

Reagents and Supplies

Get high-value clones faster and advance your BioPharma research with complete reagents and supplies



Introduction

Unlock excellence in biopharmaceutical research.

Welcome to the forefront of biopharmaceutical research and development with Molecular Devices. This guide showcases an exceptional range of reagents meticulously designed to streamline and enhance your biopharmaceutical workflow.

In the dynamic realm of biopharmaceutical production, the right clone selection can significantly elevate results. With instruments, reagents, and supplies for microbial and mammalian cell users, you can retain the integrity of growing cell colonies, pave the way for automated identification and extraction of clonal cell lines, and accelerate the journey from concept to market.

Embark on a path of excellence and innovation, where every product is a step towards reshaping the future of biopharmaceuticals.

For more information, visit:

www.moleculardevices.com/products/biologics/clone-screening.

Table of contents

01	Cell biology reagents	3
	CloneMedia CHO Growth A	3
	Recombinant CloneDetect	3
	CloneDetect	4
	CloneMatrix and CloneXL	4
	Cleaning solutions	5
	Calibeads	5
02	QPix® Chroma Filters	5
	QPix Chroma Filters	5
03	Adjustable petri dish and microplate holders	6
04	Bioassay trays	6
05	Microplates	7
	Low profile microplates	7
	384-well high sample recovery plate	7

06	Deep well plates and lids	8
	Deep well plates	8
	Lids	8
07	Adhesive and heat seals	9
	Breathe seal	9
80	Replicators	9
	QReps polypropylene replicators	9
09	Accessories	10
10	ClonePix Welcome Kits	11
	ClonePix Welcome Kit A	11
	ClonePix Welcome Kit B	11

O1 Cell biology reagents

Cell culture media, supplements, and antibody-based reagents for use in cell line development of hybridomas, CHO and HEK cells

CloneMedia CHO Growth A

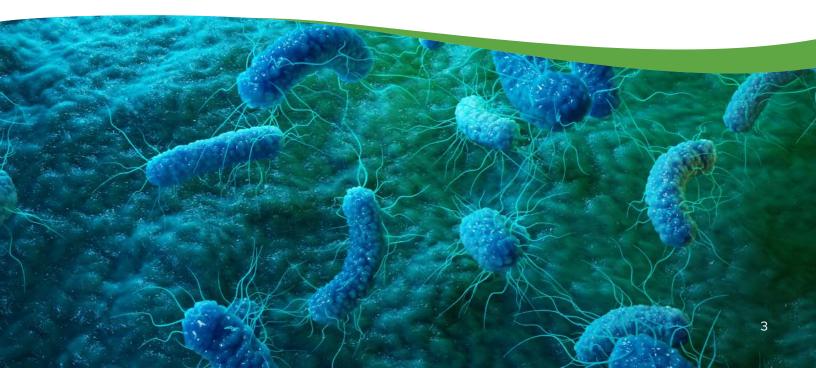
The CloneMedia CHO Growth A semi-solid media portfolio provides a fast, simple, and comprehensive solution for CHO cell line development. Each batch of K8495 is subjected to rigorous safety assessment, with mycoplasma, virus and endotoxin full certification. Optimized for recombinant protein production. Free of animal components.

Product name and description	Quantity/pack size	Part number	Compatibility
CloneMedia CHO Growth A with L-Gln	6 × 90 mL	K8800	
CloneMedia CHO Growth A with L-Gln	1 × 90 mL	K8810	ClanaDiv
CloneMedia CHO Growth A without L-Gln	6 × 90 mL	K8830	ClonePix
CloneMedia CHO Growth A without L-Gln	1 × 90 mL	K8840	

Recombinant CloneDetect

Newest detection agent free of animal origin for visualizing secreted antibodies.

Product name and description	Quantity/pack size	Part number	Compatibility
CloneDetect Human IgG (Fc) Specific, Recombinant, Fluorescein	10,000U / 1 mL	K8495	ClonePix, CloneSelect Imager FL



O1 Cell biology reagents

CloneDetect

Detection agents for visualizing secreted antibodies. All products are animal-derived, sterile, and azide-free.

Product name and description	Quantity/pack size	Part number	Compatibility
Human			
CloneDetect Human IgG (H+L) Specific, Fluorescein	10,000U / 1 mL	K8400	
CloneDetect Human IgG (H+L) Specific, Fluorescein, BSA Free	10,000U / 1 mL	K8202	ClonePix, CloneSelect Imager FL
CloneDetect Human IgG (Fc) Specific, Fluorescein	10,000U / 1 mL	K8205	_ IIIIagei i L
Mouse			
CloneDetect Mouse IgG (H+L) Specific, Fluorescein	10,000U / 1 mL	K8420	ClonePix,
CloneDetect Mouse IgG (Fc) Specific, Fluorescein	10,000U / 1 mL	K8425	CloneSelect Imager FL

CloneMatrix

A 2.5x concentrate that serves as a matrix to produce semi-solid media from 2x concentrated liquid media provided by user. Proven for hybridoma, myeloma, CHO, HEK and mESC. Supplied ready to dilute in 100 mL bottles for convenience.

Product name and description	Quantity/pack size	Part number	Compatibility
CloneMatrix, makes 6 × 100 mL final media volume	6 × 40 mL	K8500	
CloneMatrix, makes 100 mL final media volume	1 × 40 mL	K8510	
CloneMatrix-CHO: optimized for CHO cells, makes 6 × 100 mL final media volume; contains 6 vials of 2ml CloneXL and 6 bottles of 40ml CloneMatrix	6 × 40 mL	K8520	ClonePix
CloneXL reagent: 50x concentrated animal component-free cell culture supplement converts CloneMatrix to CloneMatrix-CHO	5 × 2 mL	K8525	
CloneMatrix-CHO: optimized for CHO cells, makes 100 mL final media volume; contains 1 vial of 2ml CloneXL and 1 bottle of 40ml CloneMatrix	1 × 40 mL	K8530	

O1 Cell biology reagents

Cleaning solutions

Ideally suited to clean instruments and accessories. Removes proteins and biological materials from most metals, glass, plastic, rubber, fabric surface, and is safe to use with ultrasonic cleaners/washers. Environmentally friendly — contains no phosphates, chlorofluorocarbons (CFCs) or borates.

Product name and description	Quantity/pack size	Part number	Compatibility
aQu Clean 50x concentrated pin cleaning solution, anionic and non-ionic detergent based	1L	K2505	ClonePix, QPix 400 Series

Calibeads

A must-have tool to calibrate, optimize, and keep your ClonePix instrument with optimal performance. This ready-to-use product is made from the mixture of fluorescent microspheres with 4 different wavelengths in a semi-solid substrate format. Microspheres fluoresce at four distinct wavelengths corresponding to the CFP, FITC / GFP, Rhodamine / PE and Cy5™ filter sets of ClonePix systems.

Product name and description	Quantity/pack size	Part number	Compatibility
Calibeads: Fluorescent calibration beads for ClonePix. 200 µm diameter. Pan wavelength. Provided ready-to-plate (in 100 mL of semi-solid substrate, ca. 15-25 beads / mL)	100 mL	K8150	ClonePix

QPix Chroma Filter

QPix Chroma Filters

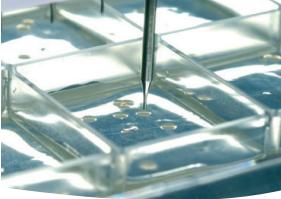
Thin film optical filters for selection of bacterial colonies with multiple ranges of color intensities. Interchangeable filters last up to 10 uses and improve recognition of colored colonies for accurate picking of desired clones. Compatible with QPix Software Version 2.0 and higher.



Product name and description	Quantity/pack size	Part number	Compatibility
QPix Chroma Filter; Thin film optical filter for Blue/White colony selection	1 filter	X1101	
QPix Chroma Filter; Thin film optical filters for Blue/White colony selection	25 filters	X1102	QPix 400 Series
QPix Chroma Colorimetric Colony Selection Software Kit; Blue/White Colony Selection Software Kit for QPix 400 Series. Includes software license and 2 QPix™ Chroma Filters (X1101)	1 license, 25 filters	X1103	







O3 Petri dishes and plate holders

Adjustable petri dish and microplate holders

Spring-loaded holders automatically adjust to the perimeter of the plates to ensure secure fit of the plastic ware with a greater size variations. Black finish enhances optical imaging and reduces light scattering. Compatible with QPix Software Version 2.0 and higher.

Product name and description	Plate diameter	Part number	Compatibility
1-Way Adjustable Petri Dish Holder	138–141 mm	X9401	
2-Way Adjustable Petri Dish Holder	Any ANSI/SLAS 1-2004 microplates	X9402	QPix 400 Series
4-Way Adjustable Petri Dish Holder	97–100.5 mm	X9403	
5-Way Adjustable Petri Dish Holder	87– 90.5 mm	X9404	

04 Bioassay trays

QTrays

Ideal for large scale culture and screening of bacterial colonies, regional picking and high throughput protein expression. The vented lid aids removal of condensation. Can be supplied with a 48-well divider to provide a cost effective alternative to petri dishes.

Product name and description	Quantity/pack size	Part number	Compatibility
Vented QTray with cover, polystyrene, 240 \times 240 \times 20 mm, sterile, individually wrapped	20 per case	X6023	ODiv 400 Sorios
Vented Qtray with cover & 48-Well Divider, polystyrene, 242 x 240 x 20 mm, sterile, individually wrapped	20 per case	X6029	QPix 400 Series

05 Microplates

A comprehensive range of 96- and 384-well microplates in a variety of profiles and materials.



Low profile microplates

Low profile design maximizes robotic plate stacker volume and storage. Dimensions: $128 \times 86 \times 10 \text{ mm}$.

Product name and description	Quantity/pack size	Part number	Compatibility
384-Well Low Profile Microplate, round well, flat bottom, sterile, with cover, working volume 50 μ L, packaged in sleeves of 10	160 per case	X6001	
384-Well Low Profile Microplate, round well, V-bottom, sterile, with cover. Individually wrapped	160 per case	X6004	- QPix 400 Series
96-Well Low Profile Microplate, polystyrene, round well, flat bottom, sterile, with cover, working volume 180 μ L, packaged in sleeves of 10	160 per case	X6011	QPIX 400 Selles
384-Well Low Profile, Large Volume Microplate, round well, flat bottom, sterile, with cover, working volume 60 μ L, packaged in sleeves of 10	160 per case	X7001	

384-well high sample recovery plate

Microplates for applications that require high sample recovery rates. Made from polypropylene with conical well profiles to maximize recovery and allow working with sample volumes as low as 5 μ L. The low profile increases robotic stack capacity and storage. Dimensions: 128 x 86 x 10 mm. Working volume: 5–65 μ L.

Product name and description	Quantity/pack size	Part number	Compatibility
384-Well High Sample Recovery Plate, polypropylene, without cover, working volume 5-65 $\mu L, packaged$ in sleeves of 10	200 per case	X7020	ODiv 400 Sovice
384-Well High Sample Recovery Plate Plate, polypropylene, with cover, working volume 5-65 μL , packaged in sleeves of 10	200 per case	X7022	QPix 400 Series

06 Deep well plates and lids

Deep well plates

Polypropylene deep well plates with elevated well walls for easier sealing. Conform to SBS standards. Non-pyrogenic, free of detectable DNase, RNase, human DNA. Ideal for storing sample materials, cell stock or cultivating bacterial, phage, yeast, suspension mammalian cell culture etc.

Product name and description	Quantity/pack size	Part number	Compatibility
96 Deep Square Well Plate, round bottom, working volume 2000 μL sterile, packed in sleeves of 6	24 plates per case	S1205	
384 Deep Square Well Plate, round bottom, working volume 400 $\mu\text{L},$ packed in sleeves of 6	24 plates per case	S1207	QPix 400 Series
24 Deep Square Well Plate, conical bottom, working volume 10 mL sterile, individually wrapped	25 plates per case	X7060	

Lids

QLid is a rigid polystyrene, close fitting lid. It is automation compatible, prevents cross-contamination among samples and reduces excessive evaporation.

Product name and description	Quantity/pack size	Part number	Compatibility
QLid, sterile, packed in sleeves of 10: fits S1205, S1207 and X7020	240 lids per case	X5101	QPix 400 Series



O7 Adhesive and heat seals

Breathe Seal

Porous film for sealing culture plates where air and gas exchange is necessary. Fits any ANSI/SLAS 1-2004 microplates. The 114 μ m thick film is easily pierced by pipette tips. Recommended for temperatures between -20 to +80°C.



Product name and description	Quantity/pack size	Part number
Breathe Seal for cell and tissue culture, non-sterile	100 seals per pack	E1004
Breathe Seal for cell and tissue culture, sterile, packed in tamper proof bags of 25	50 seals per pack	E1005

08 Replicators

Sterile QReps polypropylene replicators

Increase processing speed by eliminating flaming/autoclaving, washing and using these pre-sterilized chemically resistant replicators.



Product name and description	Quantity/pack size	Part number	Compatibility
QRep 384 Pin Replicators: polypropylene, 10 mm length (low profile plates), 1 mm diameter	2 × 64	X5050	
QRep 96 Pin Replicators: polypropylene, 10 mm length (low profile plates), 1 mm diameter	2 × 64	X5051	
QRep 96 Pin Replicators: polypropylene, 10 mm length (low profile plates), 2 mm diameter	2 × 64	X5052	QPix 400 Series
QRep 384 Pin Replicators: polypropylene, 15 mm length (standard profile plates), 1 mm diameter	2 × 56	X5053	
QRep 96 Pin Replicators: polypropylene, 15 mm length (standard profile plates), 1 mm diameter	2 × 56	X5054	



09 Accessories

Product name and description	Quantity/pack size	Part number	Compatibility
130 μL ART pipette tips, filtered, sterile	96 tips per pack (10 packs total 960 tips)	T0130	QPix 400 Series
DISPENKIT all-in-one pack for single-cell dispensing workflow, sterile	4 sensing tips, 4 cell strainers, 4 reservoirs, 5 vials of optimized medium, 1 box of 96-wide bore tips	5306092	DispenCell Single-Cell Dispenser
DISPENSTAR pack for single-cell dispensing workflow, sterile	4 sensing tips, 4 cell strainers, 4 reservoirs, 5 vials of optimized media with beads, 1 box of 96-wide bore tips	5307040	

10 ClonePix Welcome Kits

ClonePix Welcome Kit A (for CHO Users)

Are you already a CHO Cell user? The CHO Growth A media portfolio provides a fast, simple, and comprehensive solution for CHO cell line development. Combined with the CloneDetect detection agent, CloneSelect® Imager, and ClonePix® 2 System, researchers can more efficiently develop new protein-producing CHO cell lines, thus accelerating time-to-market.

- · Optimized for high recombinant protein production
- · Ready-to-use culture media and supplements are free of animal components
- · Enhanced optical quality for colony imaging
- One-step fast screening and selection of high-value clones

Kit (PN# K9000) includes:

- 1 bottle of CHO growth A without L-Glutamine
- · 1 vial of Human CloneDetect
- 1 vial of Mouse CloneDetect

- 1 vial of CloneXL 50x Concentrated Animal Component-Free Supplement
- · 1 bottle of ClonePix Calibeads



CloneMatrix[™] is an easy-to-use product that converts existing mammalian cell biology media to semi-solid form. Semi-solid media maintains the positional integrity of growing cell colonies, thus allowing the automated identification and picking of clonal cell lines.

- The use of semi-solid media combined with the unique fluorescent screening and picking technology of the ClonePix System will revolutionize clonal cell line development in biopharmaceuticals.
- CloneMatrix is a unique methylcellulose concentrate that allows users to add their own media formulation to produce an optimized semisolid media reagent.
- CloneMatrix is conveniently supplied as 40 mL in a 100 mL bottle, reducing the risk of
 contamination as there is no requirement to aliquot the concentrate. Supplied as a 2.5x
 concentrate, simply add optimized media (2x concentration) selection and supplements
 as required.
- CloneXL is a 50x-concentrated cell growth supplement proven to improve mammalian cell survival and support CHO, HEK293, stem cell clone screening, and single-cell cloning applications. It contains amino acids and thymidine family members with patented formulation.

Kit (PN# K9001) includes:

- 1 bottle of CloneMatrix
- 1 vial of Human CloneDetect
- 1 vial of Mouse CloneDetect

- 1 vial of CloneXL 50x Concentrated Animal Component-Free Supplement
- 1 bottle of ClonePix Calibeads









ClonePix® systems
Screening and selection
of mammalian clones



QPix® systemsSets the industry standard in accurate colony picking



A simple single-cell dispenser for proof of monoclonality

DispenCell[™] Single-Cell

Dispenser



CloneSelect® Imager
Objective, quantitative
assessment of cell growth



Next-generation fluorescent imaging solution for the assurance of monoclonality

CloneSelect® Imager FL

Contact Us

Phone: +1.800.635.5577

Web: www.moleculardevices.com

Email: info@moldev.com

Check our website for a current

listing of worldwide distributors.

Regional Offices

 USA and Canada
 +1.800.635.5577
 Taiwan/Hong Kong
 +886.2.2656.7585

 United Kingdom
 +44.118.944.8000
 Japan
 +81.3.6362.9109

 Europe*
 00800.665.32860
 South Korea
 +82.2.3471.9531

 China
 +86.4008203586
 India
 +91.73.8661.1198

rldwide distributors.
*Austria, Belgium, Denmark, Finland, France, Germany, Iceland, Ireland, Italy, Luxembourg, Netherlands, Portugal, Spain, Sweden, Switzerland and United Kingdom