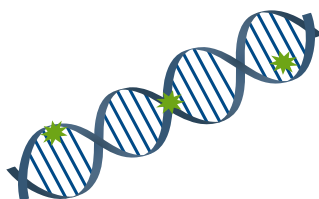


# Optimized assay kits for life science research and drug discovery



## SpectraMax Quant dsDNA Assay Kits

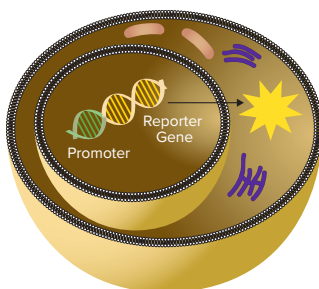
The SpectraMax® Quant™ dsDNA Assay Kits are designed for fluorescence-based dsDNA quantitation across a broad range of concentrations. Tailored to your different needs, these kits are configured and optimized for our SpectraMax® microplate readers with preconfigured protocols provided in SoftMax® Pro Software for streamlined data acquisition and analysis.



Ordering information			
Item	Size	Part #	US Price
SpectraMax® Quant™ AccuClear™ Nano dsDNA Assay Kit	Explorer Kit	R8356	\$100.00
	Bulk Kit	R8357	\$320.00
SpectraMax® Quant™ AccuBlue™ HiRange dsDNA Assay Kit	Explorer Kit	R8358	\$110.00
	Bulk Kit	R8359	\$330.00
SpectraMax® Quant™ AccuBlue™ Pico dsDNA Assay Kit	Explorer Kit	R8354	\$130.00
	Bulk Kit	R8355	\$390.00

## SpectraMax Glo Steady-Luc Reporter Assay Kit

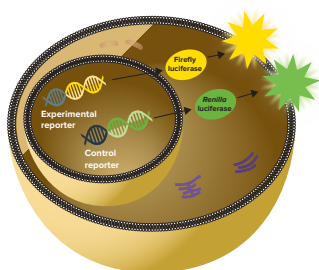
The SpectraMax® Glo Steady-Luc™ Reporter Assay Kit provides a highly sensitive assay for the quantitation of firefly luciferase expression in mammalian cells. This kit is optimized for our SpectraMax microplate readers with a preconfigured protocol provided in SoftMax Pro Software for streamlined data acquisition and analysis.



Ordering information			
Item	Size	Part #	US Price
SpectraMax Glo Steady-Luc Reporter Assay Kit	Explorer Kit	R8352	\$220.00
	Bulk Kit	R8353	\$650.00

## SpectraMax DuoLuc Reporter Assay Kit

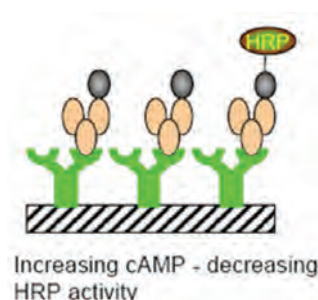
The SpectraMax® DuoLuc™ Reporter Assay Kit enables highly sensitive quantitation of both firefly and *Renilla* luciferases in mammalian cells. The assay kit is optimized for Molecular Devices SpectraMax® microplate readers with injectors and includes a preconfigured protocol in SoftMax Pro Software for simplified data acquisition and analysis.



Ordering information			
Item	Size	Part #	US Price
SpectraMax DuoLuc Reporter Assay Kit	Evaluation Kit	R8360	\$104.00
	Explorer Kit	R8361	\$256.00
	Bulk Kit	R8362	\$955.00
Passive Lysis Buffer	1 bottle	R8363	\$18.00

## CatchPoint cAMP and cGMP Fluorescent Assay Kits

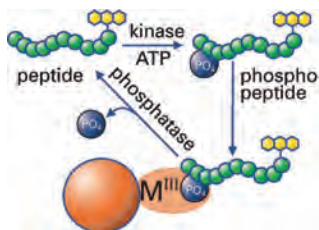
The CatchPoint® cAMP and cGMP Fluorescent Assay Kits' high-affinity reagents are optimized for sensitivity and precision in applications where cAMP and cGMP levels are low. A single wash step removes unbound material prior to the development step, so the assays are highly resistant to interference from colored or fluorescent test compounds.



Ordering information			
Item	Size	Part #	US Price
CatchPoint cAMP 96-well Fluorescent Assay Kit	Explorer Kit	R8088	\$1,949.00
	Bulk Kit	R8089	\$3,389.00
CatchPoint cAMP 384-well Fluorescent Assay Kit	Explorer Kit	R8044	\$2,159.00
	Bulk Kit	R8053	\$5,599.00
CatchPoint cGMP 96-well Fluorescent Assay Kit	Explorer Kit	R8074	\$1,949.00
	Bulk Kit	R8075	\$3,389.00
CatchPoint cGMP 384-well Fluorescent Assay Kit	Explorer Kit	R8065	\$2,159.00
	Bulk Kit	R8066	\$5,599.00

## IMAP Kinase, Phosphatase, and Phosphodiesterase Assays

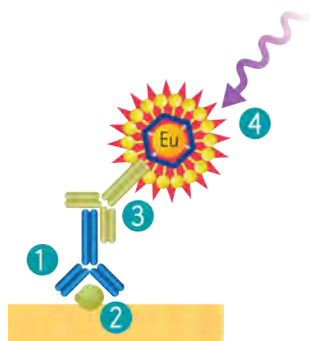
IMAP® technology provides a homogeneous assay for the assessment of kinase, phosphatase, and phosphodiesterase (PDE) activity. The assay affords a simple “mix-and-read” procedure utilizing free phosphate-binding nanoparticles directly reporting converted product, not enzyme reaction components or by-products.



Ordering information			
Item	Size	Part #	US Price
IMAP Demo Kit	Evaluation Kit	R8166	\$619.00
IMAP FP Kit with Progressive Binding System	Evaluation Kit	R8155	\$639.00
IMAP TR-FRET Kit with Progressive Binding System	Evaluation Kit	R8161	\$639.00
IMAP FP PDE Kit	Evaluation Kit	R8175	\$1,259.00
IMAP TR-FRET PDE Kit	Evaluation Kit	R8176	\$1,259.00

## ScanLater Western Blot Assay Kit

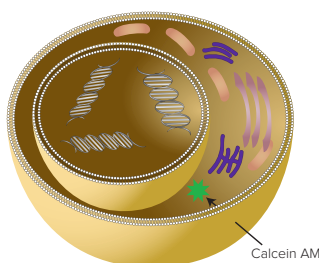
The ScanLater™ Western Blot Assay Kit enables the first western blot detection capability in a microplate reader. Utilizing this substrate-free assay, researchers can achieve equivalent sensitivity to traditional chemiluminescence and fluorescence-based assays, while consolidating western blot and ELISA applications onto a single reader.



Ordering information		
Item		
ScanLater Western Blot Assay Kit	Anti-Mouse Kit	Eu-Streptavidin Kit
	Anti-Rabbit Kit	Protein Ladder
	Anti-Goat Kit	Secondary Antibodies
	Anti-Rat Kit	
	<a href="#">See webpage for details</a>	

## EarlyTox Cell Viability Assay Kits

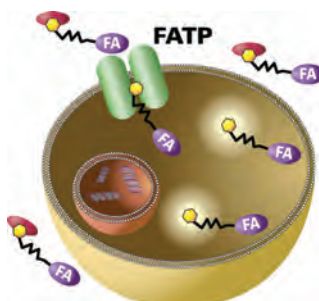
The EarlyTox™ Cell Viability Assay Kits are a family of fluorescence-based reagents for the assessment of cell viability, cell proliferation, and various apoptosis events using mammalian cells. Optimized for use with microplate readers, these assay kits employ a no-wash, homogeneous assay protocol that enables characterization of a full concentration-response profile of test compounds.



Ordering information			
Item	Size	Part #	US Price
EarlyTox™ Live/Dead Assay Kit	Explorer Kit	R8340	\$136.00
	Bulk Kit	R8341	\$293.00
EarlyTox™ Live Cell Assay Kit	Explorer Kit	R8342	\$115.00
	Bulk Kit	R8343	\$247.00
EarlyTox™ Glutathione Assay Kit	Explorer Kit	R8344	\$160.00
	Bulk Kit	R8345	\$345.00
EarlyTox™ Caspase-3/7 R110 Assay Kit	Explorer Kit	R8346	\$345.00
	Bulk Kit	R8347	\$748.00
EarlyTox™ Caspase-3/7-D NucView™ 488 Assay Kit	Explorer Kit	R8348	\$690.00
	Bulk Kit	R8349	\$2,392.00
EarlyTox™ Caspase-3/7 NucView™ 488 Assay Kit	Explorer Kit	R8350	\$690.00
	Bulk Kit	R8351	\$2,392.00

## QBT Fatty Acid Uptake Assay Kit

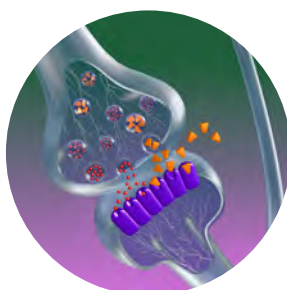
The QBT™ Fatty Acid Uptake Assay is a homogeneous assay amenable to high-throughput screening. The kits deliver pre-optimized, fluorescence-based formulations to expedite assay development and screening of fatty acid transporters.



Ordering information			
Item	Size	Part #	US Price
QBT Fatty Acid Uptake Assay Kit	Explorer Kit	R8132	\$1,059.00
	Bulk Kit	R8133	\$2,409.00

## Neurotransmitter Transporter Uptake Assay Kit

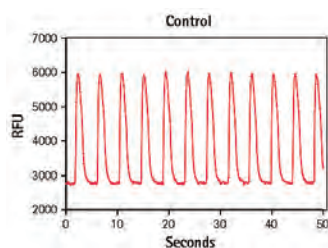
The ability to monitor serotonin, norepinephrine, and dopamine neurotransmitter uptake is key to a better understanding of diseases such as Alzheimer's and Parkinson's. With the Neurotransmitter Transporter Uptake Assay Kit, researchers now have a tool to study these three key neurotransmitters with a live-cell, fluorescent, plate reader-based assay.



Ordering information			
Item	Size	Part #	US Price
Neurotransmitter Transporter Uptake Assay Kit	Explorer Kit	R8173	\$609.00
	Bulk Kit	R8174	\$4,239.00

## EarlyTox Cardiotoxicity Kit

The EarlyTox™ Cardiotoxicity Kit provides a fast, simple, and reliable fluorescence-based method for identifying cardiotoxic compounds in a biorelevant assay. Using cultured cardiomyocytes and a kinetic plate reader, researchers can prioritize leads and direct medicinal chemistry efforts sooner, improving productivity and reducing costs associated with extensive safety testing downstream.

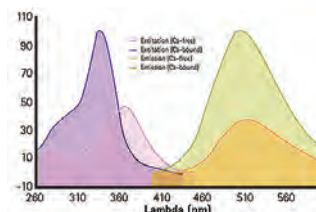


### Ordering information

Item	Size	Part #	US Price
EarlyTox Cardiotoxicity Kit	Explorer Kit	R8210	\$319.00
	Bulk Kit	R8211	\$999.00

## Fura-2 QBT Calcium Kit

The Fura-2 QBT™ Calcium Kit affords a simple, mix-and-read format that employs our proprietary masking technology with the industry-standard Fura-2 ratiometric calcium indicator. Streamline your current Fura-2 assay for calcium concentration determination, or for use with GFP-tagged proteins.

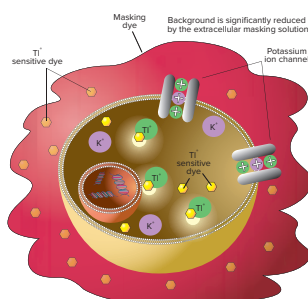


### Ordering information

Item	Size	Part #	US Price
Fura-2 QBT Calcium Kit	Explorer Kit	R8197	\$499.00
	Bulk Kit	R8198	\$3,239.00

## FLIPR Potassium Assay Kit

The FLIPR® Potassium Assay Kit measures functional activity of ligand- or voltage-gated potassium channels. The homogeneous, no-wash assay protocol also features a large signal window and high Z' values.



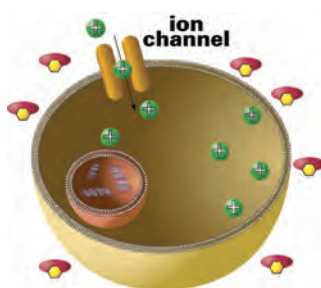
### Ordering information

Item	Size	Part #	US Price
FLIPR Potassium Assay Kit	Explorer Kit	R8222	\$809.00
	Bulk Kit	R8223	\$4,729.00



## FLIPR Membrane Potential Assay Kit

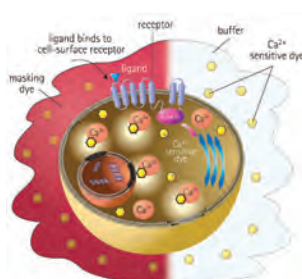
The FLIPR® Membrane Potential Assay Kit delivers homogeneous, fluorescence-based formulations for observation of real-time change of membrane potential associated with ion channel activation and ion transporter proteins.



Ordering information			
Item	Size	Part #	US Price
FLIPR Membrane Potential Red Assay Kit	Explorer Kit	R8126	\$609.00
	Bulk Kit	R8123	\$3,029.00
FLIPR Membrane Potential Blue Assay Kit	Explorer Kit	R8042	\$609.00
	Bulk Kit	R8034	\$3,029.00

## FLIPR Calcium Assay Kits

The FLIPR® Calcium Assay Kits are the most comprehensive calcium reagent portfolio on the market. These kits measure changes in intracellular calcium in response to GPCR or ion channel activities. They deliver pre-optimized, homogeneous, fluorescence-based formulations to expedite assay development and compound screening for drug discovery and basic research.



Ordering information			
Item	Size	Part #	US Price
FLIPR Calcium 6 Assay Kit	Explorer Kit	R8190	\$779.00
	Bulk Kit	R8191	\$3,929.00
FLIPR Calcium 6-QF Assay Kit	Explorer Kit	R8192	\$669.00
	Bulk Kit	R8193	\$3,679.00
FLIPR Calcium 5 Assay Kit	Explorer Kit	R8185	\$609.00
	Bulk Kit	R8186	\$3,629.00
FLIPR Calcium 4 Assay Kit	Explorer Kit	R8142	\$799.00
	Bulk Kit	R8141	\$3,929.00
FLIPR Calcium 3 Assay Kit	Explorer Kit	R8091	\$839.00
	Bulk Kit	R8090	\$4,239.00
FLIPR Calcium Assay Kit	Explorer Kit	R8041	\$619.00
	Bulk Kit	R8033	\$3,099.00

# Assay Kits and Instrument Compatibility

Select the images or text for detailed information.



**SpectraMax®  
iD3 Multi-Mode  
Microplate  
Reader**



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



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



**SpectraMax®  
L Microplate  
Reader**



**FlexStation® 3  
Multi-Mode  
Microplate  
Reader**



**FLIPR Tetra®  
System**

<b>SpectraMax Quant dsDNA Assay Kits</b>	✓	✓	✓		✓	
<b>SpectraMax Glo Steady-Luc Reporter Assay Kit</b>	✓	✓	✓	✓	✓	
<b>SpectraMax DuoLuc Reporter Assay Kit</b>	✓*	✓*		✓*		
<b>CatchPoint cAMP and cGMP Fluorescent Assay Kits</b>	✓	✓	✓		✓	
<b>IMAP Kinase, Phosphatase, and Phosphodiesterase Assays</b>	✓	✓	✓		✓	
<b>ScanLater Western Blot Assay Kit</b>		✓				
<b>EarlyTox Cell Viability Assay Kits</b>	✓	✓	✓		✓	
<b>QBT Fatty Acid Uptake Assay Kit</b>	✓	✓	✓		✓	✓
<b>Neurotransmitter Transporter Uptake Assay Kit</b>	✓	✓	✓		✓	✓
<b>EarlyTox Cardiotoxicity Kit</b>	✓	✓			✓	✓
<b>Fura-2 QBT Calcium Kit</b>					✓	✓
<b>FLIPR Potassium Assay Kit</b>					✓	✓
<b>FLIPR Membrane Potential Assay Kit</b>					✓	✓
<b>FLIPR Calcium Assay Kits</b>					✓	✓

\*Equipped with injectors



#### Contact Us

Phone: [800.635.5577](tel:800.635.5577)  
Web: [www.moleculardevices.com](http://www.moleculardevices.com)  
Email: [info@moldev.com](mailto:info@moldev.com)

Check our website for a current listing  
of worldwide distributors.