



DETECTION

GLOMAX[®] MULTI DETECTION SYSTEM

A modular, easy to use and cost effective multimode reader for Luminescence, Fluorescence, and Absorbance Life Science applications.



Promega

GLOMAX[®] - MULTI MICROPLATE MULTIMODE READER

OVERVIEW

The GloMax[®]-Multi Microplate Multimode Reader is designed for today's life science laboratory. In addition to giving performance equivalent to that of single-mode instruments, the GloMax[®]-Multi blends user-friendly operation with easy data handling and flexible purchasing options. The result of this design is an instrument with superior performance that is easy to use, is affordable, and can be customized to your laboratory's needs.

PERFORMANCE

The GloMax[®]-Multi Microplate Multimode Reader combines the superior performance expected from single-mode instruments with the functionality of multiple modes. To achieve industry-leading performance, the GloMax[®]-Multi is designed with optical channels dedicated to each individual technology. Unlike other multimode systems, readings taken with the GloMax[®]-Multi are not degraded by indirect fiber-optic transmission or crowded optical channels. Dedicated optical channeling ensures that the GloMax[®]-Multi provides sensitivity and dynamic range on par with that of the highest performing single-mode instruments.

EASE OF USE

The GloMax[®]-Multi Microplate Multimode Reader is designed to be put into use straight from the box, without the need for special training. The built-in computer eliminates the connectivity hassles, cost, and space requirements of running an external computer. Assay set-up on the GloMax[®]-Multi is effortless with several flexible set-up options to choose from.

AFFORDABLE MODULARITY

The GloMax[®]-Multi Microplate Multimode Reader is a modular instrument that easily fits into most budgets. Purchase the technology or modes that you need now and add onto the system later as your needs expand. For example, the GloMax[®]-Multi can be purchased as a luminometer, then fluorescence and/or absorbance modules can be purchased and added later.

Installations take less than ten minutes and can be done right in your lab using just the Allen wrench provided with the module. In addition, after installing a new module, you will not need to download and install new software. The GloMax[®]-Multi instantly detects the newly installed module(s) and will automatically adjust screens, protocols, and options.

INSTRUMENT FEATURES

DATA HANDLING

To transfer data to either PC or Mac, simply copy the data from the GloMax®-Multi then move the USB stick (included) to your computer. Analyze your data when and where you find it most convenient.

MICROPLATE FORMAT

Accepts 96-well plates conforming to the SBS plate standard.

COMPUTER INTERFACE (OPTIONAL)

For those labs that would prefer to operate the GloMax®-Multi through an external PC, the optional PC Software contains all of the same ease-of-use features that are available via the instrument's built-in touch screen.



INJECTORS

Single or dual auto injectors are available for flash-based luminescence applications or dual-reporter assays. Each injector system has a volume range of 25 – 200 μ l in 5 μ l increments. Injectors are controlled using the touch screen and fluidics wizard. PRIME and FLUSH commands provide easy maintenance and a REVERSE PURGE command saves valuable reagents.

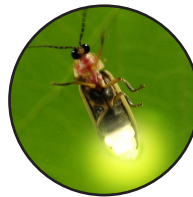
COLOR TOUCH SCREEN

The GloMax®-Multi combines a 6.6" color touch screen display with an onboard Windows-based computer.

INTUITIVE USER INTERFACE

Setting up a run and retrieving data are fast and simple while maintaining the flexibility needed for advanced or custom protocols.

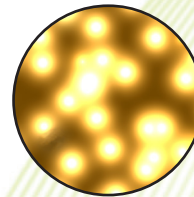
SEPARATE CHANNELS



Luminescence



Fluorescence



Absorbance

SEPARATE CHANNELS

LUMINESCENCE MODULE

(Factory Installed)

- >8 Logs Dynamic Range
- Dual-masking System Reduces Crosstalk
- 3×10^{21} moles Luciferase Detection or 1×10^{18} moles ATP



To achieve the sensitivity of a dedicated luminometer, the luminescence channel is separated from other measurement technologies and positioned directly above the sample well. These conditions maximize light capture for the best possible sensitivity. In addition, a low-noise photomultiplier tube ensures that collected light is not compromised in any way. This design makes the GloMax®-Multi Microplate Multimode Reader between 10 to 1000 times more sensitive than competing multimode luminometers. In addition to high sensitivity, the GloMax®-Multi boasts greater than 8 logs of dynamic range, eliminating the need to dilute samples or manage detector-driven gain changes. The photomultiplier tube automatically adjusts for the optimum reading of bright or dim samples. This means that the GloMax®-Multi is capable of achieving a reading range 2 - 3 logs more than that of competing multimode luminometers.

Luminescence Injector/Application Chart

Recommended Injectors	Applications
2	Dual Reporter Assays
0	Steady-Glo® Luciferase Assays
1	Flash Glow Luciferase Assays
1	Cell Viability/ATP Assays

1 Kinetics Assays

FLUORESCENCE MODULE

(User Installable)

- Epifluorescent Detection (top reading)
- Easy Optical Kit Switching
- Wavelength-matched LEDs Ensure High Sensitivity



To achieve high performance, the Fluorescence Module uses powerful light-emitting diodes (LEDs) as excitation sources. LEDs have very specific light-output profiles that closely match the excitation profiles of commonly used fluorescent molecules. LED usage increases sensitivity by fully exciting the fluorophore and reducing non-specific light leakage, a common problem when using broad-spectrum light sources. Five standard optical kits are available to measure the most popular fluorophores. In addition, custom optical kits can be readily made for non-standard applications. Optical kits can be easily exchanged in seconds and built-in software ensures that the installed optical kit matches the selected protocol. Designed as a user-installable module, you can either buy the Fluorescence Module with your GloMax®-Multi instrument or add it to your system later when you are ready to run fluorescence experiments.

Fluorescence Module Optical Kit Application Chart

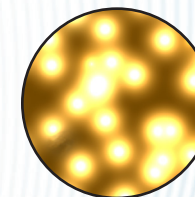
Optical Kit	Typical Applications
UV Optical Kit Ex: 365 nm Em: 410 - 460 nm	DNA Quantitation (Hoechst dye), Enzyme Activity (4-methyl-umbelliferone)
Blue Optical Kit Ex: 490 nm Em: 510 - 570 nm	DNA Quantitation, RNA Quantitation, Protein Labeling (Fluorescein), Protein Quantitation, Gene Expression (EGFP, rAcGFP)
Green Optical Kit Ex: 525 nm Em: 580 - 640 nm	Nucleotide or Protein Labeling (Rhodamine, Cy3), Enzyme Activity (Rhodamine)
Red Optical Kit Ex: 625 nm Em: 660 - 720 nm	Nucleotide or Protein Labeling (Cy5), RNA Quantitation

AFC Optical Kit
Ex: 405 nm
Em: 495 - 505 nm
Cell Viability, Cytotoxicity, and Apoptosis Multiplex Assays

ABSORBANCE MODULE

(User Installable)

- Flexible Filter System
- 0 - 5.0 OD Reading Range
- Reads Single or Dual Wavelengths



The GloMax®-Multi Microplate Multimode Reader with Absorbance Module provides measurements that are highly sensitive and cover a wide dynamic range. The absorbance channel has a reading range of 0 - 5.0 Optical Density (OD) with an accuracy that deviates less than 2%. The channel consists of a white LED with a spectral range of 450 to 750 nm positioned at the bottom of the instrument.

The Absorbance Module comes with two open filter slots and four factory-installed filters. The preinstalled filters are for 450, 560, 600, and 750 nm as these cover the most common ELISA and protein assays. With two open filter slots, you can also customize the absorbance module by adding the filter paddle of your choice.

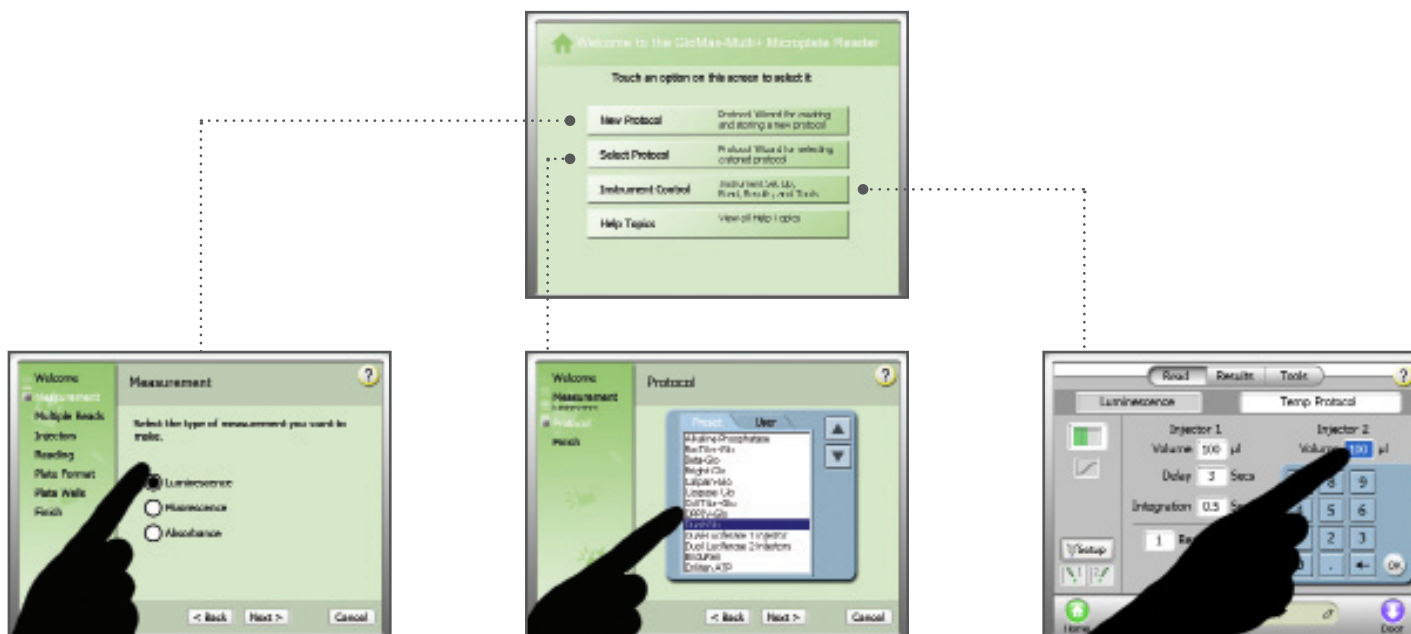
Absorbance Module Filter Application Chart

Wavelength	Applications
450	ELISA, QuantiCleave™ Protease Assay
560	BCA™ Protein Assays
600	Bradford Protein Assays, Coomassie Blue Protein Assays, PeroXOquant™ Quantitative Peroxide Assay
750	Lowry Protein Assay

INTUITIVE USER INTERFACE

TOUCH SCREEN SELECTION

The Windows-based computer built into the GloMax®-Multi Microplate Multimode Reader offers dynamic user-interface capabilities. Simply choose from three options on the instrument touch screen to start setting up a run.



CREATE A NEW PROTOCOL

Use the NEW PROTOCOL wizard to create a customized protocol. The wizard guides you step-by-step through choosing read parameters, injection methods, and plate wells to read, and in saving your protocol for future use.

CHOOSE A PRELOADED PROTOCOL

Choose SELECT PROTOCOL to access to the most popular assays from common reagent suppliers, or instantly access previously saved user-customized protocols.

TOTAL INSTRUMENT CONTROL

From the INSTRUMENT CONTROL screen, simply touch the parameter to change or customize all settings necessary to obtain the perfect read for your application. Once your parameters are set, you can save your settings for future use.

INSTRUMENT SPECIFICATIONS

GLOMAX® - MULTI

Available Detection Modes	Luminescence, Fluorescence, Visible Absorbance
Read Type	Glow, Flash, Kinetic, Repeat
Sample Format	96-well plates
User Interface	Built-in PC, touch screen navigation and operation
Data Output	PC or Mac compatible .csv file format exported to USB flash drive, or connect to PC (not included) through optional PC Connect Kit
External PC Requirements (optional)	Windows XP or Vista
Power	100 - 240 VAC, 50 - 60 Hz
Auto Shutoff	Touch screen hibernates after 15 minutes of inactivity
Dimensions	21" D x 17.3" W x 12.83" H (53 cm D x 44 cm W x 32.6 cm H)
Weight	~35 lbs (~16 kg)
Operating Temperature	60 - 85 °F (15 - 30 °C)
Warranty	One year parts and labor
Approvals	CE

LUMINESCENCE

Detector	Head-on photomultiplier tube (PMT) for photon counting
Spectral Range	350 - 650 nm
Peak Wavelength	420 nm
Detection Limit	3×10^{-21} moles of luciferase or 1×10^{-18} moles of ATP
Linear Dynamic Range	>8 logs
Crosstalk	5×10^{-5}

SINGLE INJECTOR SYSTEM

Number of Injectors	One injector
Dispense Volume Range	Selectable between 25 - 200 μ l in 5 μ l increments.
Waste Tray Volume	~50 ml

DUAL INJECTOR SYSTEM

Number of Injectors	Two injectors
Dispense Volume Range	Selectable between 25 - 200 μ l in 5 μ l increments.
Waste Tray Volume	~50 ml

FLUORESCENCE

Light Source	Wavelength-matched LED
Detector	PIN-photodiode
Read Position	Top reading
Wavelength Selection	Snap-in Fluorescence Optical Kits
Wavelengths	UV (Ex 365 nm: Em 410 - 460 nm), Blue (Ex 490 nm: Em 510 - 570 nm), Green (Ex 525 nm: Em 580 - 640 nm), Red (Ex 625 nm: Em 660 - 720 nm), AFC (Ex 405 nm: Em 495 - 505 nm)
Detection Limit	0.5 fmol/200 μ l or 1 ppt of fluorescein in 96-well plate, 30 pg/well dsDNA with DNA Quantitation Dye
Linear Dynamic Range	6 logs, assay dependent
Read Out	Relative Fluorescence units
Discrete Sample Average	Sample readings averaged over 0.5 second to improve accuracy

VISIBLE ABSORBANCE

Light Source	LED
Detector	Large-area photodiode
Spectral Range	450 - 750 nm
Wavelengths for Installed Filters	450, 560, 600, 750 nm, 2 slots for customization
Photometric Measuring Range	0 - 5.0 OD
Linear Dynamic Range	0 - 4.0 OD, assay dependent
OD Accuracy	0.01 OD \pm 3% at \leq 2.5 OD
OD Precision	0.01 OD \pm 1%
Stray Light	0.002% @ 560 nm in clear bottom, black wall plate

For research use only. Not for use in diagnostic procedures. GloMax is a trademark of Promega Corporation. All other trademarks are the sole property of their respective owners. For the most up-to-date specifications, visit www.promega.com.

