



Fluorescence Intensity (GeneBLAzer) SpectraMax Detection Cartridge

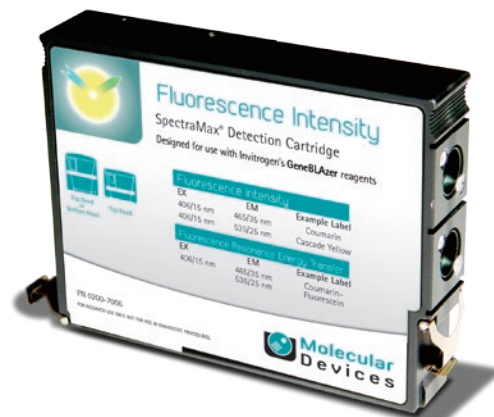
Fluorescence Intensity SpectraMax® Detection Cartridges for SpectraMax® i3 and Paradigm® Systems provide speed, sensitivity and flexibility. The cartridge design provides uncompromising performance for both top and bottom reads. In addition, the standard dual-emission design enables simple, straightforward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

Fluorescence Intensity (GeneBLAzer) SpectraMax Detection Cartridge

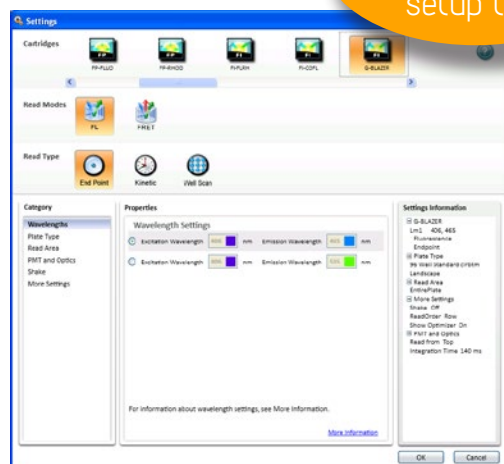
Designed to provide optimal performance when used with Life Technologies' GeneBLAzer reagents.

Description

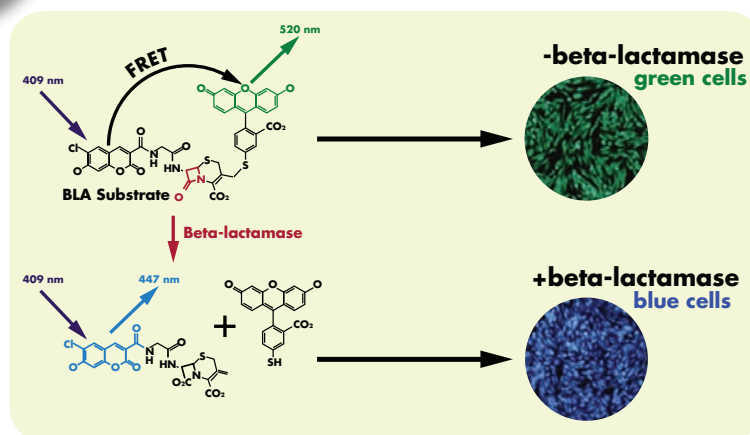
Compatible systems	SpectraMax i3 and Paradigm Readers
Part Numbers	0200-7006
Mounting	Top Read or Bottom Read
Cartridge Size	1 Position
Light Source	High-Power LED
Wavelength-Tuned Excitation	406/15 nm
Wavelength-Tuned Emission	EM1 465/35 nm
.....	EM2 535/25 nm



Fully enabled
protocols reduce
setup time



Principles of technology



Key applications

Cell-based assays

- GPCR Assays
- Ion Channel
- Nuclear and Cytokine Receptor Assays
- Signal Transduction Pathways
- Transcriptional Regulators (Kinases)

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
Brazil
China (Beijing)
China (Shanghai)
Germany

+1-800-635-5577
+55-11-3616-6607
+86-10-6410-8669
+86-21-3372-1088
00800-665-32860

Japan (Osaka)
Japan (Tokyo)
South Korea
United Kingdom

+81-6-7174-8831
+81-3-6362-5260
+82-2-3471-9531
+44-118-944-8000



**Molecular
Devices**