



## Fluorescence Intensity Dual-Label (Multi-ToxFluor)

SpectraMax Detection Cartridge

Fluorescence Intensity Dual-Label (FI-DL) (MultiTox-Fluor) SpectraMax® Detection Cartridges for SpectraMax® Paradigm® Systems are perfect for efficient multiplexing of assays utilizing two high-powered LEDs to provide fast, sensitive, simultaneous reads of two distinct labels.

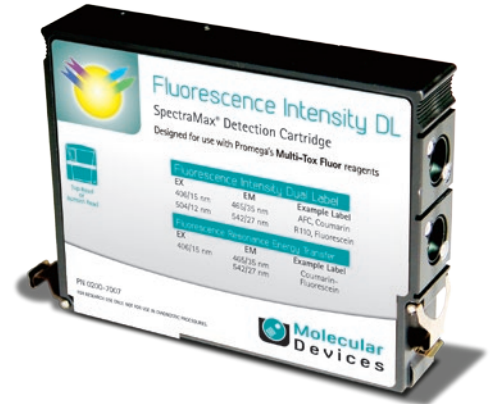
# Fluorescence Intensity Dual-Label (FI-DL) (Multitox-Fluor) SpectraMax Detection Cartridge

## Description

|                          |                                |
|--------------------------|--------------------------------|
| Compatible systems ..... | SpectraMax Paradigm Reader     |
| Part number .....        | 0200-7007                      |
| Mounting .....           | Top read or bottom read        |
| Cartridge size .....     | 1 Position                     |
| Light source .....       | 2 High-powered LEDs            |
| Excitation range .....   | EX1 406/15 nm<br>EX2 504/12 nm |
| Emission range .....     | EM1 465/35 nm<br>EM2 542/27 nm |

† Bottom mounting for SpectraMax Paradigm Readers only.

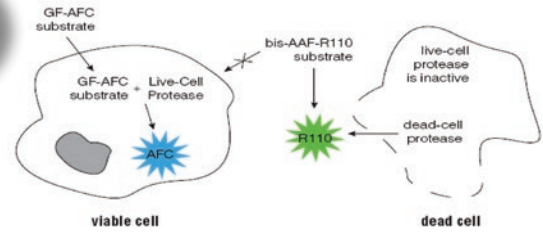
Designed to provide optimal performance when used with Promega's MultiTox-Fluor reagents. Reads both labels simultaneously. For more information and typical applications refer to [www.promega.com](http://www.promega.com).



Fully enabled protocols reduce setup time

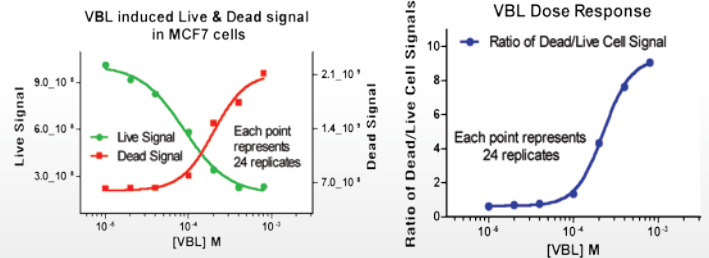


## Principles of technology



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.

## Key applications

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

## 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 website for a current listing of worldwide distributors.

## Regional Offices

|                  |                  |                |                  |
|------------------|------------------|----------------|------------------|
| USA and Canada   | +1-800-635-5577  | Japan (Osaka)  | +81-6-7174-8831  |
| Brazil           | +55-11-3616-6607 | Japan (Tokyo)  | +81-3-6362-5260  |
| China (Beijing)  | +86-10-6410-8669 | South Korea    | +82-2-3471-9531  |
| China (Shanghai) | +86-21-3372-1088 | United Kingdom | +44-118-944-8000 |
| Germany          | 00800-665-32860  |                |                  |

