



## Luminescence

### SpectraMax Detection Cartridge

Luminescence SpectraMax® Detection Cartridges for SpectraMax® i3 and Paradigm® Systems utilize a patent-pending design that isolates each well to provide optimal performance for glow-luminescence applications.

# Luminescence SpectraMax Detection Cartridge

Maximum flexibility and sensitivity

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

## Description

Compatible systems ..... SpectraMax i3 and Paradigm Readers

Part Numbers ..... 0200-7012, 0200-7014, 0200-7015

Mounting ..... Top Read or Bottom Read<sup>†</sup>

Cartridge Size ..... 1 Position

Emission Range ..... Visible to 650 nm

<sup>†</sup> Bottom mounting for SpectraMax Paradigm Readers only.

## Technical specifications

Detection limit for ATP Glow

<b>0200-7012 (LUM)</b>	<b>Optimized</b>	<b>Guaranteed*</b>
96-well (200 $\mu$ L)	1.5 pM	2.5 pM
384-well (50 $\mu$ L)	6 pM	10 pM
1536-well (8 $\mu$ L)	19 pM	25 pM

  

<b>0200-7015 (LUM 384)</b>	<b>Optimized</b>	<b>Guaranteed*</b>
96-well (200 $\mu$ L)	1.2 pM	2 pM
384-well (50 $\mu$ L)	3 pM	4 pM

  

<b>0200-7014 (LUM 96)</b>	<b>Optimized</b>	<b>Guaranteed*</b>
96-well (200 $\mu$ L)	1 pM	1.5 pM

\* For properly functioning, operated, and maintained equipment.

## Key applications\*\*

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

\*\* This list does not represent a complete list of applications or labels which can be performed by the specific detection cartridge.



### 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  
Brazil +55-11-3616-6607  
China (Beijing) +86-10-6410-8669  
China (Shanghai) +86-21-3372-1088  
Germany 00800-665-32860

Japan (Osaka) +81-6-7174-8831  
Japan (Tokyo) +81-3-6362-5260  
South Korea +82-2-3471-9531  
United Kingdom +44-118-944-8000

