

SpectraMax Paradigm

Time-Resolved Fluorescence Detection Cartridge



The **SpectraMax® Paradigm® Time-Resolved Fluorescence Detection Cartridge** is the benchmark against which all other readers will be measured. Built to measure Europium and Samarium, it is ideal for your protein-protein interaction, GPCR and enzyme activity applications.

SpectraMax Paradigm Multi-Mode Detection Platform

Time-Resolved Fluorescence Detection Cartridge



DESCRIPTION

Part Number.....	0200-7008
Mounting.....	Top Read or Bottom Read
Cartridge Size.....	1 Position
Light Source.....	High-Power LED
Wavelength-Tuned Excitation	370/80 nm
Wavelength-Tuned Emission	616/10 nm 642/10 nm

TYPICAL PERFORMANCE

	Stop-Read	On-The-Fly
Time Resolved Fluorescence (Europium)		
384-well (75 µL)	5 amol (2 min)	8 amol (1 min)
1536-well (8 µL)	1.5 amol (4 min)	3 amol (1.5 min)

TYPICAL APPLICATIONS*

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

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

TYPICAL FLUOROPHORES*

- Europium Chelates
- Samarium Chelates



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