

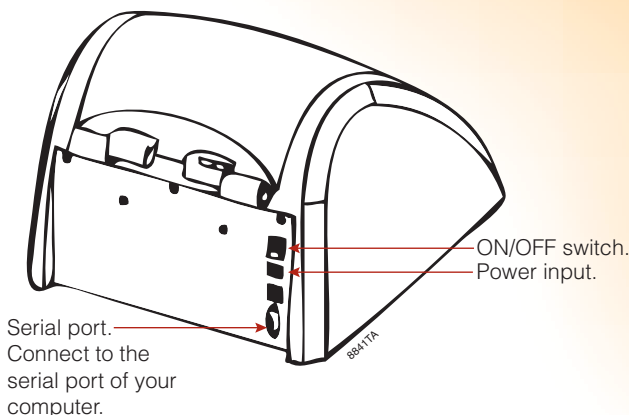
GloMax®-Multi Jr Detection System: Luminometer

INSTRUCTIONS FOR USE OF PRODUCT E6070 WITH LUMINESCENCE MODULE.

Quick
PROTOCOL

Setup

1. Connect the AC Adapter to the instrument.
2. Install the Luminescence Module.
3. Turn ON the instrument. A 60-second warm-up period optimizes performance.
4. (Optional) Connect the instrument to a PC using the 9-pin serial cable.
5. (Optional) To export data from the GloMax®-Multi Jr to an Excel® spreadsheet, you will need to install the Spreadsheet Interface Software on a PC.



Choose a Protocol

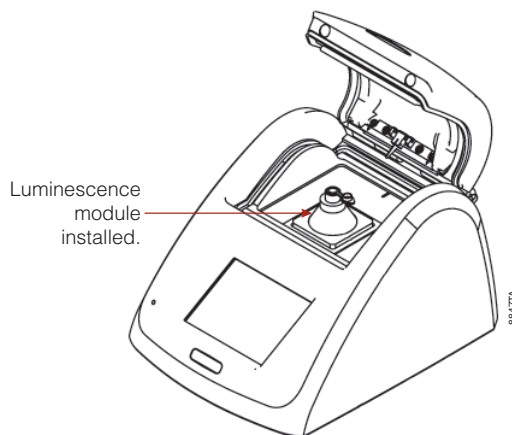
1. Touch "Protocol" to run a Promega protocol, create a new protocol, or access a saved protocol. You may also choose to run the GloMax®-Multi Jr default protocol. The default protocol provides a simple method to measure samples for an integration (measurement) time of 1 second.
2. Touch "Create New Protocol" to set up your own protocol. The Measurement Parameter screen will list the different options available.



Luminescence Module

Measure Samples

1. Insert the microcentrifuge tube containing your sample into the Luminescence Module.
2. Close the instrument lid, and touch the "Measure Luminescence" bar at the top of the touch screen. **Note:** Do not touch the screen after measurement starts. Touching any part of the screen during a protocol run will abort the measurement.
3. The data will appear on the touch screen. Use the arrow keys to scroll through the data. Only 20 data points are stored at a time.
4. To collect data from kinetic measurements, connect the GloMax®-Multi Jr to a PC or a printer.
5. Touch the "SampleID" key to use the SampleID feature. You may name samples individually using the touch screen keypad.
6. Touch the "Show Parameters" key to adjust your protocol parameters. After adjustment, touch "OK" to return to the Home screen.
7. To save a protocol for future use, touch the "Show Parameters" key, then the "Save As" key. Use the keypad to name your protocol and store it for future use.



Continued on other side

ORDERING/TECHNICAL INFORMATION:

www.promega.com • Phone 608-274-4330 or 800-356-9526 • Fax 608-277-2601

GloMax®-Multi Jr Detection System: Luminometer

INSTRUCTIONS FOR USE OF PRODUCT E6070 WITH LUMINESCENCE MODULE.

Quick
PROTOCOL

Clean the Touch Screen

1. Record all the data collected, and save the calibration or protocol.
2. Turn OFF the instrument.
3. Clean the touch screen with a delicate task wipe dampened with 70% ethanol.
4. Turn ON the instrument to resume measurements.

Help

Two Quick Ways to Access Help

1. From the instrument touch screen, touch the “?” button to view the “Help” Menu.
2. Contact Promega Technical Services at: techserv@promega.com



AF9FB120 0310FB120

ORDERING / TECHNICAL INFORMATION:

www.promega.com • Phone 608-274-4330 or 800-356-9526 • Fax 608-277-2601

©2010 Promega Corporation. All Rights Reserved.




Promega

Printed in USA. 3/10
Part #9FB120