Strasse 80634 Munich Germany
AVT EMEA CVBA	Esko	Belgium	Kortrijksesteenweg 1095 9051 Ghent Belgium
AVT, Advanced Vision Technology, Inc.	Esko	United States	1209 Orange Street Wilmington DE 19801 United States

Blue Software (UK) Limited	Esko	United Kingdom	19 Jessops Riverside 800 Brightside Lane Sheffield S9 2RX United Kingdom
Blue Software Acquisition Corporation	Esko	United States	8430 W. Bryn Mawr Ave. Suite 1100 Chicago IL 60631 United States
Blue Software Holdings Corporation	Esko	United States	1209 Orange Street Corporation Trust Center Wilmington DE 19801 United States
Blue Software, LLC	Esko	United States	1209 Orange Street Corporation Trust Center Wilmington DE 19801 United States
Enfocus BV	Esko	Belgium	Raymonde de Larochelaan 13 9051 Gent Belgium
Esko BV	Esko	Belgium	Raymonde de Larochelaan 13 9051 Gent Belgium
Esko Finance BV	Esko	Belgium	Raymonde de Larochelaan 13 9051 Gent Belgium
Esko Graphics Mexico, S. De R.L. de C.V.	Esko	Mexico	Avenida Coyoacan 1213, Colonia del Valle, Benito Juarez Distrito Federal, MX 03340 Mexico
Esko Graphics UK Limited	Esko	United Kingdom	19 Jessops Riverside 800 Brightside Lane Sheffield S9 2RX United Kingdom
Esko Iberia S.L	Esko	Spain	Ronda General Mitre, 28-30 08017 Barcelona Spain
ESko Software BV	Esko	Belgium	Raymonde de Larochelaan 13 9051 Gent Belgium
Esko Stonecube Limited	Esko	United Kingdom	19 Jessops Riverside 800 Brightside Lane Sheffield S9 2RX United Kingdom
Esko-Graphics (Shanghai) Trading Company Ltd.	Esko	China	Room 325 Xinxing Building No. 2005 North Yanggao Road Waigaoqiao District Shanghai China 200131 China
Esko-Graphics BV	Esko	Belgium	Raymonde de Larochelaan 13 9051 Gent Belgium

Esko-Graphics Co. Ltd.	Esko	Japan	West Wing 6F, Telecom Center Bldg. 2-5-10 Aomi, Koto-ku Tokyo 135-0064 Japan
Esko-Graphics Danmark A-S	Esko	Denmark	Akandevej 21 KD-2700 Bronshoj Denmark
Esko-Graphics do Brasil Comércio de Produtos Tecnológicos Ltda.	Esko	Brazil	Rua Professor Aprigio Gonzaga 78 Andar 10 CJTS 101 102 103 E 104 São Judas São Paulo 04303-000 Brazil
Esko-Graphics GesmbH	Esko	Austria	Lehmanngasse 7 1230 Vienna Austria
ESKO-GRAPHICS GmbH	Esko	Switzerland	c/o GM Rechtsanwälte St. Jakob-Strasse 37 9000 St. Gallen Switzerland
Esko-Graphics GmbH [Germany]	Esko	Germany	Heerskamp 6 25524 Itzehoe Germany
Esko-Graphics Imaging GmbH	Esko	Germany	Heerskamp 6 25524 Itzehoe Germany
Esko-Graphics Inc.	Esko	United States	C T Corporation System 289 S. Culver St. Lawrenceville GA 30046
Esko-Graphics India Private Limited	Esko	India	Esteem Tower, No. 71, Railway Parallel Road Kumara Park West Bangalore 560 020 India
Esko-Graphics Pte Ltd.	Esko	Singapore	1 Kaki Bukit View #04-15/17 Techview 415941 Singapore
Esko-Graphics S.A.S.	Esko	France	ALLEE DU MOULIN PINARD ZI NORD 87100 Limoges France
Esko-Graphics S.r.l	Esko	Italy	Viale Sondrio, 2 20124 Milan Italy

Esko-Graphics s.r.o.	Esko	Czech Republic	Bieblova 19/1227 Praha 5 Smichov 150 00 Czech Republic
Graphics Microsystems, LLC	Esko	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
Viki Solutions Inc.	Esko	Canada	B701-810 Humbolt Victoria BC V8V5B1 Canada
Arista Molecular, Inc.	BEC - Diagnostics	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
BC Distribution B.V.	BEC - Diagnostics	Netherlands	Pelmolenlaan 15 3447GW Woerden Netherlands
BC Japan LLC	BEC - Diagnostics	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
BCK Holdings S.à.r.l.	BEC - Diagnostics	Luxembourg	1B Heienhaff, L-1736 Senningerberg Grand-Duchy of Luxembourg Luxembourg
Beckman Coulter (Israel) Ltd.	BEC - Diagnostics	Israel	2 Pekeris St Park Tamar Rehovot 7670202 Israel
Beckman Coulter (U.K.) Pension Trustees Limited	BEC - Diagnostics	United Kingdom	Oakley Court, Kingsmead Business Park London Road High Wycombe HP11 1JU United Kingdom
Beckman Coulter AB	BEC - Diagnostics	Sweden	Archimedesvägen 7 Box 11156 S-168 6 161 11 Bromma Sweden
Beckman Coulter Argentina S.A.	BEC - Diagnostics	Argentina	Guemes 4168 Villa Martelli Buenos Aires B1603BEN Argentina
Beckman Coulter Australia Pty Ltd	BEC - Diagnostics	Australia	23-27 Chaplin Drive Lane Cove West NSW 2066 Australia
Beckman Coulter Biomedical GmbH	BEC - Diagnostics	Germany	Sauerbruch Strasse 50 81377 Munich Germany

Beckman Coulter Biyomedikal Ürünler Sanayi ve Ticaret Limited Şirketi	BEC - Diagnostics	Turkey	Askent Sokak No:3/A Kosifler İş Merkezi A Blok K:18 İçerenköy Atasehir Istanbul 34752 Turkey
Beckman Coulter Canada LP	BEC - Diagnostics	Canada	PwC Law LLP 18 York Street, Suite 2600 Toronto ON M5J 0B2 Canada
Beckman Coulter Ceska republika s.r.o.	BEC - Diagnostics	Czech Republic	Praha 10 Radiová 1122/1 PŠC 10200 Prague Czech Republic
Beckman Coulter Chile Limitada	BEC - Diagnostics	Chile	Piso 11 Floor Isidora Goyenechea 2939 Las Condes Santiago Chile
Beckman Coulter Colombia S.A.S.	BEC - Diagnostics	Colombia	CL 13 NO. 60 49/61, BOGOTA D.C. Colombia
Beckman Coulter Commercial Enterprise (China) Co., Ltd.	BEC - Diagnostics	China	4/F, 518 North Fu Quan Road Chang Ning District 450 Fu Shan Shanghai City China
Beckman Coulter d.o.o.	BEC - Diagnostics	Croatia	Av. V. Holjevca 40 Zagreb 10020 Croatia
Beckman Coulter Danmark ApS	BEC - Diagnostics	Denmark	Åbenrå 23, Ground floor DK-1124 Copenhagen K Denmark
Beckman Coulter de Mexico, S.A. de C.V.	BEC - Diagnostics	Mexico	Suite 396 Ave. Popocatepetl Mexico City 03340 Mexico
Beckman Coulter de Venezuela, S.A.	BEC - Diagnostics	Venezuela, Bolivarian Republic of	Avenida Anauco, Quinta Yumare #23-11 San Bernardino Caracas 1010 DF Venezuela
Beckman Coulter Diagnostics Limited	BEC - Diagnostics	Ireland	Lismeehan, O' Callaghan Mills, Co. Clare Ireland
Beckman Coulter do Brasil Comércio e Importação de Produtos de Laboratório Ltda.	BEC - Diagnostics	Brazil	Calçada Aldebarã 39 Centro de Apoio 2 Santa de Parnaíba São Paulo 06541-055 Brazil

Beckman Coulter España, S.A.U.	BEC - Diagnostics	Spain	Plaza Europa 41-43 L'Hospitalet de Llobregat 08902 Barcelona Spain
Beckman Coulter Eurocenter S.A.	BEC - Diagnostics	Switzerland	22 Rue Juste-Olivier P.O. Box 1044 Nyon-1 1260 Switzerland
Beckman Coulter Finland Oy	BEC - Diagnostics	Finland	c/o Tilitoimisto Hörhammer Oy Henry Fordin katu 5 C Helsinki 00150 Finland
Beckman Coulter Foundation	BEC - Diagnostics	United States	C T Corporation System 818 West 7th Street Los Angeles CA 90017 United States
Beckman Coulter France S.A.S.	BEC - Diagnostics	France	22 avenue des Nations Immeuble le Rimbaud 93420 Villepinte France
Beckman Coulter Genomics Inc.	BEC - Diagnostics	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
Beckman Coulter GmbH	BEC - Diagnostics	Austria	Giefinggasse 6, Top 8 1210 Vienna Austria
Beckman Coulter GmbH	BEC - Diagnostics	Germany	Europark Fichtenhain B 13 47807 Krefeld Germany
Beckman Coulter Holdings G.K.	BEC - Diagnostics	Japan	TOC Ariake West Tower 3-5-7 Ariake Tokyo, Koto-ku 135-0063 Japan
Beckman Coulter Hong Kong Limited	BEC - Diagnostics	Hong Kong	22/F. AXA Tower, Landmark East 100 How Ming Street Kwun Tong Kowloon Hong Kong
Beckman Coulter India Private Limited	BEC - Diagnostics	India	Unit Nos.TF-B-07(A) to15, Third Floor, B Wing Art Guild House, Phoenix Market City LBS Road, Kurla (W) Mumbai Maharashtra 400070 India

Beckman Coulter International S.A.	BEC - Diagnostics	Switzerland	22 Rue Juste-Olivier P.O. Box 1059 CH-1260 Nyon 1 Switzerland
Beckman Coulter International Trading (Shanghai) Co., Ltd.	BEC - Diagnostics	China	379 Debao Road East A Part of 1st/F, factory building No. 4 Waigaoqiao Bonded Area Shanghai China
Beckman Coulter Ireland Inc.	BEC - Diagnostics	Ireland	Lismeehan, O' Callaghan Mills, Co. Clare Ireland
Beckman Coulter K.K.	BEC - Diagnostics	Japan	TOC Ariake West Tower 3-5-7 Ariake Tokyo, Koto-ku 135-0063 Japan
Beckman Coulter Korea Ltd.	BEC - Diagnostics	Korea, Republic of	3rd Floor Suseo-dong Hyosung Sudeo Building, 281 Gwangpyeong-ro, Gangnam- gu Seoul 06349 Korea, Republic of
Beckman Coulter Laboratory Systems (Suzhou) Co. Ltd.	BEC - Diagnostics	China	181 Su Hong Road West Sino Singapore Suzhou Industrial Park Suzhou CN215021 China
Beckman Coulter Limited Liability Company	BEC - Diagnostics	Russian Federation	Stanislavskogo Street 21, Bldg. 3 Bus Center Fabrika Stanis Moscow 109004 Russia
Beckman Coulter Magyarország Kft	BEC - Diagnostics	Hungary	1038 Budapest Papírgyár u. 58-59 Budapest Hungary
Beckman Coulter Malaysia Sdn Bhd	BEC - Diagnostics	Malaysia	No. 18, Jalan Tandang 51/205A Seksyen 51, 46050 Petaling Jaya Malaysia
Beckman Coulter Mishima K.K.	BEC - Diagnostics	Japan	454-32 Higashino, Nagaizumi- cho Sunto-gun Shizuoka-ken 411-0931 Japan
Beckman Coulter Nederland B.V.	BEC - Diagnostics	Netherlands	Pelmolenlaan 15 3447GW Woerden Netherlands

Beckman Coulter Nippon G.K.	BEC - Diagnostics	Japan	TOC Ariake West Tower 3-5-7 Ariake Tokyo, Koto-ku 135-0063 Japan
Beckman Coulter NZ	BEC - Diagnostics	New Zealand	Unit 3, 33 Spartan Road Takanini Takanini 2105 New Zealand
Beckman Coulter Polska sp. z.o.o.	BEC - Diagnostics	Poland	Al. Jerozolimskie 181 A 02-222 Warsaw Poland
Beckman Coulter Portugal Unipessoal Lda.	BEC - Diagnostics	Portugal	Avenida do Forte n.03 Edificio Suecia III Pisa 0, Sala 0.35 Carnaxide, Portugal Portugal
Beckman Coulter Puerto Rico Inc.	BEC - Diagnostics	United States	C T Corporation System 208 South LaSalle Street, Suite 814 Chicago IL 60604 United States
Beckman Coulter Saudi Arabia Co. Ltd.	BEC - Diagnostics	Saudi Arabia	OLIA ST. Raseel Building office # 8 Riyadh 11335 Saudi Arabia
Beckman Coulter Servicios, S. de R.L. de C.V.	BEC - Diagnostics	Mexico	Suite 396 Av. Popocatepetl, Col Gral Pedro Ma. Anauya Delg Mexico City 03340 Mexico
Beckman Coulter Singapore Pte. Ltd.	BEC - Diagnostics	Singapore	7 Straits View #12-00 Marina One East Tower Singapore 018936 Singapore
Beckman Coulter Slovenska republika s.r.o.	BEC - Diagnostics	Slovakia	Digital Park II, Einsteinova 23 Bratislava SK85101 Slovakia
Beckman Coulter South Africa (Proprietary) Limited	BEC - Diagnostics	South Africa	Stand 1A Growth Point Business Park Tonetti Street Halfway House Johannesburg ZA 1685 South Africa
Beckman Coulter Srl	BEC - Diagnostics	Italy	Via Roma 108 - Palazzo F/1 20060 Cassina De' Pecchi Milan Italy

Beckman Coulter Taiwan Inc.	BEC - Diagnostics	United States	C T Corporation System 818 West 7th Street Los Angeles CA 90017 United States
Beckman Coulter United Kingdom Limited	BEC - Diagnostics	United Kingdom	Oakley Court, Kingsmead Business Park London Road High Wyco Buckinghamshire GB HP111JU United Kingdom
Beckman Coulter, Inc.	BEC - Diagnostics	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
Beckman Coulter, S.L.U.	BEC - Diagnostics	Spain	Plaza Europa 41-43 L'Hospitalet de Llobregat 08908 Barcelona Spain
Beckman Holdings Limited	BEC - Diagnostics	United Kingdom	Oakley Court, Kingsmead Business Park London Road High Wycombe Buckinghamshire HP11 1JU United Kingdom
Coulter Electronics Pty. Limited	BEC - Diagnostics	Australia	23-27 Chaplin Drive Lane Cove NSW 2066 Australia
DH II Luxembourg Sarl	BEC - Diagnostics	Luxembourg	1B Heienhaff, L-1736 Senningerberg Grand-Duchy of Luxembourg Luxembourg
DHR Holding Spain S.L.	BEC - Diagnostics	Spain	L'Hospitalet de Llobregat Plaza de Europa 41-34 Torre Realia 4a planta 08908 Barcelona Spain
Diagnostic Systems Laboratories Australia Pty Ltd	BEC - Diagnostics	Australia	23-27 Chaplin Drive Lane Cove West NSW 2066 Australia
DTIL Ireland Holdings Ltd	BEC - Diagnostics	Ireland	Lismeehan, O' Callaghan Mills, Co. Clare Ireland
DX Laboratuvar Cozumleri A.S.	BEC - Diagnostics	Turkey	Askent Sokak No: 3 A Kosifler İş Merkezi A Blok K 8 34752 İçerenköy Istanbul Turkey Turkey
Fullerton LLC	BEC - Diagnostics	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
Hybrigenetics Cancer Research Incorporated	BEC - Diagnostics	United States	C T Corporation System 818 West 7th Street Los Angeles CA 90017 United States

Hybrigenetics Cancer Research Limited	BEC - Diagnostics	United States	C T Corporation System 818 West 7th Street Los Angeles CA 90017 United States
Hybritech Incorporated	BEC - Diagnostics	United States	C T Corporation System 818 West 7th Street Los Angeles CA 90017 United States
Hybritech International, Incorporated	BEC - Diagnostics	United States	C T Corporation System 818 West 7th Street Los Angeles CA 90017 United States
IAF Corporation Pty Limited	BEC - Diagnostics	Australia	23-27 Chaplin Drive Lane Cove West NSW 2066 Australia
Immunotech Sro	BEC - Diagnostics	Czech Republic	Radiova 1 Prague 10 Czech Republic
Inhibin Pty Limited	BEC - Diagnostics	Australia	c/o Scott Associates Suite 203 Level 2 379-383 Pitt Street Sydney NSW 2000 Australia
Iris International Europe GmbH	BEC - Diagnostics	Switzerland	Hardturmstrasse 161 8005 Zurich Switzerland
IRIS International, Inc.	BEC - Diagnostics	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
KFL Finance Ltd	BEC - Diagnostics	Ireland	Lismeehan, O' Callaghan Mills, Co. Clare Ireland
KHL Ireland Investment Holdings Ltd	BEC - Diagnostics	Ireland	Lismeehan, O' Callaghan Mills, Co. Clare Ireland
Lumigen, Inc.	BEC - Diagnostics	United States	The Corporation Company 40600 Ann Arbor Road E Suite 201 Plymouth MI 48170 United States
Normand Info SAS	BEC - Diagnostics	France	7 rue Frederic Degeorge, Bat B, CS 10451 62028 Arras Cedex France
QHC Ireland Finance Limited	BEC - Diagnostics	Ireland	Lismeehan, O' Callaghan Mills, Co. Clare Ireland

Beckman Coulter (Dalian) Software Co., Ltd.	BEC - Life Sciences	China	No. 1406, Building A, No. 32 Huoju Road, Dalian High-Tech Industrial Zone, Dalian, The People's Republic of China (PRC China
Beckman Coulter Biotechnology (Suzhou) Co., Ltd.	BEC - Life Sciences	China	Unit 101, 201, 301, Building 4 218 Sang Tian Street Suzhou Industry Park Suzhou City China
Immunotech SAS	BEC - Life Sciences	France	130 Avenue Jean de Lattre de Tassigny - BP 177 13276 Marseille Ce France
Labcyte Europe Ltd.	BEC - Life Sciences	Ireland	Unit 12, Enterprise Centre Pearse Street Dublin 2 Ireland
Labcyte GmbH	BEC - Life Sciences	Germany	c/o LM Leinauer Müller & Partner Paul-Gerhardt-Allee 50 81245 München Germany
Labcyte Inc.	BEC - Life Sciences	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
Labcyte Ltd.	BEC - Life Sciences	United Kingdom	19 Jessops Riverside 800 Brightside Lane Sheffield S9 2RX United Kingdom
m2p-labs Limited	BEC - Life Sciences	Hong Kong	UNIT 1213, 12/F, WING ON PLAZA, 62 MODY ROAD, TSMIT SHA TSUI EAST, KOWLOON Hong Kong
Pantone Japan Co, Ltd.	X-Rite	Japan	West wing 6F, Telecom Center Bldg. 2-5-10, Aomi, Koto-ku, Tokyo 135-0064 Japan
Pantone Japan, Inc.	X-Rite	United States	The Corporation Trust Company 820 Bear Tavern Road West Trenton NJ 08628 United States
Pantone LLC	X-Rite	United States	The Corporation Trust Company 820 Bear Tavern Road West Trenton NJ 08628 United States

X-Rite (Shanghai) Color Management Col, Ltd.	X-Rite	China	6/F, Building#1 , IBP Phase 2 518 Fuquan Road North Changning District Shanghai 200335 China
X-Rite Asia Pacific Limited	X-Rite	Hong Kong	Suite 2801, 28th Floor, AXA Tower, Landmark East 100 How Ming Street Kwun Tong Kowloon Hong Kong
X-Rite Europe GmbH	X-Rite	Switzerland	Althardstrasse 70 CH--8105 Regensdorf Switzerland
X-Rite Global, Incorporated	X-Rite	United States	The Corporation Company 40600 Ann Arbor Road E Suite 201 Plymouth MI 48170 United States
X-Rite GmbH	X-Rite	Germany	Fraunhoferstr 14 82152 Martinsried Germany
X-Rite Holdings, Inc.	X-Rite	United States	The Corporation Company 40600 Ann Arbor Road E Suite 201 Plymouth MI 48170 United States
X-Rite India Private Limited	X-Rite	India	115, Free Press House, 11th Floor, Free Press Journal Marg, Nariman Point Mumbai India 400 021 India
X-Rite Italy S.R.L.	X-Rite	Italy	Via Luigi Bralle 69-69 I-59100 Prato Italy
X-Rite Limited	X-Rite	United Kingdom	19 Jessops Riverside 800 Brightside Lane Sheffield S9 2RX United Kingdom
X-Rite Mediterranee SARL	X-Rite	France	43 rue du Saule Trapu 91300 Massy France
X-Rite, Incorporated	X-Rite	United States	The Corporation Company 30600 Telegraph Road, Suite 2345 Bingham Farms MI 48025 United States

DH Life Sciences Switzerland Sàrl	Pall	Switzerland	Avenue de Tivoli 3 CH-1700 Fribourg Switzerland
Filter Specialists Inc. de Mexico S. de R.L. de C.V.	Pall	Mexico	Camino Temixco a Emiliano Zapata km3 interior 4a y 4b, exterior S/N, Colonia Palo Escrito CP 62760, Emiliano Zapata, Morelos MEXICO D.F. MX Mexico
Filter Specialists India Private Limited	Pall	India	C/O Kalyaniwalla & Misty Plant 13, Extension Off, Eastern Express Highway Pirojsha Ngr, Vikroli (E) Mumbai Maharashtra 400079 India
FSI Leasing, LLC	Pall	United States	2200 Pennsylvania Ave., NW, Suite 800, Washington, DC 20037
Gelman Sciences Inc.	Pall	United States	The Corporation Company 40600 Ann Arbor Road E Suite 201 Plymouth MI 48170 United States
Nihon Pall Ltd.	Pall	Japan	SEE MAIN OFFICE
Nihon Pall Manufacturing Limited	Pall	Japan	SEE MAIN OFFICE
PALL (CANADA) ULC	Pall	Canada	2800 Park Place 666 Burrard Street Vancouver BC V6C2Z7 Canada
Pall (China) Co., Ltd.	Pall	China	No 12 Hongda Nanlu Beijing Economic Technological Development Area Beijing 100176 China
Pall (Malaysia) Sdn Bhd	Pall	Malaysia	Level 11, 1 Sentral Jalan Rakyat Kuala Lumpur Sentral 50470 Kuala Lumpur Malaysia
Pall (Philippines) Inc.	Pall	Philippines	16th Floor, Tower II, Insular Life Corporate Centre, Insular Drive, Filinvest Corporate City, Alabang 1781 Muntinlupa City Philippines Philippines
Pall (Schweiz) GmbH	Pall	Switzerland	Schaeferweg 16 4057 Basel Switzerland

Pall Aeropower Corporation	Pall	United States	The Corporation Trust Company Corporation Trust Center 1209 Orange Street Wilmington DE 19801 United States
PALL Arabia for Trading	Pall	Saudi Arabia	Al-Turki Business Park (2nd Floor) 7244 King Saud Road, Ad Dawhah Al Janubiyah Dist., Unit No. 55 Dhahran 34455-3955 Saudi Arabia
Pall Artelis BV	Pall	Belgium	Reugelstraat 2 3320 Hoegaarden Belgium Belgium
Pall Asia Holdings, Inc.	Pall	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
Pall Asia International Ltd.	Pall	Hong Kong	21st Floor Edinburgh Tower The Landmark 15 Queen's Road Central Hong Kong Hong Kong
PALL AUSTAR Lifesciences Limited	Pall	Hong Kong	c/o 14/F, San Toi Building, 137 - 139 Connaught Road Central Hong Kong Hong Kong
Pall Australia Pty Ltd	Pall	Australia	1-2 Wandarri Court Cheltenham, Victoria Melbourne VIC 3192 Australia
Pall Austria Filter Ges.m.b.h.	Pall	Austria	Handelskai 94-96 / Top 422 A-1200 Vienna Austria
Pall Belgium BV	Pall	Belgium	Reugelstraat 2 3320 Hoegaarden Belgium
Pall Biotech (Beijing) Co., Ltd.	Pall	China	Yanqi Road, No 15 Yanqi Economic Development Zone Huairou Beijing China
Pall Corporation	Pall	United States	C T Corporation System 111 Eighth Avenue New York NY 10011 United States
Pall Corporation Filtration & Separations (Thailand) Ltd	Pall	Thailand	SEE MAIN OFFICE

Pall do Brasil Ltda.	Pall	Brazil	Avenue Luigi Papaiz no. 239 Administrative Block, 1st Floor, Part B, Industrial and Commercial Center, Campanario Diadema Sao Paulo 09.931-610 Brazil
Pall Espana, S.A.	Pall	Spain	Isla De Corcega 2 28100 Alcobendas Spain
Pall Eurasia L.L.C.	Pall	Russian Federation	Stanislavskogo Street 21 Buld 3 Room 1 109004 Moscow Russian Federation
Pall Europe Limited	Pall	United Kingdom	5 Harbourgate Business Park Southampton Road Portsmouth Hampshire PO6 4BQ United Kingdom
Pall European Holdings Limited	Pall	United Kingdom	5 Harbourgate Business Park Southampton Road Portsmouth Hampshire PO6 4BQ United Kingdom
Pall Exekia S.A.S.	Pall	France	Usine de bazet 65460 Bazet France
Pall Export Sales Ltd	Pall	Jamaica	c/o Myers, Fletcher & Gordon Park Place, 21 East St, PO Box 162 Kingston Jamaica
Pall Filter Specialists, Inc.	Pall	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
Pall Filtersystems GmbH	Pall	Germany	Planiger Strasse 137 55543 Bad Kreuznach Germany
Pall Filtration and Separations Group Inc.	Pall	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
Pall Filtration Pte Ltd	Pall	Singapore	7 Straits View #12-00 Marina One East Tower Singapore 018936 Singapore
Pall France SAS	Pall	France	3 Rue des Gaudines 78100 Saint Germain en Laye France

Pall GeneDisc Technologies SAS	Pall	France	1 rue du Courtil Parc d"Affaires Cicea Bat 1 35170 Bruz France
Pall GmbH	Pall	Germany	Philipp-Reis-Str. 6 D-63303 Dreieich Germany
Pall India Pvt Ltd	Pall	India	Unit No. TF-B-08 Art Guild House, CTS No. 124-A Off LBS Marg, Near Pheonix Mall Kurla (West) Mumbai 400070 India
Pall International Holdings, Inc.	Pall	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
Pall International Sarl	Pall	Switzerland	Avenue de Tivoli 3 CH-1700 Fribourg Switzerland
Pall Italia S.R.L.	Pall	Italy	Via Emilia, 26 20090 Buccinasco MI Italy
Pall Korea Ltd	Pall	Korea, Republic of	Dongwon Building 333 4th Flr Youngdong-daero Gangnam-gu, Seoul 06188 Korea, Republic of
Pall Life Sciences Belgium BV	Pall	Belgium	Reugelstraat 2 3320 Hoegaarden Belgium
Pall Life Sciences Mexico, S. de R.L. de C.V.	Pall	Mexico	Colinas 11731 Parque Industrial El Florido Baja California C.P. 22224 Mexico
Pall Life Sciences Puerto Rico, LLC	Pall	United States	CT Corporation System 361 San Francisco St. 4th Floor San Juan PR 00901 United States
Pall Manufacturing UK Limited	Pall	United Kingdom	5 Harbourgate Business Park Southampton Road Portsmouth Hampshire PO6 4BQ United Kingdom
Pall Medistad B.V.	Pall	Netherlands	Nijverheidsweg 1 1671GC Medemblik Netherlands

Pall Middle East FZE	Pall	United Arab Emirates	Office No. LB14330 Jebel Ali Dubai United Arab Emirates
Pall Modultechnik GmbH	Pall	Germany	Planiger Strasse 137 55543 Bad Kreuznach Germany
Pall Netherlands B.V.	Pall	Netherlands	Nijverheidsweg 1 1671GC Medemblik Netherlands
Pall New Zealand Ltd	Pall	New Zealand	188 Quay Street Auckland New Zealand
Pall Norden AB	Pall	Sweden	Pall Norden AB c/o Medicon Village AB Scheelevagen 2 223 81 Lund Sweden Sweden
Pall Norge AS	Pall	Norway	Skjærdalen 2 3533 Tyrstrand 0605 Ringerike Norway
Pall Poland Sp. z.o.o.	Pall	Poland	ul. Wenecka 12 03-244 Warsaw Warsaw Poland
Pall S. A. R. L. A. U.	Pall	Morocco	Sidi Maarouf Business Center Casablanca Nearshore Park 1100 Bvd. Al Quods 20190 Sidi Maarouf Casablanca Morocco
Pall Saudi International Limited	Pall	Saudi Arabia	Al-Turki Business Park (Loft 1) Dhahran 34455 - 4088 King Saud Road Doha 7448 Unit No. 3 Saudi Arabia
Pall Singapore Taiwan Branch Holding Company Pte Ltd	Pall	Singapore	1 Science Park Road, #05 - 09 East Wing The Capricorn, Singapore Science Park II Singapore 117528 Singapore
Pall South Africa (Pty)Ltd	Pall	South Africa	16th Floor South Tower The Towers Heerengracht Cape Town Western Cape 8001 South Africa

Pall Technologies S. A.	Pall	Argentina	Riobamba 1236, 8C Buenos Aires C1116ABJ Argentina
Pall Technology UK Limited	Pall	United Kingdom	5 Harbourgate Business Park Southampton Road Portsmouth Hampshire PO6 4BQ United Kingdom
Pall UK Holdings Limited	Pall	United Kingdom	19 Jessops Riverside 800 Brightside Lane Sheffield S9 2RX United Kingdom
PLLN Inc.	Pall	United States	1209 Orange Street Corporation Trust Center Wilmington DE 19801 United States
poromembrane GmbH	Pall	Germany	Alfred-Schefenacker-Strasse 1 71409 Schwaikheim Germany
Precision Nanosystems Singapore Pte. Ltd.	Pall	Singapore	7 Straits View #12-00 Marina One East Tower Singapore 018936 Singapore
Precision Nanosystems UK Limited	Pall	United Kingdom	19 Jessops Riverside 800 Brightside Lane Sheffield S9 2RX United Kingdom
Precision Nanosystems ULC	Pall	Canada	Suite 1400, 250 Howe Street Vancouver BC V6C 357 Canada
Precision Nanosystems US, Inc.	Pall	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
PT Pall Filtration Indonesia	Pall	Indonesia	SEE MAIN OFFICE
Cepheid	Cepheid	United States	C T Corporation System 818 West 7th Street Los Angeles CA 90017 United States
Cepheid (Shanghai) Commerce and Trade Co., Ltd.	Cepheid	China	R1806-1807, No. 1602 West Zhongshan Road, Hongwell International Plaza, Shanghai 200235 China
Cepheid (Shanghai) Diagnostics Co., Ltd.	Cepheid	China	Room A02, Floor 4 No. 207 North Fu Te Road Shanghai China China

Cepheid (Suzhou) Medical Science Co. Ltd.	Cepheid	China	Unit 201, Building 27, Phase II of BioBay No.218 Sangtian Street Suzhou Industrial Park Jiangsu Pilot Free Trade Zone Suzhou Are 215123 China
Cepheid AB	Cepheid	Sweden	Box 1427, 171 27 Solna Sweden
Cepheid Benelux BV	Cepheid	Belgium	Bellevue 5, box 1001 9050 Ghent Belgium
Cepheid Brasil Exportacao e Comercio de Produtos de Diagnosticos Ltda.	Cepheid	Brazil	Rua Luigi Galvani, 200 - conjunto 51 5.andar 04575-020 - São Paulo – SP Brasil Brazil
Cepheid Canada ULC	Cepheid	Canada	PwC - Vancouver 250 Howe Street Suite 1400 Vancouver BC V6C3S7 Canada
Cepheid Denmark ApS	Cepheid	Denmark	2 Oslo Plads, 2100 Copenhagen DK Denmark
Cepheid Diagnostics & Lifesciences (Private) Limited	Cepheid	Zimbabwe	CEP Zimbabwe 203 Josiah Chinamano Avenue Harare Zimbabwe
Cepheid Diagnostics And LifeSciences Limited	Cepheid	Kenya	Nine Riverside Building, Riverside Drive, P.O. Box 12342 - 00100 – G.P.O. Nairobi Kenya
Cepheid Europe SAS	Cepheid	France	Vira Solelh, 81470 Maurens-Scopont France
Cepheid Finland Oy	Cepheid	Finland	Fabianinkatu 23 Helsinki 00130 Finland
Cepheid GK	Cepheid	Japan	703 Burex Toranomom, 2-7-5 Toranomom Minato-ku, Tokyo 105-0001 Japan Japan
Cepheid GmbH	Cepheid	Germany	Europark Fichtenhain A 4, 47807 Krefeld Germany
Cepheid HBDC	Cepheid	United States	C T Corporation System 818 West 7th Street Los Angeles CA 90017 United States
Cepheid HBDC SAS	Cepheid	France	Vira Solelh, 81470 Maurens-Scopont France

Cepheid Holdings Pty Limited	Cepheid	Australia	495 Blackburn Road Mount Waverly VIC 3149 Australia
Cepheid Iberia SLU	Cepheid	Spain	C/ Caleruega, 81 28033 Madrid Spain
Cepheid India Private Limited	Cepheid	India	Unit no: DSM 214 & 215, 2nd Floor, DLF Towers Shivaji Marg, Najafgarh Road, New Delhi-10015 India
Cepheid Limited	Cepheid	Hong Kong	21st Floor, Edinburgh Tower, The Landmark 15 Queen's Road Central Hong Kong Hong Kong
Cepheid Netherlands BV	Cepheid	Netherlands	1e Tochtweg 11 2913LN Nieuwerkerk aan den IJssel Netherlands
Cepheid Norway AS	Cepheid	Norway	c/o Pall Norge AS Skjaerdalen 2 3533 TYRISTRAND 3007 RINGERIKE Norway
Cepheid Proprietary Limited	Cepheid	South Africa	Famous Grouse House, Kildrummy Office Park, Cnr. Witkoppen & Umhlanga Avenue, Paulshof 2191 South Africa
Cepheid Rus LLC	Cepheid	Russian Federation	Stanislavskogo Street 21, Bldg. 3 Bus Center Fabrika Stanis Moscow 109004 Russia
Cepheid S.r.l.	Cepheid	Italy	Attn: DR EMILIANO LEONE Via Emilia 26 20090 BUCCINASCO 20090 Milan Italy Italy
Cepheid Singapore Pte. Ltd.	Cepheid	Singapore	138 Market Street #05-06 CapitaGreen Singapore 048946 Singapore
Cepheid Switzerland GmbH	Cepheid	Switzerland	Zurcherstrasse 66 Postfach 124, Thalwil CH-8800 Switzerland

Cepheid UK Limited	Cepheid	United Kingdom	Oakley Court, Kingsmead Business Park Frederick Place High Wycombe HP11 1JU United Kingdom
InSilixa, Inc.	Cepheid	United States	1201 Orange Street Wilmington DE 19801 United States
Oncomark Limited	Cepheid	Ireland	Lismeehan, O' Callaghan Mills, Co. Clare Ireland
Devicor Medical Canada, Inc.	Devicor	Canada	250 HOWE STREET SUITE 1400 VANCOUVER BC V6C 3S7 Canada
Devicor Medical Device (Shanghai) Co., Ltd.	Devicor	China	Room 301, Building 2 No. 518, Fuquan North Road Changning District Shanghai China
Devicor Medical Europe GmbH	Devicor	Germany	Südportal 5, Nordport Towers, Erdgeschoss, 22848 Norderstedt Germany
Devicor Medical France SAS	Devicor	France	8 avenue de l'Île Saint Martin 92737 Nanterre Cedex France
Devicor Medical H.K. Limited	Devicor	Hong Kong	21st Floor Edinburgh Tower The Landmark 15 Queen's Road Central Hong Kong Hong Kong
Devicor Medical Holdings LLC	Devicor	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
Devicor Medical Italy S.r.l.	Devicor	Italy	Milan, Corso Magenta. 84, Italy Italy
Devicor Medical Japan K.K.	Devicor	Japan	DKK-TOA CORPORATION Annex Office Building 7th Floor, 29-9 Takadanobaba 1-chome Shinjuku-ku, Tokyo Japan
Devicor Medical Korea Limited	Devicor	Korea, Republic of	5F, 741(Eunsung Building, Chungdam-dong), Youngdong-daero, Gangnam-gu, Seoul, Korea, Republic of Korea, Republic of
Devicor Medical Poland sp. z o.o.	Devicor	Poland	31-121 Krakow, ul. Karmelicka 27 Poland

Devicor Medical Products de Mexico, S. de R.L. de C.V.	Devicor	Mexico	Avenida Angel Urruza 314 Del Valle Mexico 03100 Mexico
Devicor Medical Products Holdings, Inc.	Devicor	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
Devicor Medical Products Servicios, S. de R.L. de C.V.	Devicor	Mexico	Avenida Angel Urruza 314 Del Valle Distrito Federal 03100 Mexico
Devicor Medical Products, Inc.	Devicor	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
Devicor Medical Puerto Rico LLC	Devicor	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
Mammotome Medical Device Manufacturing (Shanghai), Co, Ltd.	Devicor	China	Floor 3B, 4B and % Building 1 T20 #258, Jin Znag Road Shanghai China
DH Sweden Holdings AB	Hemocue	Sweden	Box 1204, SE-262 23 Ängelholm Sweden
DHK Holding Company AB	Hemocue	Sweden	c/o Hemocue AB, P.O. Box 1204 262 23 Ängelholm 26223 Ängelholm Sweden
DHKAB Company AB	Hemocue	Sweden	c/o Hemocue AB, P.O. Box 1204 262 23 Ängelholm 26223 Ängelholm Sweden
FHAB Company Sweden AB	Hemocue	Sweden	c/o Hemocue AB, P.O. Box 1204 262 23 Ängelholm 26223 Ängelholm Sweden
HemoCue (Suzhou) Medical Technology Co., Ltd.	Hemocue	China	Room 207, No. 11 Modern Industrial Lane, No. 333 Xingpu Road, Suzhou Industrial Park Suzhou Free Trade Zone China
HemoCue AB	Hemocue	Sweden	Box 1204, SE-262 23 Ängelholm Sweden
HemoCue South Africa Pty Ltd	Hemocue	South Africa	P.O.Box 508, 2123, Pinegowrie, South Africa South Africa

RADCUE Holding AB	Hemocue	Sweden	Box 1204, SE-262 23 Ängelholm Sweden
IDT (Shanghai) Biotechnology Co., Ltd.	IDT	China	Shanghai One Site Room 602, Bldg 2, 518 North Fuquan Road IBP, Changning District Shanghai 200335 China
Integrated DNA Technologies Australia, Pty Ltd.	IDT	Australia	Level 19 181 William Street Melbourne VIC 3000 Australia
Integrated DNA Technologies Canada, Inc.	IDT	Canada	TD Tower, 77 Westmorland Street Suite 300 Fredericton NB E3B 6Z3 Canada
Integrated DNA Technologies Germany, GmbH	IDT	Germany	Terminal Straße Mitte 18 85356 München-Flughafen Germany
Integrated DNA Technologies K.K	IDT	Japan	Mita Kokusai Bldg. 24F 1-4-28 Mita Minato-ku Tokyo 108-0073 Japan
Integrated DNA Technologies Pte Ltd.	IDT	Singapore	7 Straits View, Marina One East Tower 018936 Singapore
Integrated DNA Technologies Spain, S.L.	IDT	Spain	Calle Prim 19, Bajo-Izquierda, 28004 Madrid Spain
Integrated DNA Technologies UK, Ltd	IDT	United Kingdom	19 Jessops Riverside 800 Brightside Lane Sheffield S9 2RX United Kingdom
Integrated DNA Technologies, B.V.	IDT	Belgium	Interleuvenlaan 12A B-3001 Leuven Belgium
Integrated DNA Technologies, Inc.	IDT	United States	c/o CT Corporation 1209 Orange Street Wilmington DE United States
Messenger Biotech, Inc.	IDT	Korea, Republic of	I-Teco 942-ho, 150, Jojeong- daero, Hanam-si, Gyeonggi- do 12930 Korea, Republic of

DH (Shanghai) Life Sciences Co., Ltd.	Life Sciences Group Staff	China	Shanghai One Site Room 602, Bldg 2, 518 North Fuquan Road IBP, Changning District Shanghai 200335 China
DH Life Sciences LLC	Life Sciences Group Staff	United States	1209 Orange Street Wilmington DE 19801 United States
Asymptote Limited	Cytiva	United Kingdom	Sovereign House, Vision Park Chivers Way Histon Cambridge CB24 9BZ United Kingdom
Biacore International AB	Cytiva	Sweden	Bjorkgatan 30 751 84 Uppsala Sweden
Biosafe America Inc.	Cytiva	United States	1999 Bryan Street Dallas TX 75201 United States
Biosafe Asia-Pacific Ltd.	Cytiva	Hong Kong	8/F., Lee Garden One, 33 Hysan Avenue, Causeway Bay Hong Kong
Biosafe Brasil Distribuidora Ltda	Cytiva	Brazil	Rua Luis Gama, 971, Bonfim Campinas 13070-717 Brazil
Biosafe do Brasil Participações e Representações Ltda	Cytiva	Brazil	Av. Magalhães de Castro, 4800, 11th floor, room 156, conj. 1 São Paulo 05676-120 Brazil
Biosafe Group SA	Cytiva	Switzerland	1-3 Route du Petit Eysins 1262 Eysins Switzerland
Biosafe Medical Device International Trading (Shanghai) Co Ltd.	Cytiva	China	Room 3804-H, No. 1601 West Nanjing Road Jingan District Shanghai China
Biosafe SA	Cytiva	Switzerland	1-3 Route du Petit Eysins 1262 Eysins Switzerland
Cytiva Belgium BV	Cytiva	Belgium	Reugelstraat 2 3320 Hoegaarden Belgium
Cytiva BioProcess R&D AB	Cytiva	Sweden	Bjorkgatan 30 751 84 Uppsala Sweden

Cytiva Bioscience Holding Limited	Cytiva	United Kingdom	Amersham Place, Little Chalfont Buckinghamshire HP7 9NA United Kingdom
Cytiva Biotechnology (Guangzhou) Co., Ltd.	Cytiva	China	Room 633, No. 333 Jian She Road, Jiu Fo Sino-Singapore Guang Guangzhou China
Cytiva Biotechnology (Hang Zhou) Co., Ltd.	Cytiva	China	1568 East Chunjiang Road Tonglu Hangzhou 311500 China
Cytiva CTH Holding Limited	Cytiva	United Kingdom	Amersham Place, Little Chalfont Buckinghamshire HP7 9NA Buckinghamshire HP7 9NA United Kingdom
Cytiva Denmark APS	Cytiva	Denmark	Akandevvej 21 DK-2700 Bronshoj Denmark
Cytiva do Brasil Comércio e Serviços para Biotecnologia Ltda.	Cytiva	Brazil	Av. Magalhaes de Castro, 4.800 Continental Tower Sao Paulo 305676-120 Brazil
Cytiva Europe GmbH	Cytiva	Germany	Munzinger Straße 5 79111 Freiburg Germany
Cytiva Finland Oy	Cytiva	Finland	Biolinja 12 Turku FI-20750 Finland
Cytiva Ireland Limited	Cytiva	Ireland	Crosshaven Road, Connemara Business Park Carrigaline Industrial Estate Carrigaline Cork Ireland
Cytiva Netherlands BV	Cytiva	Netherlands	Nijverheidsweg 1 1671 GC Medemblik Netherlands
Cytiva RUS LLC	Cytiva	Russian Federation	Stanislavsky str., 21 Building 3, Premise I Office 57 109004 Moscow Tagansky municipal district Russian Federation
Cytiva Sweden AB	Cytiva	Sweden	Bjorkgatan 30 751 84 Uppsala Sweden

Cytiva Sweden Holding AB	Cytiva	Sweden	Bjorkgatan 30 751 84 Uppsala Sweden
Cytiva Testa Center AB	Cytiva	Sweden	Bjorkgatan 30 753 23 Uppsala Sweden
Cytiva US LLC	Cytiva	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
Global Life Sciences IP Holdco LLC	Cytiva	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
Global Life Sciences Solutions Australia Pty Ltd	Cytiva	Australia	495 Blackburn Road Mount Waverly VIC 3149 Australia
Global Life Sciences Solutions Austria GmbH & Co KG	Cytiva	Austria	Kremplstrasse 5 4061 A Pasching Austria
Global Life Sciences Solutions Austria Verwaltungs GmbH	Cytiva	Austria	Kremplstraße 5, 4061 Pasching Austria
Global Life Sciences Solutions Canada ULC	Cytiva	Canada	PwC - Vancouver 250 Howe Street Suite 1400 Vancouver BC V6C3S7 Canada
Global Life Sciences Solutions Germany GmbH	Cytiva	Germany	Hahnestrasse 3 DE-37586 Dassel Germany
Global Life Sciences Solutions Korea Limited	Cytiva	Korea, Republic of	202, R&D Bldg., No. 2, 9, Songdomirae-ro Yeonsu-gu Incheon Korea, Republic of
Global Life Sciences Solutions Manufacturing UK Ltd	Cytiva	United Kingdom	19 Jessops Riverside 800 Brightside Lane Sheffield S9 2RX United Kingdom
Global Life Sciences Solutions New Zealand	Cytiva	New Zealand	Buddle Findlay, level 18, PwC Tower, 188 Quay Street Auckland 1010 New Zealand
Global Life Sciences Solutions Operations UK Ltd	Cytiva	United Kingdom	19 Jessops Riverside 800 Brightside Lane Sheffield S9 2RX United Kingdom

Global Life Sciences Solutions Poland Sp. z o.o.	Cytiva	Poland	Al. Jerozolimskie 181A 02-222 Mazowieckie Warsaw Poland
Global Life Sciences Solutions Singapore Pte. Ltd.	Cytiva	Singapore	25 Tuas South Street 1 Singapore 638034 Singapore
Global Life Sciences Solutions USA LLC	Cytiva	United States	1209 Orange Street Wilmington DE 19801 United States
Global Life Sciences Technologies (Shanghai) Co., Ltd.	Cytiva	China	Room A-738, 739, 188 Yesheng Road Pilot Free Trade Zone Shanghai China
Global Life Sciences Technologies Japan K.K.	Cytiva	Japan	3-25-1 Hyakunincho Sanken Building, Shinjuku-ku Tokyo 169-0073 Japan
Global Life Sciences Technologies Sweden Holdings AB	Cytiva	Sweden	Bjorkgatan 30 751 84 Uppsala Sweden
GLS Sweden Holdings AB	Cytiva	Sweden	Bjorkgatan 30 751 84 Uppsala Sweden
HyClone Laboratories, LLC	Cytiva	United States	C T Corporation System 1108 East South Union Avenue Midvale UT 84047 United States
Hyclone Life Sciences Solutions India Private Limited	Cytiva	India	No. 4 Kadugodi, Industrial Area, Whitefield Karnataka Bangalore 560067 India
Life Sciences IP Holdings Corporation	Cytiva	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
Puridify Ltd.	Cytiva	United Kingdom	Stevenage BioScience Catalyst Gunnels Wood Road Stevenage Hertfordshire SG1 2FX United Kingdom
Vanrx Pharmsystems Inc.	Cytiva	Canada	PwC - Vancouver 250 Howe Street Suite 1400 Vancouver BC V6C3S7 Canada

Bonna-Agela India Pvt. Ltd.	Phenomenex	India	2nd Floor, G-212 Sector 63 Noida 201301 India
BONNA-AGELA TECHNOLOGIES INC.	Phenomenex	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
FFM Holdings Pvt. Ltd.	Phenomenex	Mauritius	c/o Kross Border Corporate Corporate Services limited St Lous Business Centre, CNR Desroches & St Lous Streets, Port Louis, Mauritius India
Orange Holding Corp.	Phenomenex	United States	818 West 7th Street Los Angeles CA 90017 United States
Phenomenex ApS	Phenomenex	Denmark	Birketofte 13, Hareskov, 3500 Vaerloese Denmark
Phenomenex Australia PTY Ltd.	Phenomenex	Australia	Unit 35 2-6 Chaplin Drive Lane Cove NSW 2066 Australia
Phenomenex B.V.	Phenomenex	Netherlands	Maarssebroeksedijk 13-D, 3542DL Utrecht Netherlands
Phenomenex Distribution Centre GmbH	Phenomenex	Germany	Zeppelinstrasse 5, 63741 Aschaffenburg Germany
Phenomenex Espana, S.L.U.	Phenomenex	Spain	Valgrande 6, Edificio Thanworth II, 2, 1 B 28108 Alcobendas Spain
Phenomenex Helvetia GmbH	Phenomenex	Switzerland	Schaeferweg 16 4057 Basel Switzerland
Phenomenex India Holdings LLC	Phenomenex	United States	818 West 7th Street Los Angeles CA 90017 United States
Phenomenex India Pvt. Ltd.	Phenomenex	India	2-91/14/11 to 14/B, Laxmi Cyber City, B-Block Kondapur,, Hyderabad- 500084, Andhra Pradesh India
Phenomenex Ltd.	Phenomenex	United Kingdom	Melville House Queens Avenue, Hurdsfield Industrial Estate Macclesfield, Cheshire, SK10 2BN United Kingdom
Phenomenex N.Z. Ltd.	Phenomenex	New Zealand	Suite 5, Bld G, 4 Orbit Drive, Rosedale, North Shore City 0757 New Zealand
Phenomenex S.r.l.	Phenomenex	Italy	Milan, Via Montenapoleone nr. 8. Italy

Phenomenex SAS Ltd.	Phenomenex	France	60 Route De Sartrouville Parc Des Grillons Batiment 3 782390 Le Pecq France
Phenomenex, Inc.	Phenomenex	United States	818 West 7th Street Los Angeles CA 90017 United States
Phenova, Inc.	Phenomenex	United States	818 West 7th Street Los Angeles CA 90017 United States
Tianjin Bojieke Technologies Co. Ltd.	Phenomenex	China	No. 179, South Avenue, West District Tianjin Economy and Technology Development Zone China
Tianjin Bonna-Agela Technologies Co., Ltd.	Phenomenex	China	No. 179, South Avenue, West District Tianjin Economy and Technology Development Zone China
Hach DHR India Private Limited	OTT	India	D-128-129, First Floor, Okhla Industrial Area Phase-I, New Delhi, Delhi, India 110020 India
Lufft China Limited	OTT	Hong Kong	21st Floor, Edinburgh Tower, The Landmark 15 Queen's Road Central Hong Kong
OTT HydroMet B.V.	OTT	Netherlands	Delftechpark 36 2628 XH Delft Netherlands
OTT Hydromet Corp	OTT	United States	CT Corporation System 4701 Cox Road, Suite 301 Glen Allen VA 23060 United States
OTT HydroMet Fellbach GmbH	OTT	Germany	Gutenbergstr. 20, 70736 Fellbach Germany
OTT Hydromet GmbH	OTT	Germany	Ludwigstrasse 16 87437 Kempten Germany
OTT HydroMet Limited	OTT	United Kingdom	19 Jessops Riverside 800 Brightside Lane Sheffield S9 2RX United Kingdom
Ott Hydromet Pte. Ltd.	OTT	Singapore	10 Ubi Crescent #06-55 Lobby C Ubi Techpark 408564 Singapore

OTT HydroMet SARL	OTT	France	Immeuble le CLAMAR – Bât B 240 rue René Descartes 13799 Aix-en-Provence France
Alaska Blocker Corp.	Aldevron	United States	1209 Orange Street Wilmington DE 19801 United States
Aldevron Madison, LLC	Aldevron	United States	c/o CT Corporation System 120 W Sweet Ave Bismarck ND 58504 United States
Aldevron, L.L.C.	Aldevron	United States	c/o CT Corporation System 120 W Sweet Ave Bismarck ND 58504
Bison Holdco, LLC	Aldevron	United States	1209 Orange Street Wilmington DE 19801 United States
Nature Technology Corporation	Aldevron	United States	1201 Orange Street Wilmington DE 19801 United States
DH (Shanghai) Diagnostics Co., Ltd.	Diagnostics Group Staff	China	Shanghai One Site Room 602, Bldg 2, 518 North Fuquan Road IBP, Changning District Shanghai 200335 China
DH Diagnostics LLC	Diagnostics Group Staff	United States	1209 Orange Street Wilmington DE 19801 United States
DH Product Identification LLC	PID Group Staff	United States	1209 Orange Street Wilmington DE 19801 United States
AC INTERMEDIATE COMPANY	Corporate Entity	United States	C T Corporation System 4400 Easton Commons Way Suite 125 Columbus OH 43219 United States
ACME-Cleveland LLC	Corporate Entity	United States	C T Corporation System 4400 Easton Commons Way Suite 125 Columbus OH 43219 United States
Danaher Corporation	Corporate Entity	United States	CSC 251 Little Falls Drive Wilmington DE 19808 United States
Danaher Digital LLC	Corporate Entity	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
Danaher Finance Company	Corporate Entity	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States

Danaher Finance Company, LLC	Corporate Entity	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
Danaher Foundation	Corporate Entity	United States	C T Corporation System 208 South LaSalle Street, Suite 814 Chicago IL 60604 United States
Danaher Innovation Center LLC	Corporate Entity	United States	1209 Orange Street Wilmington DE 19801 United States
Danaher S&T Innovation Group LLC	Corporate Entity	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
DH Cooperative, Inc.	Corporate Entity	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
DH Holdings Corp.	Corporate Entity	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
FJ 900, Inc.	Corporate Entity	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
Industrial Fasteners, Inc.	Corporate Entity	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
Jennings Land Co.	Corporate Entity	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
Joslyn Electronic Systems Company, LLC	Corporate Entity	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
Joslyn Holding Company, LLC	Corporate Entity	United States	1209 Orange Street Corporation Trust Center Wilmington DE 19801 United States
Joslyn Manufacturing Company, LLC	Corporate Entity	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
KIIC Investment Company	Corporate Entity	Mauritius	Sanne House Bank Street Twentyeight Cybercity Ebene 72201 Mauritius

Premier Global Search, Inc.	Corporate Entity	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
Voluntary Employee Benefit Association	Corporate Entity	United States	2200 Pennsylvania Ave., NW, Suite 800, Washington, DC 20037
Alpha Lux Holdings S.à.r.l.	Corporate Tax Entity	Luxembourg	1B, Heienhaff, L-1736 Senningerberg Luxembourg
ARGO (DH) CANADA HOLDING ULC	Corporate Tax Entity	Canada	250 HOWE STREET SUITE 1400 VANCOUVER BC V6C 3S7 Canada
BCII Lux S.à.r.l.	Corporate Tax Entity	Luxembourg	1B Heienhaff, L-1736 Senningerberg Grand-Duchy of Luxembourg Luxembourg
BIAB Finance Corp.	Corporate Tax Entity	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
Danaher Australia Finance Pty Ltd	Corporate Tax Entity	Australia	495 Blackburn Road Mount Waverly VIC 3149 Australia
Danaher Australia Holding Pty Ltd	Corporate Tax Entity	Australia	495 Blackburn Road Mount Waverly VIC 3149 Australia
Danaher Canada Partners Inc.	Corporate Tax Entity	Canada	PwC Law LLP 18 York Street, Suite 2500-C Toronto ON M5J 0B2 Canada
Danaher China Holdings Limited	Corporate Tax Entity	Hong Kong	21st Floor Edinburgh Tower The Landmark 15 Queen's Road Central Hong Kong Hong Kong
Danaher Earnings Protection Insurance Company	Corporate Tax Entity	United States	Paul Frank + Collins P.C. One Church Street P.O. Box 1307 Burlington VT 05402-1307 United States
Danaher European Finance Company EHF	Corporate Tax Entity	Iceland	c/o Deloitte & Touche Smáratorgi 3 201 Kópavogur Iceland
Danaher European Finance S.a.r.l.	Corporate Tax Entity	Luxembourg	1B Heienhaff, L-1736 Senningerberg Grand-Duchy of Luxembourg Luxembourg
Danaher Finance Company Limited	Corporate Tax Entity	Cayman Islands	c/o International Corporation Services Ltd. PO Box 472 H George Town KY1-1106 Cayman Islands

Danaher General Partner Company Inc.	Corporate Tax Entity	Canada	PwC Law LLP 18 York Street, Suite 2600 Toronto ON M5J 0B2 Canada
Danaher HK Holdings Limited	Corporate Tax Entity	Hong Kong	21st Floor Edinburgh Tower The Landmark 15 Queen's Road Central Hong Kong Hong Kong
Danaher Iceland Finance Company EHF	Corporate Tax Entity	Iceland	c/o Deloitte & Touche Storhofdi 23 110 Reykjavik Iceland
Danaher Insurance Company	Corporate Tax Entity	United States	C T Corporation System 400 Cornerstone Drive, Suite 240 Williston VT 05495 United States
Danaher Luxembourg Finance S.à.r.l.	Corporate Tax Entity	Luxembourg	22, rue de l'Industrie, L-8399 Windhof, Grandy Duchy of Luxembourg Luxembourg
Danaher Participações Ltda	Corporate Tax Entity	Brazil	Rua Júlio González, 132 14th floor, conjunto 81 Memorial Office Building, Barra Funda Sao Paulo 01156-060 Brazil
DH Business Services (Shanghai) Co., Ltd.	Corporate Tax Entity	China	Address: 8 Floor Min Run Mansion, No.1388 Yishan Road, Xuhui China
DH Business Services LLC	Corporate Tax Entity	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
DH China Limited	Corporate Tax Entity	Cayman Islands	c/o International Corporation Services Ltd. PO Box 472 H George Town KY1-1106 Cayman Islands
DH Egypt	Corporate Tax Entity	Egypt	M002-M003 Al Raya building Plot 133, BAnking Sector New Cairo Cairo Egypt
DH Europe Finance II Sarl	Corporate Tax Entity	Luxembourg	1B, Heienhaff, L-1736 Senningerberg Luxembourg
DH Europe Finance S.à.r.l.	Corporate Tax Entity	Luxembourg	1B Heienhaff, L-1736 Senningerberg Grand-Duchy of Luxembourg Luxembourg

DH Finance Corp.	Corporate Tax Entity	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
DH France Sarl	Corporate Tax Entity	France	SILIC 703, 16 Avenue du Québec, Villebon sur Yvette 91961 Courtaboeuf Cedex France
DH Holding Vietnam Limited	Corporate Tax Entity	Vietnam	Room 037 (037D) Centec Business Center, 4th Floor, 72-74 Nguyen Thi Minh Vietnam
DH Holdings Italia S.r.l.	Corporate Tax Entity	Italy	Corso Venezia 61 20121 Milano Italy
DH Japan Finance S.à.r.l.	Corporate Tax Entity	Luxembourg	1B Heienhaff, L-1736 Senningerberg Grand-Duchy of Luxembourg Luxembourg
DH Japan Holding G.K.	Corporate Tax Entity	Japan	6-5-1 Nishishinjyuku, Shinjyuku-ku Tokyo 163-1325 Japan
DH Japan LLC	Corporate Tax Entity	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
DH Luxemburg S.a.r.l.	Corporate Tax Entity	Luxembourg	1B Heienhaff, L-1736 Senningerberg Grand-Duchy of Luxembourg Luxembourg
DH Omega Holdings Inc	Corporate Tax Entity	United States	1201 Orange Street Wilmington DE 19801 United States
DH Ontario Finance Limited	Corporate Tax Entity	Canada	PwC Canada - Ontario 18 York Street Suite 2600 Toronto ON M5J 0B2 Canada
DH Rus Services LLC	Corporate Tax Entity	Russian Federation	Malookhtinsky 64 bldg 3 Saint-Petersburg Russia 195112 Russian Federation
DH SB Holdings, LLC	Corporate Tax Entity	United States	1209 Orange Street Corporation Trust Center Wilmington DE 19801 United States
DH Surety LLC	Corporate Tax Entity	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
DH Switzerland Finance S.à.r.l.	Corporate Tax Entity	Luxembourg	1B Heienhaff, L-1736 Senningerberg Grand-Duchy of Luxembourg Luxembourg

DH UK Partners	Corporate Tax Entity	United Kingdom	c/o Shepherd & Wedderburn Condor House 10 St. Paul's Churchyard London England EC4M 8AL United Kingdom
DHCDAN Australia Pty Ltd	Corporate Tax Entity	Australia	495 Blackburn Road Mount Waverly VIC 3149 Australia
DHCY Limited	Corporate Tax Entity	Cayman Islands	c/o International Corporation Services Ltd. PO Box 472 H George Town KY1-1106 Cayman Islands
DJ Acquisition Pty Ltd	Corporate Tax Entity	Australia	495 Blackburn Road Mount Waverly VIC 3149 Australia
DJAPL Investments LLC	Corporate Tax Entity	United States	c/o The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
DLFSA Lux Holdings S.à.r.l.	Corporate Tax Entity	Luxembourg	1B Heienhaff, L-1736 Senningerberg Grand-Duchy of Luxembourg Luxembourg
DMG Partners	Corporate Tax Entity	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
DMG Plastics, Inc.	Corporate Tax Entity	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
Dolan-Jenner Industries, Inc.	Corporate Tax Entity	United States	C T Corporation System 155 Federal Street, Suite 700 Boston MA 02110 United States
Easco Cayman Holdings Ltd.	Corporate Tax Entity	Cayman Islands	c/o International Corporation Services Ltd. PO Box 472 H George Town KY1-1106 Cayman Islands
Easco Lux S.à.r.l.	Corporate Tax Entity	Luxembourg	1B Heienhaff, L-1736 Senningerberg Grand-Duchy of Luxembourg Luxembourg
FCL Hong Kong Holding Limited	Corporate Tax Entity	Hong Kong	21st Floor Edinburgh Tower The Landmark 15 Queen's Road Central Hong Kong Hong Kong

FFCL Lux Sàrl	Corporate Tax Entity	Luxembourg	1B Heienhaff, L-1736 Senningerberg Grand-Duchy of Luxembourg Luxembourg
Joslyn LLC	Corporate Tax Entity	United States	Corporation Trust Center 1209 Orange Street Wilmington DE 19801 United States
KAIC Investment Company	Corporate Tax Entity	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
Launchchange, Inc.	Corporate Tax Entity	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
Omega Lux Finance S.à.r.l.	Corporate Tax Entity	Luxembourg	1B Heienhaff, L-1736 Senningerberg Grand-Duchy of Luxembourg Luxembourg
PN UK Finance Limited	Corporate Tax Entity	Cayman Islands	c/o International Corporation Services Ltd. PO Box 472 H George Town KY1-1106 Cayman Islands
PN UK Holding Limited	Corporate Tax Entity	Cayman Islands	c/o International Corporation Services Ltd. PO Box 472 H George Town KY1-1106 Cayman Islands
QHC Holding Company	Corporate Tax Entity	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
Radiometer Finance Corp.	Corporate Tax Entity	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
Radiometer Medical Sales Pty Limited	Corporate Tax Entity	Australia	1st floor, 96 Ricketts Rd Mount Waverly VIC 3149 Australia
Seneca Realty Inc.	Corporate Tax Entity	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
Utica Holding Company	Corporate Tax Entity	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
VSE Holdings Pty Ltd	Corporate Tax Entity	Australia	495 Blackburn Road Mount Waverly VIC 3149 Australia

X-Rite Holdings S.à.r.l.	Corporate Tax Entity	Luxembourg	1B Heienhaff, L-1736 Senningerberg Grand-Duchy of Luxembourg Luxembourg
Aegean Partner LLC	Legacy Tools Division	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
Armstrong Tools, LLC	Legacy Tools Division	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
Cispus Hong Kong Holding Limited 塞司香港控股有限公司	Legacy Tools Division	Hong Kong	21st Floor Edinburgh Tower The Landmark 15 Queen's Road Central Hong Kong Hong Kong
Danaher (Shanghai) Management Co. Ltd.	Legacy Tools Division	China	Room 2723, 450 Fu Shan Road Shanghai pilot free trade zone , Shanghai China
Danaher Hong Kong Finance Limited 丹納赫香港財務有限公司	Legacy Tools Division	Hong Kong	21st Floor Edinburgh Tower The Landmark 15 Queen's Road Central Hong Kong Hong Kong
Danaher Hong Kong Limited	Legacy Tools Division	Hong Kong	21st Floor Edinburgh Tower The Landmark 15 Queen's Road Central Hong Kong Hong Kong
Danaher Innovation Centre (Shanghai) Co., Ltd. 丹纳赫创思(上海)管理有限公司	Legacy Tools Division	China	Building C, No.888 Huang Hu North No. 2 Road, Lingang Free Trade Zone Shanghai China
DCI Consolidated Industries, Inc.	Legacy Tools Division	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
Easco Hand Tools, Inc.	Legacy Tools Division	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
H-K Connecticut Holding Company, LLC	Legacy Tools Division	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
JS Technology, LLC	Legacy Tools Division	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States

Quality Wire Processing, Inc.	Legacy Tools Division	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
Terryville Realty Inc.	Legacy Tools Division	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
Yukon Hong Kong Holding Limited	Legacy Tools Division	Hong Kong	21st Floor Edinburgh Tower The Landmark 15 Queen's Road Central Hong Kong Hong Kong
DHR Holding India Pvt. Ltd.	Corporate - India	India	3rd floor, B Wing, Art Guild House Phoenix Marketcity, Kurla West Mumbai 400070 India
PT DHH Trading Indonesia	Legacy PSA	Indonesia	10F World Trade Centre, WTC 5, 10th Floor Jl. Jend. Sudirman Kav. 29 Jakarta 12920 Indonesia
Life Sciences Holdings France SAS	Legacy PSA	France	16 avenue du Quebec 703-91961 Courtaboeuf Cedex France
WA 10 SAS	Legacy PSA	France	16 avenue du Quebec 703-91961 Courtaboeuf Cedex France
WA 11 SAS	Legacy PSA	France	16 avenue du Quebec 703-91961 Courtaboeuf Cedex France
WA Mistral SAS	Legacy PSA	France	SILIC 703, 16 Avenue du Québec, Villebon sur Yvette 91961 Courtaboeuf Cedex France
WA9 SAS	Legacy PSA	France	SILIC 703, 16 Avenue du Québec, Villebon sur Yvette 91961 Courtaboeuf Cedex France
TCIL Ireland Finance Ltd	Legacy Comms Enterprise Group	Ireland	Lismeehan, O' Callaghan Mills, Co. Clare Ireland
Aristotle Brazil Finance Limited	Fortive Legacy	United Kingdom	19 Jessops Riverside 800 Brightside Lane Sheffield S9 2RX United Kingdom

Danaher Malta Limited	Fortive Legacy	Malta	Level 1 LM Complex Brewery Street Mrieħel Birkirkara BKR 3000 Malta
DH Alpha Limited	Fortive Legacy	United Kingdom	19 Jessops Riverside 800 Brightside Lane Sheffield S9 2RX United Kingdom
DH Holdings US II Inc.	Fortive Legacy	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
DH Omega Limited	Fortive Legacy	United Kingdom	19 Jessops Riverside 800 Brightside Lane Sheffield S9 2RX United Kingdom
DH TM US LLC	Fortive Legacy	United States	1209 Orange Street Wilmington DE 19801 United States
DH UK Finance Limited	Fortive Legacy	United Kingdom	19 Jessops Riverside 800 Brightside Lane Sheffield S9 2RX United Kingdom
DHR MENA FZ-LLC	Fortive Legacy	United Arab Emirates	301, BUILDING NO 34, AL ZAHRAWI DUBAI HEALTHCARE CITY PO BOX 71569 DUBAI United Arab Emirates
FHC Holding Corp.	Fortive Legacy	United States	1209 Orange Street Wilmington DE 19801 United States
FIC International Corporation	Fortive Legacy	United States	The Corporation Trust Company 1209 Orange Street Wilmington DE 19801 United States
FSIL Finance Ltd	Fortive Legacy	Ireland	Lismeehan, O' Callaghan Mills, Co. Clare Ireland
GCC Canada Corporation	Fortive Legacy	Canada	Stewart McKelvey Lawyers 900-1959 Upper Water Street HALIFAX NS Canada B3J 3N2 Halifax NS B3J 3N2 Canada
GHC Holding Corporation	Fortive Legacy	United States	1209 Orange Street Wilmington DE 19801 United States

GHKH Hong Kong Limited	Fortive Legacy	Hong Kong	21st Floor Edinburgh Tower The Landmark 15 Queen's Road Central Hong Kong Hong Kong
Gilzoni Holdings Italia S.R.L.	Fortive Legacy	Italy	SCAFATI (SA) VIA PASSANTI 174 CAP 84018 Italy
Gilzoni Limited	Fortive Legacy	United Kingdom	19 Jessops Riverside 800 Brightside Lane Sheffield S9 2RX United Kingdom
GL UK Holdings Limited	Fortive Legacy	United Kingdom	19 Jessops Riverside 800 Brightside Lane Sheffield S9 2RX United Kingdom
Launchchange Finance Limited	Fortive Legacy	United Kingdom	19 Jessops Riverside 800 Brightside Lane Sheffield S9 2RX United Kingdom
Launchchange Holding Company	Fortive Legacy	United Kingdom	19 Jessops Riverside 800 Brightside Lane Sheffield S9 2RX United Kingdom
Launchchange Holdings Limited	Fortive Legacy	United Kingdom	19 Jessops Riverside 800 Brightside Lane Sheffield S9 2RX United Kingdom
Launchchange Instrumentation Limited	Fortive Legacy	United Kingdom	19 Jessops Riverside 800 Brightside Lane Sheffield S9 2RX United Kingdom
Launchchange Limited	Fortive Legacy	United Kingdom	19 Jessops Riverside 800 Brightside Lane Sheffield S9 2RX United Kingdom
Launchchange Operations Limited	Fortive Legacy	United Kingdom	19 Jessops Riverside 800 Brightside Lane Sheffield S9 2RX United Kingdom
Launchchange UK Holdco Limited	Fortive Legacy	United Kingdom	19 Jessops Riverside 800 Brightside Lane Sheffield S9 2RX United Kingdom

LIL Holdco UK Limited	Fortive Legacy	United Kingdom	19 Jessops Riverside 800 Brightside Lane Sheffield S9 2RX United Kingdom
Namco Controls Corporation	Fortive Legacy	United States	C T Corporation System 4400 Easton Commons Way Suite 125 Columbus OH 43219 United States
PallBio Holdings Limited	Fortive Legacy	United Kingdom	19 Jessops Riverside 800 Brightside Lane Sheffield S9 2RX United Kingdom
Radiometer Holdco Limited	Fortive Legacy	United Kingdom	19 Jessops Riverside 800 Brightside Lane Sheffield S9 2RX United Kingdom
Sarbel UK Holding Company	Fortive Legacy	United Kingdom	19 Jessops Riverside 800 Brightside Lane Sheffield S9 2RX United Kingdom
Stampede Acquisition Limited	Fortive Legacy	United Kingdom	19 Jessops Riverside 800 Brightside Lane Sheffield S9 2RX United Kingdom
VJ HoldCo UK Limited	Fortive Legacy	United Kingdom	19 Jessops Riverside 800 Brightside Lane Sheffield S9 2RX United Kingdom
VJH Holding Co.	Fortive Legacy	United States	1209 Orange Street Wilmington DE 19801 United States
VR Finance Company	Fortive Legacy	United Kingdom	19 Jessops Riverside 800 Brightside Lane Sheffield S9 2RX United Kingdom
VRC Holding Corporation	Fortive Legacy	United States	1209 Orange Street Wilmington DE 19801 United States