

DATA SHEET

Fluorescence Intensity Detection Cartridges



Fluorescence Intensity Detection Cartridges for SpectraMax® i3x and Paradigm® Multi-Mode Microplate Readers 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.

Maximum flexibility and sensitivity

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

Key fluorophores**

0200-7002

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

KEY APPLICATIONS**

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

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

Description

Compatible systems	SpectraMax i3x and Paradigm Multi-Mode Microplate Readers
Mounting	Top read or bottom read*
Cartridge size	1 position
Light source	High-power LED

Part number

0200-7002	Cartridge name	Coumarin–fluorescein
	Wavelength-tuned excitation	360/35 nm
	Wavelength-tuned emission	465/35 nm 535/25 nm
0200-7005	Cartridge name	CFP-FP
	Wavelength-tuned excitation	485/xx nm
	Wavelength-tuned emission	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	

* Bottom mounting for SpectraMax Paradigm readers only.

* For properly functioning, operated, and maintained equipment.



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	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

