

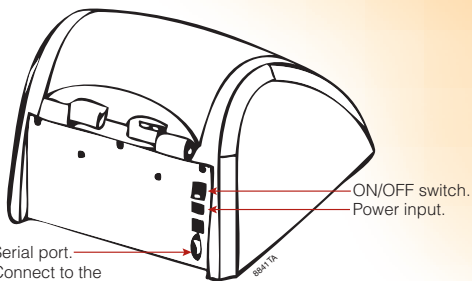
GloMax®-Multi Jr Detection System: Photometer

INSTRUCTIONS FOR USE OF PRODUCT E6070 WITH ABSORBANCE MODULE.

Quick
PROTOCOL

Setup

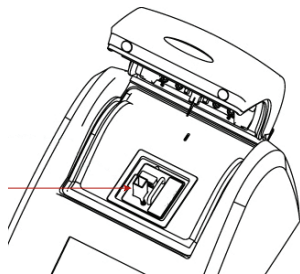
1. Connect the AC Adapter to the instrument.
2. Install the Absorbance Optical Kit.
3. Turn ON the instrument.
4. (Optional) Connect the instrument to a PC using the 9-pin serial cable.
5. (Optional) To export data from the GloMax®-Multi Jr to an Excel® spreadsheet, you will need to install the Spreadsheet Interface Software on a PC.



Optical Kit Installation/Removal

1. Turn OFF the instrument.
2. Hold the handle of the Optical Kit, and align it with the sample compartment. Press down to lock the Optical Kit in place.
3. Close the lid and turn ON the instrument.
4. Select "Photometer" from the list of options on the instrument touch screen.
5. Install the filter paddle that corresponds to the wavelength of absorbance for the assay.
6. Before removing an Optical Kit, turn OFF the instrument. Grasp the handle, and gently pull up to remove the kit from the sample compartment.

Serial port.
Connect to the serial port of your computer.

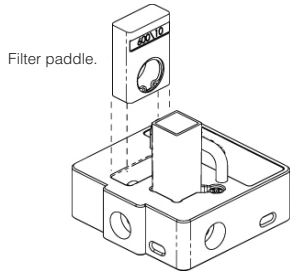


Measure Samples

Blank Instrument

Blank the GloMax®-Multi Jr after powering ON and changing filters. For best results, set a blank immediately before reading a series of samples using a buffer that is equivalent to your sample buffer. Blank values are not stored between power cycles.

1. Touch "Blank Instrument" to update the blank.
2. Insert a 10 × 10mm cuvette containing the appropriate buffer (blank).
3. Touch "OK" to complete the update.
4. Touch the "Cancel" key to abort the calibration.



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: Photometer

INSTRUCTIONS FOR USE OF PRODUCT E6070 WITH ABSORBANCE MODULE.

Quick
PROTOCOL

Measure Samples

1. Open the lid, and insert the cuvette. Close the lid.
2. Touch "Sample ID" to name your sample (optional). Using the keypad, enter the sample name into the name field. Touch "Save" to save the sample ID.
3. Touch "Measure Absorbance" to begin measurement. The GloMax®-Multi Jr will measure the sample for 5 seconds and report the average reading for the sample.

Note: The GloMax®-Multi Jr does not store more than 20 measurements at one time. Measurements are not stored between power cycles.

Save Calibrations

1. You may save calibrations for future use. After the calibration is complete, enter a name for the calibration and touch "Save As".
2. Touch the "Calibrate" key on the navigation bar to retrieve a saved calibration.

Clean the Touch Screen

1. Turn OFF the instrument.
2. Clean the touch screen with a delicate task wipe dampened with 70% ethanol.
3. 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



AF9FB122 0310FB122

ORDERING/TECHNICAL INFORMATION:

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