



## Tunable Wavelength SpectraMax Detection Cartridge

TUNE SpectraMax<sup>®</sup> Detection Cartridges for the SpectraMax<sup>®</sup> Paradigm<sup>®</sup> System offer unmatched sensitivity in wavelength tunable detection. Designed to work with top or bottom read, it supports detection in:

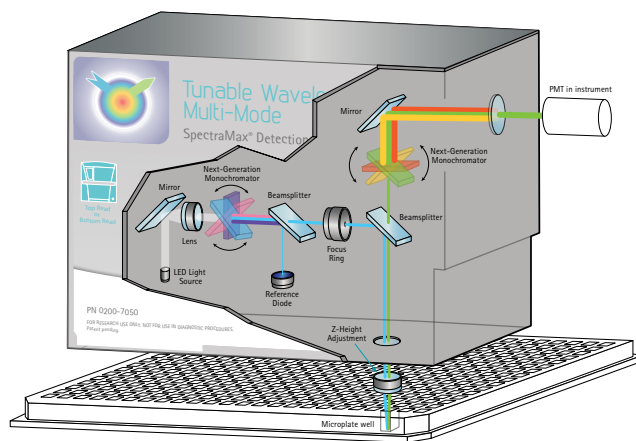
- Fluorescence Intensity (FI)
- Time-Resolved Fluorescence (TRF)
- Luminescence (LUM)

# Tunable Wavelength SpectraMax Detection Cartridge

## Description

The TUNE technology combines the wavelength flexibility of a monochromator-based system with the signal strength of a filter design. High-powered LEDs provide an adjustable light source, while the TUNE technology selects the excitation and emission wavelengths of your choice. The optimization wizard in SoftMax® Pro Software quickly scans excitation and emission spectra to select the ideal settings for the user's assay.

Compatible systems .....	SpectraMax Paradigm Reader
Part Number .....	0200-7050
Mounting .....	Top or Bottom Read
Cartridge Size .....	3 Positions
Light Source.....	High-Power LEDs
Wavelength Scanning Range .....	EX 360-790 nm EM 400-850 nm
Wavelength Selection .....	1 nm increments
Wavelength Accuracy .....	± 2 nm
Wavelength Repeatability.....	1 nm
Bandwidth .....	EX/EM - 20 nm



## Technical specifications

	Optimized	Guaranteed*	
Fluorescence Intensity (Fluorescein)			
96-well	0.2 pM	1.5 pM	
384-well	0.5 pM	2.5 pM	
1536-well	1.5 pM	6 pM	
Time Resolved Fluorescence (Europium)			
96-well	0.1 pM	0.43 pM	
384-well	0.1 pM	0.50 pM	
1536-well	0.3 pM	2.50 pM	
Luminescence (ATP)			
96-well	20 pM	40 pM	
384-well	50 pM	100 pM	
Optimized speed	<b>96-Well</b>	<b>384-Well</b>	<b>1536-Well</b>
Read times (seconds)	12	24	54

\* For properly functioning, operated, and maintained equipment

## Typical applications\*\*

- cAMP quantitation
- Cell migration
- Cell viability and proliferation
- Cytotoxicity
- DNA/RNA quantitation (fluorescent)
- ELISA (fluorescent/luminescent)
- Enzyme kinetics
- GPCR
- Immunoassay
- Neurotransmitter transporter uptake
- Protease
- QBT Fatty Acid Uptake
- Reporter assays

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

### 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	+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		

