










# How to upgrade your SpectraMax® i3x Multi-Mode Microplate Reader with additional detection modes



Cartridge	Description	Part number	Specifications	Optimized sensitivity	Guaranteed sensitivity	Slots used
 Dual Auto Injector	<ul style="list-style-type: none"> <li>SpectraMax i3x Injector Cartridge with SmartInject™ Technology</li> <li>Expand your research capabilities to include flash-based applications, including dual luciferase and ATP assays.</li> </ul>	0200-7029	Wavelength range (LUM): Visible to 650 nm Dead volume: < 10 µL with Reverse Prime function	20 amol ATP ("flash" luminescence using Promega Enliten Assay System)	50 amol ATP (250 fM @ 0.2mL/well, "Flash" luminescence using Promega Enliten Assay System) 3 fMol ATP (15 pM @ 0.2mL/well, "Glow" luminescence using PerkinElmer ATPlite 1step Luminescence Assay System)	2
 ScanLater	<ul style="list-style-type: none"> <li>Western blot detection using ScanLater Western Blot Assay Kit</li> <li>TRF-based with 340/80 nm EX and 616/10 nm EM</li> </ul>	0200-7027	EX range: 340/80 nm EM range: 616/10 nm	High fg levels of Streptavidin	High fg levels of Streptavidin	2
 AlphaScreen	<ul style="list-style-type: none"> <li>AlphaScreen and AlphaLISA detection using 1 W 680 nm EX laser diode and a 570 nm (100) EM filter</li> <li>Pick best speed, sensitivity, and price for your needs</li> <li>Guaranteed sensitivity: &lt;100 amol of biotinylated-LCK-P peptide, 25 µL/well in a 384-well plate, using AlphaScreen Phosphotyrosine (PT66) Assay Kit</li> </ul>	0200-7017	Alpha 384 STD 96- and 384-well plates		< 100 amol (384-well)	1
		0200-7018	Alpha 384 HTS 96- and 384-well plates		< 100 amol (384-well)	1
		0200-7019	Alpha 1536 HTS 96, 384, and 1536-well plates		< 100 amol (384-well)	1
 HTRF	<ul style="list-style-type: none"> <li>Cisbio HTRF detection with optimized Xenon light source and 616, 665 nm EM filters</li> </ul>	0200-7011	6- to 1536-well plates	Exceeds Cisbio certification requirements		2
 TRF	<ul style="list-style-type: none"> <li>LED light source EX: 370/80 nm EM: 616/10 nm, 642/10 nm</li> <li>Suitable for assays using europium and samarium chelate labels</li> </ul>	0200-7008	6- to 1536-well plates	96-well: 0.03 pM 384-well: 0.03 pM 1536-well: 0.125 pM *Specifications listed are for europium	96-well: 0.1 pM 384-well: 0.1 pM 1536-well: 0.375 pM	1
 FP	<ul style="list-style-type: none"> <li>Fluorescence Polarization detection for fluorescein- or rhodamine-like labels, using specific LED and EX/EM filters</li> <li>6- to 1536-well plates</li> </ul>	0200-7009	Fluorescein FP EX 485 nm, EM 535P and 535S nm	96-well: 1.0 mP 384-well: 1.5 mP 1536-well: 2.0 mP	96-well: 3 mP 384-well: 4 mP 1536-well: 6 mP	1
		0200-7010	Rhodamine FP EX 535 nm, EM 595P and 595S nm			

Cartridge	Description	Part number	Specifications	Optimized sensitivity	Guaranteed sensitivity	Slots used
 Fluorescence Intensity (FI)	<ul style="list-style-type: none"> <li>Fluorescence Intensity (FI) detection for coumarin-, fluorescein-, or rhodamine-like labels, and FRET</li> <li>Ultra-high-power LEDs for the excitation of various fluorescent labels</li> </ul>	0200-7002	EX: 360/35 nm EM1: 465/35 nm EM2: 535/25 nm	384-well plate (75 µL): 10 fmol/well 1536-well plate (8 µL): 3 fmol/well		1
 Glow Luminescence (LUM) Detection Cartridge	<ul style="list-style-type: none"> <li>Glow Luminescence (LUM) Detection</li> </ul>	0200-7012	96-, 384- and 1536-well plates Wavelength range: Visible to 650 nm	96-well (200 µL) 3 pM 384-well (50 µL) 7 pM 1536-well (8 µL) 20 pM *Specifications listed are for ATP	96-well (200 µL) 15 pM 384-well (50 µL) 30 pM 1536-well (8 µL) 60 pM	1
		0200-7015	384-well plates Wavelength range: Visible to 650 nm	96-well (200 µL) 2 pM 384-well (50 µL) 3 pM	96-well (200 µL) 10 pM 384-well (50 µL) 15 pM	1
		0200-7014	96-well plates Wavelength range: Visible to 650 nm	96-well (200 µL) 2 pM	96-well (200 µL) 8 pM	1
Dual Color Luminescence (LUM) (BRET2) Detection Cartridge	<ul style="list-style-type: none"> <li>Dual Color Luminescence (LUM) (BRET2™)</li> <li>Detection Cartridge that isolates each well to provide optimal performance for chemiluminescence applications</li> </ul>	0200-7016	EM1: 410/35 nm EM2: 515/35 nm			1
Transmitted Light (TL) Detection Cartridge (for use with the SpectraMax MiniMax 300 Imaging Cytometer)	<ul style="list-style-type: none"> <li>Transmitted Light (TL) Detection provides white LED illumination for transmitted-light (brightfield) imaging with the SpectraMax MiniMax 300 Imaging Cytometer</li> </ul>	5022671	Proprietary solid-state illumination, white LED			1
 Custom Solutions	<ul style="list-style-type: none"> <li>Custom cartridges are available and designed to meet your specific research needs</li> </ul>	Ask your local sales representative for more information				

## Contact Us

Phone: +1.800.635.5577  
 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.

## 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

