



Assay Kits, Reagents, and Consumables

Simplifying life science and biopharmaceutical
research with robust, easy-to-use solutions

Unlock the full potential of your research with our optimized and validated assay kits, reagents and consumables



ImageXpress
High-Content
Imaging Systems



**ImageXpress
Pico**
Automated Cell
Imaging Systems



**SpectraMax®
i3/i3x**
Multi-Mode
Microplate
Readers



**SpectraMax®
Mini**
Multi-Mode
Microplate
Reader



**SpectraMax®
Paradigm®**
Multi-Mode
Microplate
Reader



**SpectraMax®
iD3/iD5**
Multi-Mode
Microplate
Readers



**SpectraMax®
M Series**
Multi-Mode
Microplate
Readers

Cell Viability Assay Kits

- Apoptosis
- Cardiotoxicity
- Cell Integrity
- Live Cell
- Live/Dead

Cell Signaling Assay Kits

- CatchPoint cAMP
- CatchPoint cGMP

Plate Washer Accessories

- AquaMax Sterilant Kit

dsDNA Quantitation Assay Kits

- Quant HiRange
- Quant Nano
- Quant Pico

DNA/RNA/Protein Quantitation Microplate

- SpectraDrop Starter
- SpectraDrop HTS
- SpectraDrop Refill

3D Ready Patient-Derived Organoids

- Healthy Controls
- Colorectal Cancer
- Breast Cancer

Enzyme Assay Kits

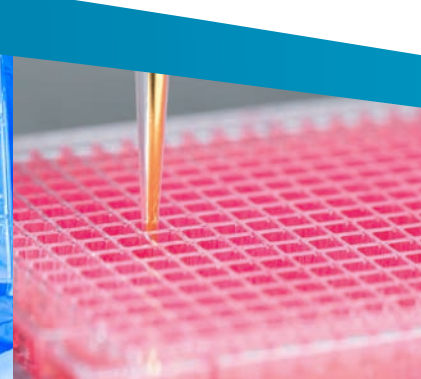
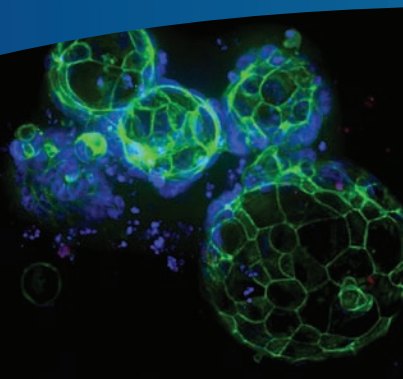
- IMAP Kinase
- IMAP Phosphatase
- IMAP Phosphodiesterase

Protein Expression Assay Kits

- CatchPoint SimpleStep ELISA
- Western Blot

Reporter Gene Assay Kit

- Dual Luciferase
- Luciferase



Tailored specifically for your instruments, our solutions come with pre-configured software protocols and streamlined analysis tools, allowing you to swiftly transform samples into insightful answers. Enhance your drug discovery, development, and bioassay processes with unmatched efficiency and reliability.



FlexStation® 3
Multi-Mode
Microplate Reader



FLIPR® Penta
High-Throughput
Cellular Screening
System



DispenseCell™
Single-Cell
Dispenser



CloneSelect®
Series
Cell Imaging
Systems



ClonePix®
Series
Mammalian Colony
Pickers



QPix®
400 Series
Microbial Colony
Pickers

Calcium Mobilization Assay Kits

Calcium
Fura-2 QBT Calcium

Ion Channels/Transporter Assay Kits

Membrane Potential
Neurotransmitter
Transporter Uptake
Potassium
QBT Fatty Acid Uptake

Pipette Tips

FLEX Station Tips
FLIPR Tips

Accessories

DISPENSTAR
DISPENKIT

Calibration

Calibeads

Pin Cleaning

aQu Clean

CloneMedia Kits

CHO Growth A
CloneMatrix
CloneXL

CloneDetect Assay Kits

Animal-free human IgG
Human IgG
Mouse IgG

QPix Accessories

Heads, Pins, Springs
Petri Dish Holders
Chroma Filters
QReps Replicators
QTips
QTrays
QBlocks
Deep Well Plates
High Recovery Plates
Low Profile Plates
Breathe Seal
QLids



Learn more



Assay kits

Easy-to-use, robust assay kits for life science research, drug discovery and development, and bioassays. All assay kits are optimized for use on our instruments.

- Fluorescent and luminescent microplate reader assay kits
- High throughput screening assay kits for FLIPR instruments
- Unique western blot detection enabled via ScanLater system
- Assess cell viability at any time with homogenous fluorescent no-wash assay kits via ImageXpress systems



eBook: Mastering optimal plate reader assays

Master the use of plate reader assays in your drug discovery research with our comprehensive eBook. Gain expert insights, step-by-step guidance, and valuable tips to enhance your assay optimization process. Explore real-world case studies and access detailed instructions on using our assay kits, instruments, and software. Don't miss this opportunity to revolutionize your research methodology and accelerate groundbreaking discoveries.



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

This guide showcases an exceptional range of reagents meticulously designed to streamline and enhance your biopharmaceutical workflow. 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.

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

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

