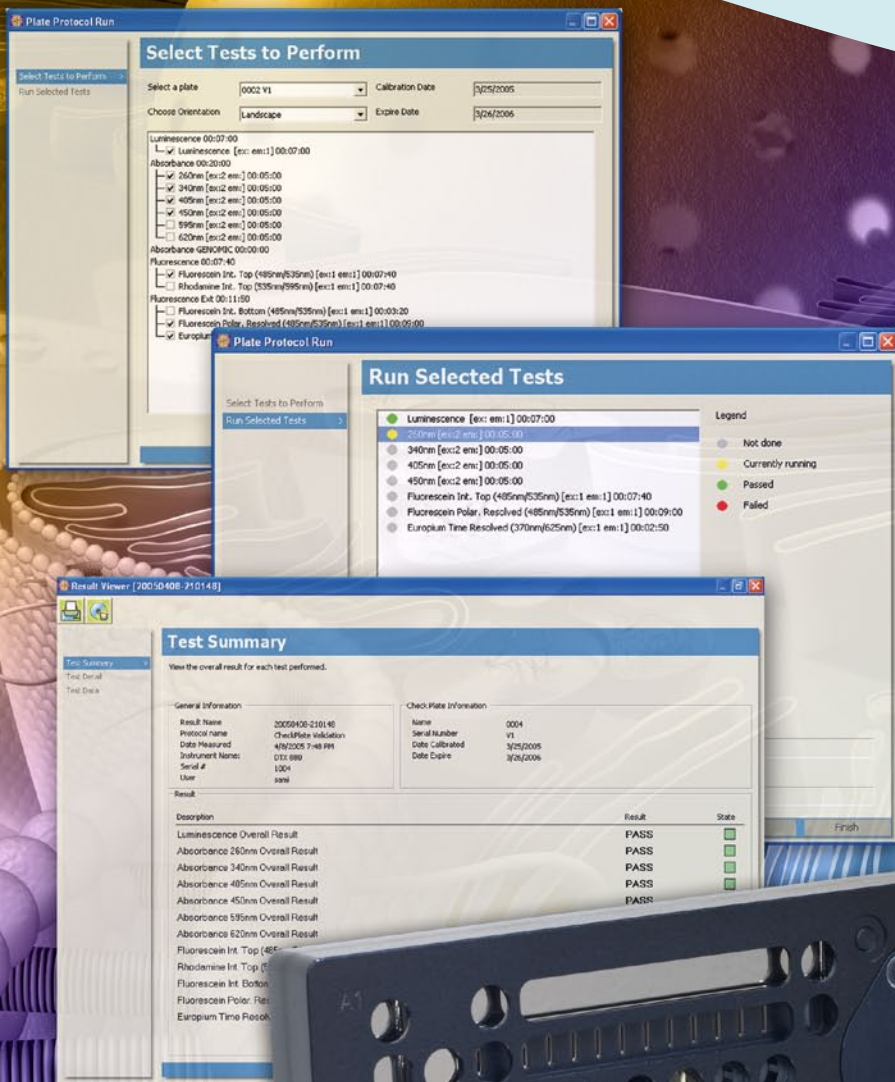


# Multi-Mode Validation Plate



for the SpectraMax Paradigm Multi-Mode Detection Platform and FilterMax Multi-Mode Microplate Readers

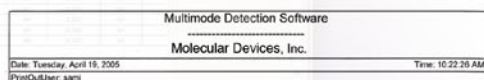
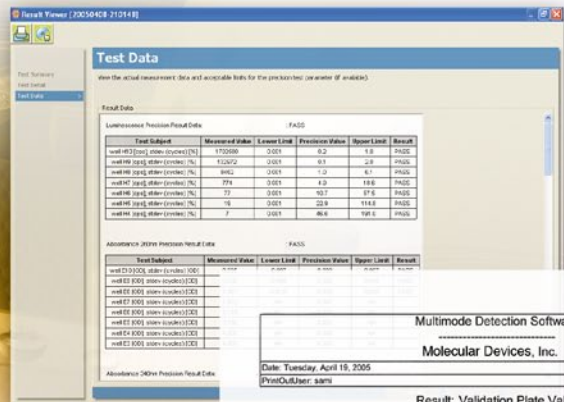
*Confident Assurance of  
Instrument Performance*



The innovative Multi-Mode Validation Plate, featuring a reference test plate and simple software, enables users to easily verify the performance of their Molecular Devices SpectraMax® Paradigm® Multi-Mode Detection Platform and FilterMax™ Multi-Mode Microplate Readers. The Multi-Mode Validation Plate provides testing for all available SpectraMax Paradigm System and FilterMax Reader detection modes on a single plate through an easy-to-use software interface. There is no programming of protocols or wet chemistry involved—just select the mode(s) you want to test, insert the plate and go. A standard summary test report delivers basic “Pass/Fail” results, or specific test data can be viewed if desired. The Multi-Mode Validation Plate is an ideal solution for any customer who wants confident assurance of their SpectraMax Paradigm System or FilterMax Multi-Mode Reader performance.

## COMPREHENSIVE PERFORMANCE TESTING – ALL ON ONE PLATE

- Ability to test one or a selection of the following detection modes:
  - Absorbance (ABS)
  - Fluorescence Intensity (FI) (top and bottom)
  - Luminescence (LUM)
  - Fluorescence Polarization (FP)
  - Time-Resolved Fluorescence (TRF)
- Tests for a variety of performance parameters, including:
  - Linearity
  - Precision
  - Accuracy
  - Dynamic Range
  - Alignment
  - Signal to Noise
  - Signal to Background
  - Filter Checks



Result Name : 20050408-194351  
 Protocol Name : ValPlate Validation  
 Application Type : Validation  
 Date Measured : Friday, April 08, 2005, 6:27 PM  
 User Measured : samf  
 Check Plate Name : 0001  
 Serial Number : V1  
 Date Calibrated : Friday, March 25, 2005, 12:00 AM  
 Date Expire : Sunday, March 28, 2005, 12:00 AM  
 Instrument Name : Test 1  
 Serial Number : 1019

#### Test Summary

Luminescence Overall Result : PASS  
 Absorbance 200nm Overall Result : PASS  
 Absorbance 340nm Overall Result : PASS  
 Absorbance 405nm Overall Result : PASS  
 Absorbance 450nm Overall Result : PASS  
 Absorbance 595nm Overall Result : PASS  
 Absorbance 620nm Overall Result : PASS  
 Fluorescein Int. Top (485nm/535nm) Overall Result : PASS

Beckman Coulter, Inc.

Page 1

## RESULTS YOU CAN TRUST

- Plate uses National Institute of Standards and Technology (NIST) traceable standards for absorbance testing
- Each plate comes with its own unique certification and quality documentation
- Annual re-certification program ensures plate and reference value reliability

## EASY OPERATION

- User chooses which mode(s) to test and how often to run the plate
- Simple software interface – no programming required
- No time-consuming pipetting or wet chemistry involved

## STRAIGHTFORWARD RESULTS REPORTING

- Standard 'pass/fail' results report printout
- View additional results data detail if desired
- Save test results in software database



## ORDERING INFORMATION

PART NO.	DESCRIPTION
0200-7200	Multi-Mode Validation Plate for FilterMax/SpectraMax Paradigm Systems
0200-7201	Multi-Mode Validation Plate Recalibration Program for FilterMax/SpectraMax Paradigm Systems



#### 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