

# SpectraMax Paradigm

## AlphaScreen Detection Cartridges



**SpectraMax® Paradigm® AlphaScreen® Detection Cartridges** utilize a 680 nm laser diode to provide a sensitive reading system for AlphaScreen assays. In addition, a patent pending design that isolates each well enables optimal performance for AlphaScreen assays.

# SpectraMax Paradigm Multi-Mode Detection Platform

## AlphaScreen Detection Cartridges

### DESCRIPTION

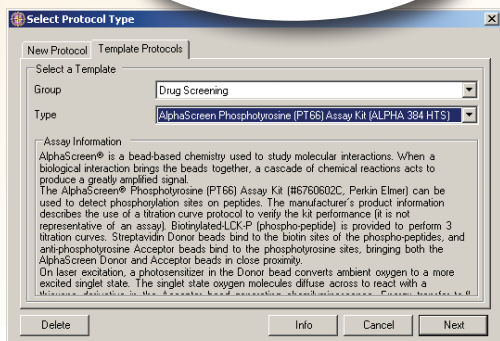
Part Numbers ..... 0200-7017, 0200-7018, 0200-7019  
 Mounting ..... Top Read or Bottom Read  
 Cartridge Size ..... 1 Position  
 Light Source ..... 680 nm Laser Diode  
 Excitation Range ..... 680 nm  
 Emission Range ..... 570/100 nm

### TYPICAL PERFORMANCE

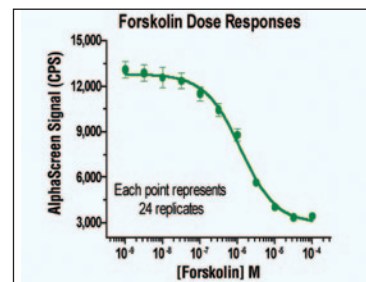
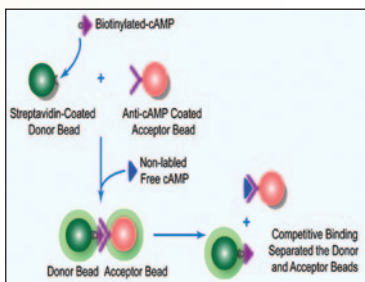
Detection limit <100 amol of biotinylated-LCK-P peptide, 25 µL/well in 384-well plate (with 0200-7017, 0200-7018), using AlphaScreen Phosphotyrosine (PT66) Assay Kit.

PART NUMBER:	0200-7017 (384 Std)	0200-7018 (384 HTS)	0200-7019 (1536 HTS)
Plate formats	96, 384	96, 384	96, 384, 1536
Typical Throughput	4 min/ 384 plate	2 min/384 plate	8 min/1536 plate

Fully enabled protocols = reduced setup time



### Principles of Technology



Source: "AlphaScreen cAMP Assay for the Human Epithelial Carcinoma Cell Line, A431, Using the BioRAPTR FRD Microfluidic Workstation and SpectraMax Paradigm Multi-Mode Detection Platform"

### TYPICAL APPLICATIONS

The SpectraMax Paradigm AlphaScreen Detection Cartridge provides optimal performance for key applications in drug discovery:

- GPCR Assays
- Cytokine Assays
- cAMP Quantitation
- Kinase Assays



#### 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