



Fluorescence Intensity SpectraMax Detection Cartridge

Fluorescence Intensity SpectraMax® Detection Cartridges for SpectraMax® i3 and Paradigm® Systems provide speed, sensitivity and flexibility. The cartridge design provides uncompromising performance for both top and bottom reads. In addition, the standard dual-emission design enables simple, straightforward FRET (fluorescence resonance energy transfer) measurement.

Fluorescence Intensity SpectraMax Detection Cartridge

Maximum flexibility and sensitivity

- Three cartridges; each optimized for popular fluorophores
- Choose one or have all three

Description

Compatible systemsSpectraMax i3 and Paradigm Readers

Part numbers.....0200-7002, 0200-7003, 0200-7004

MountingTop read or bottom read†

Cartridge size1 position

Light sourceHigh-power LED

† Bottom mounting for SpectraMax Paradigm Readers only.



Part Number	0200-7002	0200-7003	0200-7004	0200-7005
Cartridge Name	Coumarin-Fluorescein	Fluorescein-Rhodamine	Cy3-Cy5	CFP-FP
Wavelength-Tuned Excitation	360/35 nm	485/20 nm	535/25 nm	485/xx nm
Wavelength-Tuned Emission	465/35 nm 535/25 nm	535/25 nm 595/35 nm	595/35 nm LP 655 nm	485/xx nm 535/xx nm

Technical specifications

Detection limit for fluorescein (read time/plate)	Optimized	Guaranteed*
96-well	0.10 pM	0.75 pM
384-well	0.15 pM	1.3 pM
1536-well	0.4 pM	3.7 pM

Linear dynamic range > 6 decades

* For properly functioning, operated, and maintained equipment.

Key applications**

- Nucleic Acid Quantitation
- Molecular Interaction
- Ion Channel
- Protein Quantitation
- Reporter Gene Assays
- Apoptosis
- Proliferation/Viability

Key fluorophores**

0200-7002

- Coumarin
- Hoechst
- DAPI
- AMCA
- ANS
- Coumarin-Fluorescein (FRET)
- Alexa 350

0200-7003

- Fluorescein
- OliGreen
- PicoGreen
- RiboGreen
- Fluo-3,4
- GFP
- Fluorescein-Rhodamine (FRET)
- GFP-RFP (FRET) Alexa 488

0200-7004

- Rhodamine
- TAMRA, TMR
- Cy 3-Cy 5 (FRET)
- Cy 3-Cy 5.5 (FRET)
- Resorufin
- Alexa 532, 546, 555

0200-7005

- xxxxxxxxxxxx
- xxxxxxxxxxxx
- xxxxxxxxxxxx
- xxxxxxxxxxxx

** 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
 Email: 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

