

# SpectraMax Gemini XPS Microplate Reader

A dual-monochromator spectrofluorometer system

## **BENEFITS**

- · No filters needed
- · Higher sensitivity
- Validation tools
- · Robotics compatible

#### Introduction

The SpectraMax® Gemini™ XPS Microplate Spectrofluorometer from Molecular Devices® provides a flexible environment to determine the optimal excitation and emission settings for most fluorescent intensity assays. The Gemini XPS Reader has dual monochromators that allow users to utilize new and novel dyes without purchasing expensive filter sets. SoftMax® Pro Microplate Data Acquisition and Analysis Software, included with every Gemini XPS Reader, provides convenient data analysis. The optional SpectraTest® FL1 Validation Plate automatically validates the performance of the reader. Software validation, IQ/OQ/PQ and FDA 21 CFR Part 11 compliance tools are also available.

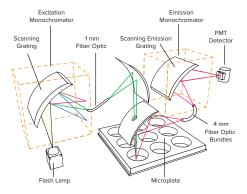
#### **Dual monochromators**

With the Gemini XPS Reader, users don't have to worry about having the right set of filters. The system uses two scanning monochromators to determine the optimal excitation and emission settings. Assays vary depending on temperature or pH conditions; the monochromators allow users to select the optimal wavelength for any assay. Alternatively, literature values for the monochromator can be easily entered. When methods or fluorophores change, it takes only a few software commands to adjust the reader.

#### Well scanning

With well scanning, multiple points within each well can be read, providing a high level of sensitivity for cell-based assays. Endpoint, kinetic and spectrum scanning assays can also be run.

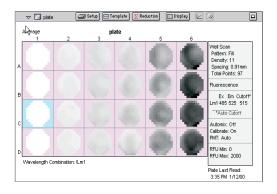




**Gemini XPS Reader optics.** The optics of the Gemini XPS Reader are engineered for superior performance and reliability.



**Streamline GLP/GMP compliance.** The SpectraTest FL1 Validation Test Plate provides a complete solution for validating optical performance of the Gemini XPS reader.



**Gemini XPS Reader scans.** Well scans from the Gemini XPS Reader illustrating serial dilution of CHO cells using the Live Cell assay from Molecular Probes Live/Dead (Calcein AM /Ethidium Homodimer-1) assay kit. A Costar 24-well tissue culture plate with col1 = blanks and doubling cell concentration left to right starting at Col. 2 with 1000 cells.

### Plate stacker and robot integration

The Gemini XPS Reader can be integrated with Molecular Devices' StakMax® Microplate Stacker in a matter of minutes and begin reading microplates with seven mouse clicks. For a higher degree of automation, the Automation Vendor Partners Program has streamlined the integration of our microplate reader systems with all leading partner robots. The "out-of-the-box" automation solution saves up-front integration time and resources.

#### Leading data analysis software

SoftMax Pro Microplate Microplate Data Acquisition and Analysis Software provides flexibility in experimental design, setup, analysis and reporting, providing the opportunity to customize assays. Choose from nine different curve-fitting routines and use default data reduction, or set up custom formulas for analysis. Data can be analyzed and combined from different plates.

#### **Applications**

- Live/Dead viability/cytotoxicity assays
- · Detection of nucleic acids
- · Green fluorescent protein
- NanoOrange protein quantitation
- · PicoGreen DNA detection
- · Molecular beacons
- · Caspase-3
- Fluorometric protease assays
- · cAMP detection

#### **Specifications** Fluorescence photometric performance specifications Dual monochromators 1 nm increment selection EX wavelength range 250-850 nm EM wavelength range 360-850 nm Wavelength bandwidth (EX, EM) 9 nm Detection limit 3.0 fmol/well FITC 200 µL in 96 wells (signal 3X SD of baseline) Time-resolved fluorescence specifications (secondary mode) 360-850 nm Wavelength range Data collection 50-1450 µsec., 200 µsec. increments 0.5 fmol/well Eu-chelate (obtained Sensitivity with DELFIA reagent from Perkin Elmer by using a 384-well plate) Luminescence specifications (secondary mode) 10 amol/well alkaline phosphatase $200 \, \mu L/well$ (obtained with Emerald II Detection limit reagent from Applied Biosystems) General photometric performance specifications 6, 12, 24, 48, 96, 384 wells Plate formats Light source Xenon flash lamp (1 joule/flash) Detector Photomultiplier (R-3896) Read time\* 96 wells < 15 sec., 384 wells < 45 sec. Shaker time 0 to 999 seconds Ambient +4°C to 45°C Temperature range **General specifications** Dimensions (in.) 8.6 (H) x 22.8 (W) x 15 (D) Dimensions (cm) 22 (H) x 58 (W) x 38 (D) Weight 35 lbs. (15.9 kg) Power consumption < 125 watts 100-240 Vac, 3 A, 50/60 Hz Power source Robot compatible

<sup>\*</sup> Measurement type may extend read time.

## Ordering information

Contact your Molecular Devices sales representative for configuration options.

#### 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.

