

Ordering Information

Catalog # Description

Lumimark Microplate Luminescence and Absorbance Systems

170-7170	Lumimark Microplate Luminescence and Absorbance System, 100–240 V, 50/60 Hz, power cord, serial cable, Lumimark software, instructions
170-7172	Lumimark System, with 2 liquid dispensers
170-7175	Lumimark Plus Microplate Luminescence and Absorbance System, with onboard software control and data reduction
170-7177	Lumimark Plus System, with 2 liquid dispensers
170-7173	One Dispenser Upgrade Kit, Lumimark or Lumimark Plus systems
170-7174	Two Dispensers Upgrade Kit, Lumimark or Lumimark Plus systems
170-7178	Lumi Test Plate Kit, for Lumimark or Lumimark Plus systems

Lumimark Software Requirements

Minimum PC requirements include a Pentium® 166 PC computer, 64 MB RAM, 240+ MB hard drive, 256-color monitor, serial port, compatible 256-gray-level graphics card or equivalent local bus, expansion slot, and Windows® 95, 98, NT, or 2000 platforms.

Lumimark System Specifications

Luminescence

Detection system:	Photon-counting photomultiplier tube
Spectral sensitivity range:	300–700 nm
Sensitivity at 545 nm:	50 amol ATP
Dynamic range:	8 orders
Filters:	Filterwheel for up to 15 luminescence and absorbance filters
Sample formats:	Opaque and transparent microplates, filter-bottom microplates, or other sample arrays of similar dimension
Dispensers:	Optional 1- or 2-dispenser systems Standard volume/stroke: 50 µl (metal free, exchangeable tubing)

Absorbance

Spectral range:	405–690 nm
Filter capacity:	Up to 15 filters
Indication range:	0–4.5 OD
Resolution:	0.001 OD
Accuracy:	±1% and 0.005 OD at 1 OD

Physical

Width:	50 cm
Depth:	34 cm
Height:	26 cm
Weight:	~14 kg
Computer interface:	RS232 port

Excel, Windows, and Windows 95, 98, NT, and 2000 are trademarks of Microsoft Corp. L2Win is a trademark of Anthos Labtec Instruments. Pentium is a trademark of Intel Corp.



The Lumimark Plus system with software onboard.

Luminescence
for the sensitive
types



For Glow- or Flash-Type Chemistries,
Lumimark Microplate Luminescence and Absorbance Systems
Go to the Extremes.

BIO-RAD

**Bio-Rad
Laboratories**

Life Science
Group

Web site www.bio-rad.com USA (800) 4BIORAD Australia 02 9914 2800 Austria (01)-877 89 01 Belgium 09-385 55 11 Brazil 55 21 507 6191
Canada (905) 712-2771 China 86-10-8201-1366/68 Denmark 45 44 52-1000 Finland 358 (0)9 804 2200 France 01 47 95 69 65 Germany 089 318 84-177
Hong Kong 852-2789-3300 India (91-124) 6398112/113/114 Israel 03 951 4124 Italy 34 91 590 5200 Japan 03-5811-6270 Korea 82-2-3473-4460
Latin America 305-894-5950 Mexico 52 5 534 2552 to 54 The Netherlands 0318-540666 New Zealand 64-9-4152280 Norway 47-23-38-41-30
Russia 7 095 979 98 00 Singapore 65-2729877 Spain 34-91-590-5200 Sweden 46 (0)8-55 51 27 00 Switzerland 061-717-9555 United Kingdom 0800-181134

BIO-RAD

Lumimark systems are sensitive flexible to light to your needs

Lumimark systems offer a high level of flexibility and value. High-sensitivity Lumimark systems are ideal for long-lasting glow-type luminescence, and can be optimized for flash- or glow-type luminescence with optional sample injectors. In addition to all types of luminescence detection, Lumimark can perform standard absorbance assays. The Lumimark Plus system has software onboard.

Features	Benefits
Absorbance capabilities	Ideal for use with multiple detection methods
Superior sensitivity	The instrument is so sensitive that a black plate can be used to improve the signal-to-noise ratio; the same assay on another instrument needs a white plate
Optional liquid dispenser	System is ideal for both flash- and glow-type chemistries
Optional onboard control	Saves space when necessary
RS232 port for computer control	Easy setup
Customizable lag time and wait time	Ideal for kinetic and low-background readings



Lumi Test Plate kit

The Lumi test plate kit, which includes a luminescence test plate and its accessories, is used to verify performance of Lumimark or Lumimark Plus systems.

The following performance parameters can be verified with the Lumi test plate:

Dynamic Range

The dynamic range covers 5.5 orders of magnitude, and can be checked by taking the reading from row A.

Cross-Talk

Cross-talk is checked by measuring a high-signal well and the 8 surrounding wells, which have no luminescence sources. The ratio of the highest value of the 8 surrounding wells to the high-signal well must be $\leq 5 \times 10^{-5}$.

Reproducibility

The Lumi test plate kit checks reproducibility by calculating the coefficient of variation, which must be $\leq 1\%$, for 8 measurements on one well.



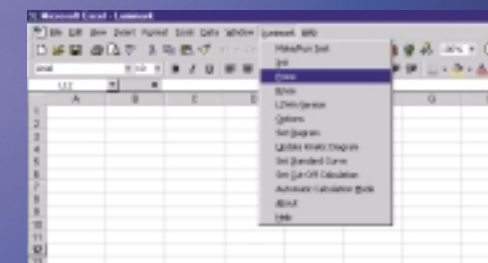
Lumimark Software

You can define tests to be executed on the Lumimark system, run them on the instrument, and receive the related measurements to evaluate the results.

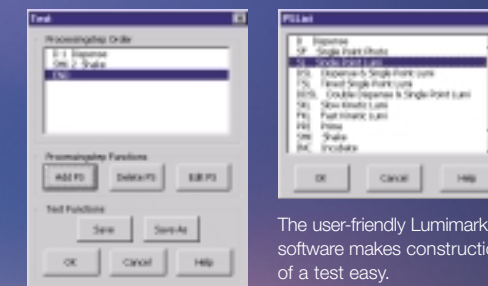
Lumimark software is designed as an Excel® macro that leads you through the steps to

set up templates and tests, and to have them performed by the instrument. Through a menu structure, it's simple to transfer your specific applications into a test.

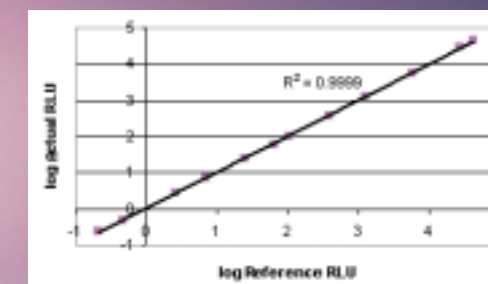
L2Win™ software is used to communicate with the instrument. Whenever the software establishes a connection, a window appears on the screen, providing information on the current communication between the computer and the instrument. This window closes after the command has been processed.



Lumimark software allows you to control the instrument with an Excel-based application, so data reduction can be accomplished in a familiar environment.



The user-friendly Lumimark software makes construction of a test easy.



Accurate linear dynamic luminescent reading with the Lumi test plate kit.