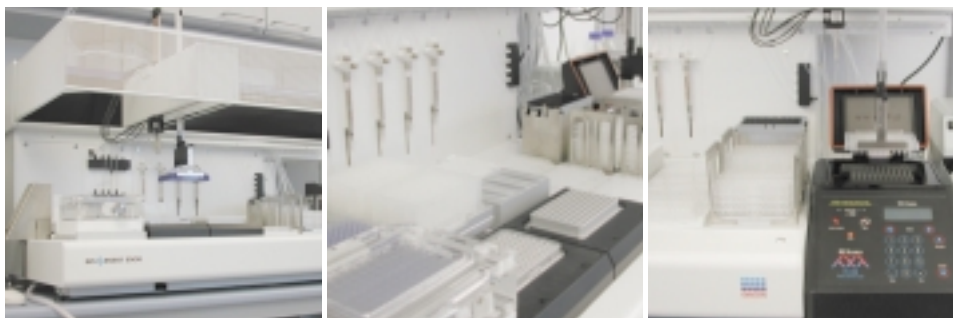




QIAGEN® BioRobot® Extended-Arm Systems



The QIAGEN® BioRobot® extended-arm systems allow the robotic handling system to move labware and accessories between the BioRobot 3000 worktable and external instruments (Figure 1, next page). The configuration (i.e., arm extending from either the left or right side of the workstation) is chosen based on the external instruments to be accessed (Table 1). Protocols that include processing or analysis performed by other instruments, such as thermal cycling or spectrophotometry, can be fully automated with this system.

The QIAsoft™ Operating System allows data to be exchanged between the workstation and external instruments for sample and data tracking. For example, dilutions of DNA samples prepared on the workstation can be automatically transferred to a spectrophotometer. After analysis, the spectrophotometric data is imported into a QIAsoft table and dilution calculations performed. The workstation then uses the calculated data to either dilute the original samples to uniform concentrations or pipette a specified quantity of DNA for reaction setup.

Table 1. Examples of instruments accessed by extended-arm workstations

Instrument	Example
Spectrophotometer	SPECTRAmax® from Molecular Devices
Thermal cycler	PTC200 from MJ Research
Heating device	Heating system from QIAGEN Instruments

This table indicates instruments with which extended-arm workstations have already been used. Other instruments can be also used with the workstation. Contact QIAGEN for the most current information.



Area Accessible by Extended-Arm BioRobot Configurations

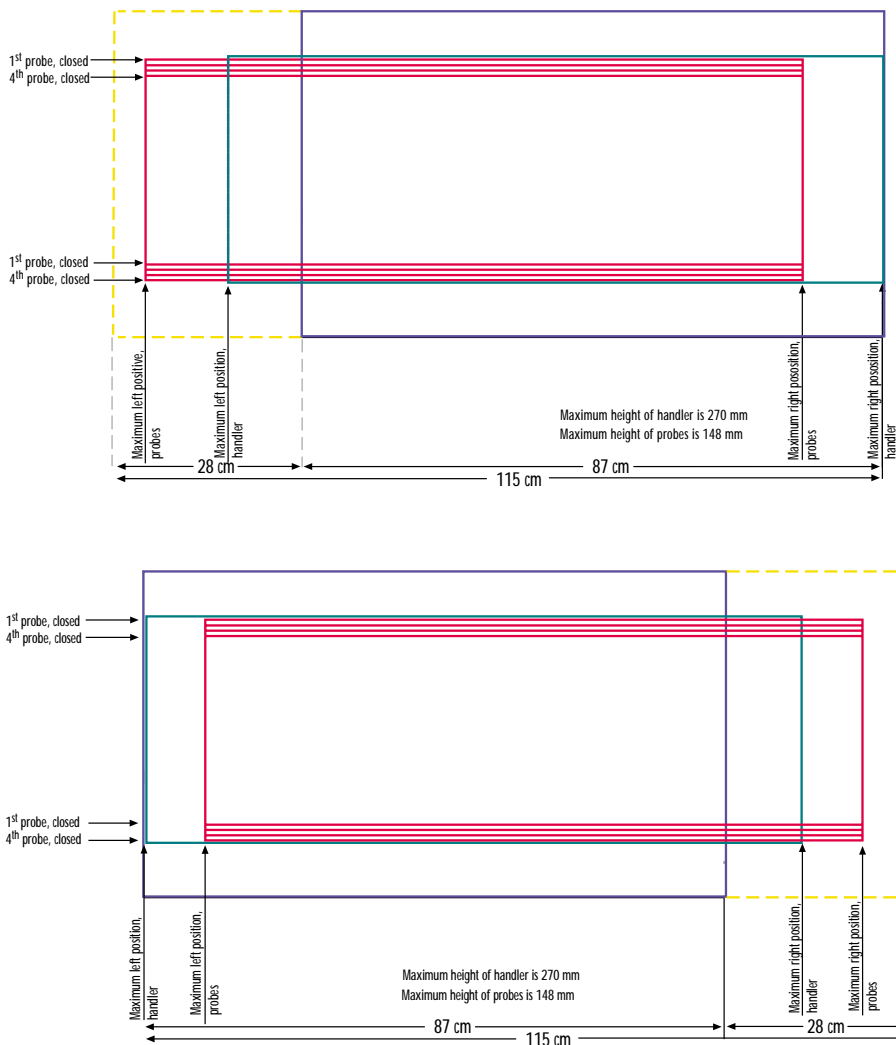


Figure 1. This diagram show the areas, both on the worktable and outside of the instrument, which can be accessed by the left and right extended-arm configurations. **Blue lines**: worktable surface; **red lines**: area accessible by the pipetting probes; **green lines**: area accessible by the handler.

Contact QIAGEN for more information about the products mentioned in this Technical Note.

QIAGEN robotic workstations are not available in all countries; please inquire. Patented or patent-pending technology and/or registered or registration-pending trademarks of QIAGEN: QIAGEN®, QIASoft™, BioRobot®. SPECTRAMax is a registered trademark of Molecular Diagnostics Corporation. Registered names, trademarks, etc. used in this document, even when not specifically marked as such, are not to be considered unprotected by law.

© 2001 QIAGEN, all rights reserved.

www.qiagen.com

QIAGEN:

- Australia** Tel. 03-9489-3666 Fax 03-9489-3888
- Canada** Tel. 800-572-9613 Fax 800-713-5951
- France** Tel. 01-60-920-930 Fax 01-60-920-925
- Germany** Tel. 02103-29-12400 Fax 02103-29-22000
- Italy** Tel. 02-33430411 Fax 02-33430426
- Japan** Tel. 03-5547-0811 Fax 03-5547-0818
- Switzerland** Tel. 061-319-30-31 Fax 061-319-30-33
- UK** Tel. 01293-422-999 Fax 01293-422-922
- USA** Tel. 800-426-8157 Fax 800-718-2056

Distributors:

- Argentina** Tecnolab S.A. (011) 4555 0010
- Austria/Hungary/Slovenia** R. u. P. MARGARITELLA Austria (01) 889 18 19
- Belgium/Luxemburg** Westburg b.v. 0800-1-9815
- Brazil** Uniscience do Brasil 011 3622 2320
- China** Gene Company Limited (852)2896-6283
- Cyprus** Scientronics Ltd (02) 765 416
- Czech Republic** BIO-CONSULT spol. s.r.o. (02)4447 1239
- Denmark** Merck EuroLab A/S 43 86 87 88
- Egypt** Clinilab 525 7212
- Finland** Merck EuroLab Oy (09)804 551
- Greece** BioAnalytica S.A. (01)640 03 18
- India** Genetix (011)542 1714 or (011)515 9346
- Israel** Westburg (Israel) Ltd. 07. 66 50 814 or 1-800 20 22 20
- Korea** LRS Laboratories, Inc. (02) 924-86 97
- Malaysia** Research Biolabs Sdn. Bhd. (03)7312099
- Mexico** Quimica Valaner S.A. de C.V. (5) 525 57 25
- The Netherlands** Westburg b.v. (033)4950094
- New Zealand** Biolab Scientific Ltd. (09) 980 6700 or 0800 933 966
- Norway** Merck EuroLab AS 22 90 00 00
- Poland** Syngen Biotech Sp.z.o.o. (071) 351 41 06 or 0601 70 60 07
- Portugal** IZASA PORTUGAL, LDA (1)424 73 64
- Singapore** Research Biolabs Pte Ltd. 445 7927
- Slovak Republic** BIO-CONSULT Slovakia spol. s.r.o. (07) 50221 336
- South Africa** Southern Cross Biotechnology (Pty) Ltd (021) 671 5166
- Spain** IZASA, S.A. (93) 902.20.30.90
- Sweden** Merck EuroLab AB (08) 621 34 00
- Taiwan** TAIGEN Bioscience Corporation (02) 2880 2913
- Thailand** Theera Trading Co. Ltd. (02) 412-5672
- In other countries** contact: QIAGEN, Germany

