

SpectraMax Paradigm



Fluorescence Intensity Dual-Label (MultiTox-Fluor) Detection Cartridge



The innovative **SpectraMax® Paradigm®** **Fluorescence Intensity Dual-Label (FI-DL) (MultiTox-Fluor) Detection Cartridge** is perfect for efficient multiplexing of assays. It utilizes two high-powered LEDs to provide fast, sensitive, simultaneous reads of two distinct labels.

SpectraMax Paradigm Multi-Mode Detection Platform

Fluorescence Intensity Dual-Label (FI-DL) (MultiTox-Fluor) Detection Cartridge

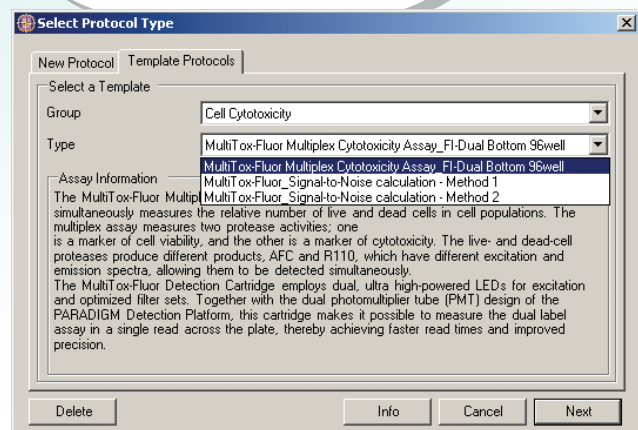
DESCRIPTION

Part Number.....	0200-7007
Mounting.....	Top Read or Bottom Read
Cartridge Size.....	1 Position
Light Source.....	2 High-Powered LEDs
Excitation Range.....	EX1 406/15nm
	EX2 504/12nm
Emission Range.....	EM1 465/35nm
	EM2 542/27nm

Designed to provide optimal performance when used with Promega's MultiTox-Fluor reagents. Reads both labels simultaneously.



**Fully enabled protocols=
reduced setup time**

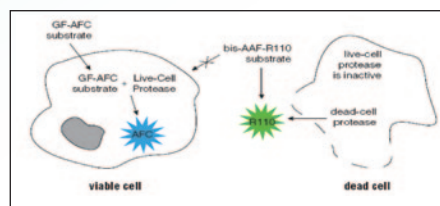


TYPICAL APPLICATIONS

The Fluorescence Intensity Dual-Label Detection Cartridge provides optimal performance for key applications for Drug Discovery, Cancer Research, ADME-Tox studies and others.

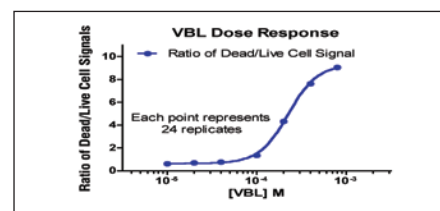
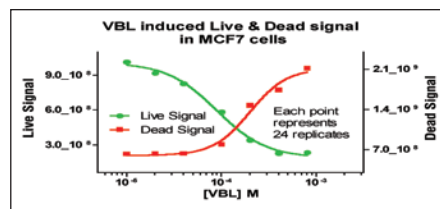
For more information and typical applications refer to www.promega.com

Principles of Technology



Source: <http://www.promega.com/tbs/tb348/tb348.pdf>

Vinblastine dose responses in MCF7 cells



"MultiTox-Fluor Multiplex Cytotoxicity Assay for the Breast Cancer Cell line, MCF7, using the BioRAPTR FRD Microfluidic Workstation and SpectraMax Paradigm Multi-Mode Detection Platform"



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