










# 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<br>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)<br>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<br>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<br/>EX: 370/80 nm<br/>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<br>384-well: 0.03 pM<br>1536-well: 0.125 pM<br>*Specifications listed are for europium | 96-well: 0.1 pM<br>384-well: 0.1 pM<br>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<br>EX 485 nm,<br>EM 535P and 535S nm   | 96-well: 1.0 mP<br>384-well: 1.5 mP<br>1536-well: 2.0 mP  | 96-well: 3 mP<br>384-well: 4 mP<br>1536-well: 6 mP   | 1          |
|  |  | 0200-7010   | Rhodamine FP<br>EX 535 nm,<br>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<br>EM1: 465/35 nm<br>EM2: 535/25 nm                     | 384-well plate (75 µL): 10 fmol/well<br>1536-well plate (8 µL): 3 fmol/well                                    |  | 1          |
|   |   | 0200-7003  | EX: 485/20 nm<br>EM1: 535/20 nm<br>EM2: 595/25 nm                     | 384-well plate (75 µL): 0.10 fmol/well<br>1536-well plate (8 µL): 0.03 fmol/well                               |  | 1          |
|   |   | 0200-7004  | EX: 535/25 nm<br>EM1: 595/35 nm<br>EM2: LP 655 nm                     | 384-well plate (75 µL): 0.15 fmol/well<br>1536-well plate (8 µL): 0.06 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<br>Wavelength range: Visible to 650 nm | 96-well (200 µL) 3 pM<br>384-well (50 µL) 7 pM<br>1536-well (8 µL) 20 pM<br>*Specifications listed are for ATP | 96-well (200 µL) 15 pM<br>384-well (50 µL) 30 pM<br>1536-well (8 µL) 60 pM | 1          |
|   |   | 0200-7015  | 384-well plates<br>Wavelength range: Visible to 650 nm                | 96-well (200 µL) 2 pM<br>384-well (50 µL) 3 pM   | 96-well (200 µL) 10 pM<br>384-well (50 µL) 15 pM                           | 1          |
|   |   | 0200-7014  | 96-well plates<br>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<br>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    China (Beijing) +86.10.6410.8669    Japan +81.120.993.656  
 United Kingdom +44.118.944.8000    China (Shanghai) +86.21.3372.1088    South Korea +82.2.3471.9531  
 Europe\* 00800.665.32860    Hong Kong +852.3971.3530

\*Austria, Belgium, Denmark, Finland, France, Germany, Ireland, Netherlands, Spain, Sweden and Switzerland