

# SpectraMax Paradigm

## Luminescence Detection Cartridges



**SpectraMax® Paradigm® Luminescence Detection Cartridges** utilize a patent pending design that isolates each well to provide optimal performance for glow-luminescence applications.

# SpectraMax Paradigm Multi-Mode Detection Platform

## Luminescence Detection Cartridges

### Maximum flexibility and sensitivity

- Three cartridges; each optimized for a popular plate format
- Choose one or have all three



### DESCRIPTION

Part Numbers ..... 0200-7012, 0200-7014, 0200-7015  
 Mounting ..... Top Read or Bottom Read  
 Cartridge Size ..... 1 Position  
 Emission Range ..... Visible to 800 nm

### TYPICAL PERFORMANCE

Detection limit for ATP (typical throughput /plate)

	0200-7012 (LUM)	0200-7014 (LUM 96)	0200-7015 (LUM 384)
<b>96 Well Plate (200 µl)</b>	0.5 fmol/well (2.5 min/plate)	<b>0.3 fmol/well</b> (2.5 min/plate)*	0.4 fmol/well (2.5 min/plate)
<b>384 Well Plate (50 µl)</b>	0.5 fmol/well (4 min)	NA	<b>0.2 fmol/well</b> (4min/plate)*
<b>1536 Well Plate (8 µl)</b>	<b>0.2 fmol/well</b> (10 min/plate)*	NA	NA

\*The cartridge's optical pickup is optimized for this plate format.

### TYPICAL APPLICATIONS\*\*

- Apoptosis
- cAMP Quantitation
- GPCR Ligand Binding
- GTP Binding
- Immunoassay

\*\*This list does not represent a complete list of applications that can be performed by the specific detection cartridges.



#### Contact Us

Phone: +1-800-635-5577  
 Web: www.moleculardevices.com  
 Email: info@moldev.com  
 Check our web site for a current listing of worldwide distributors.

#### Regional Offices

USA and Canada +1-800-635-5577  
 Brazil +55-11-3616-6607  
 China (Beijing) +86-10-6410-8669  
 China (Shanghai) +86-21-6887-8820  
 Germany +49-89/96-05-88-0

Japan (Osaka) +81-6-6399-8211  
 Japan (Tokyo) +81-3-5282-5261  
 South Korea +82-2-3471-9531  
 United Kingdom +44-118-944-8000