

# SpectraMax Paradigm

## Fluorescence Intensity (GeneBLAzer) Detection Cartridge



The **SpectraMax® Paradigm® Fluorescence Intensity Detection Cartridge** provides speed, sensitivity and exhibility. The cartridge design provides uncompromised performance for both top and bottom reads. In addition, the standard dual-emission design enables simple, straightforward FRET (uorescence resonance energy transfer) measurement.

# SpectraMax Paradigm Multi-Mode Detection Platform Fluorescence Intensity (GeneBLAzer) Detection Cartridge

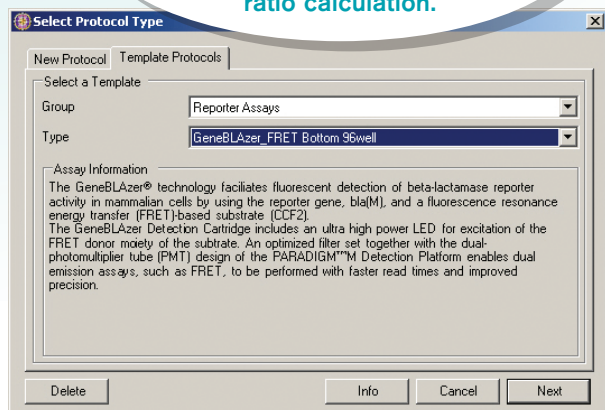
Designed to provide optimal performance when used with  
Invitrogen's GeneBLAzer® reagents.

## DESCRIPTION

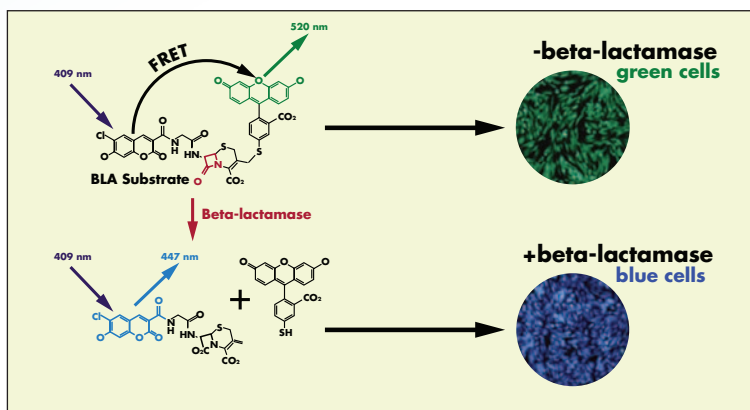
Part Numbers .....	0200-7006
Mounting.....	Top Read or Bottom Read
Cartridge Size.....	1 Position
Light Source.....	High-Power LED
Wavelength-Tuned Excitation .....	406/15 nm
Wavelength-Tuned Emission .....	EM1 465/35 nm EM2 535/25 nm



When inserted,  
template protocols are enabled  
with predened settings for a ratiometric  
readout, and data reductions such  
as blank subtraction and blue/green  
ratio calculation.



## Principles of Technology



## TYPICAL APPLICATIONS

Cell Based assays:

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



### 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 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