



GloMax[®]-Multi+ Detection System with Instinct[™] Software

Setting Up Your Instrument. It's Easy!

1. Unpacking the instrument

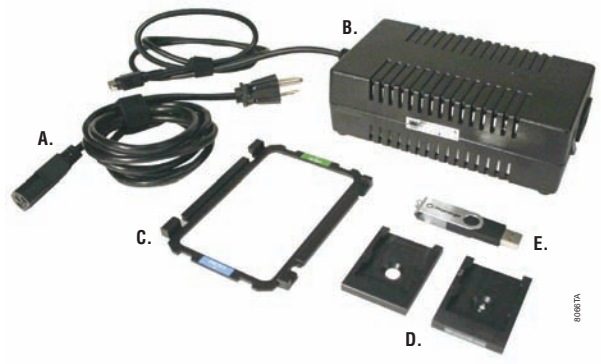


For the complete protocol see Technical Manual, TM340, on the CD provided with your instrument or go to: www.promega.com/protocols and search for TM340.

CAUTION! The GloMax[®]-Multi+ Detection System weighs approximately 35lb (16kg) and should be handled by two people.

1. Set the shipping box on a table, and remove the accessories and literature.
2. Carefully remove the GloMax[®]-Multi+ Detection System from the box and place on a flat, level, solid surface in a dust-free location with good air circulation. Leave at least 7.5 inches (19cm) of clearance in front of the instrument to allow the instrument door to open without hindrance.
3. Install the Instinct Software[™] on a PC/Windows[®] computer.

2. What's in the box?



The GloMax[®]-Multi+ Detection System with Instinct Software Accessories:

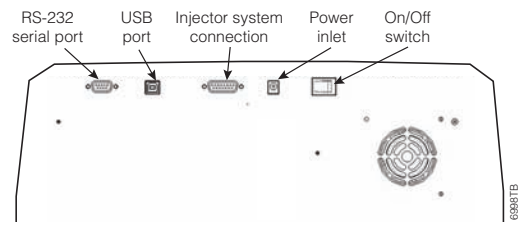
A. Power Cord. **B.** AC Power Adapter. **C.** Microplate Adapter, located inside the instrument. **D.** Two Optical Well Crosstalk Masks, one for 96-well plates and one for 384-well plates. The 96-well Optical Crosstalk Mask will be located inside the instrument. **E.** USB Flash Drive (appearance may vary from Figure E).

Included but not shown: CD containing GloMax[®]-Multi+ Detection System Technical Manual (TM340) and Instinct software, one spare red Breakaway Tab for the plate mechanism and a small screwdriver.

Check to make sure all accessories are included with the instrument and are undamaged.

3. Setting up the instrument

1. Open the accessories box included with your shipment, and locate the instrument power cable.
2. Connect the AC Power Adapter to the power inlet on the back of the Instrument. Connect the power cord to the AC Power Adapter.
3. Plug the power cord into a grounded outlet.

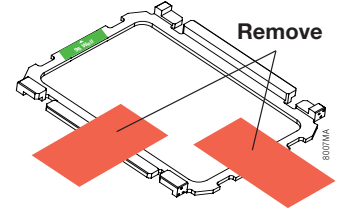


4. Starting up the instrument



1. Turn On the GloMax[®]-Multi+ Detection System. The power switch is located on the back panel next to the power connection. An orange LED located left of the touch screen will illuminate.
2. The instrument will initialize and calibrate for about one minute, at which point the touch screen will activate and the Home screen will appear.
3. In order to prevent damage to the instrument and loss of sample it is important to have the Microplate Adapter in the correct orientation. Please refer to Technical Manual TM340 for complete instructions.

Important: When using the instrument for the first time, you must press the "Door" button and remove the red labels attaching the Microplate Adapter to the Microplate Sample Tray.



Please refer to Technical Manual TM340 for detailed instructions on your specific application.

Continued on other side.





GloMax[®]-Multi+ Detection System with Instinct[™] Software

5. Setting the read parameters

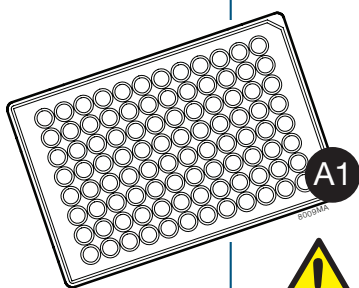
From the Home screen:

- For a wizard-guided protocol setup, select the “New Protocol” button.
- Alternatively, select a preprogrammed/saved protocol by selecting the “Select Protocol” button.
- Or, select the “Instrument Control” button to set new parameters.
- Select “Help Topics” to view all help topics.

6. Starting a run

Please refer to Technical Manual TM340 for Touch Screen Basics.


1. Touch the “Plate Mapping” button and highlight the wells to be read (green is selected and gray is deselected). For information about using the built-in curve fitting and well labeling features of the Instinct software see Technical Manual TM340.
2. Touch the “Door” button to open the instrument. Place the microplate in the sample tray. Make sure the A1 well is at the top right corner.
3. If installed, do not use force or push on the Microplate Sample Tray Cover to close it. The cover will close automatically.
4. Touch the “Start” button.
5. Touch the “Edit” button, and use the keypad to enter a file name.
6. Touch the “OK” button to start reading.




Important: To prevent evaporation and damage to the detector head, always remove the microplate when finished reading.

7. Results

To view Results:

1. From the Instrument Control screen, touch the “Results” button.
2. Touch the  button, and choose a saved file to view.

To Transfer Results File(s):

1. Insert a USB flash drive into the USB port on the front of the instrument to the left of the touch screen.
2. Touch the  button on the upper right corner of the screen.
3. Touch the “OK” button to export only the selected files, or touch the “Copy all Files” button to export all displayed and undisplayed results files.
4. To copy more than one file, touch the “Files” button, touch the Multi Select button and highlight multiple files to copy.
5. Touch the “USB” button to save to USB.
6. Move the USB flash drive to a PC/Windows[®] computer where the Instinct software has been installed to download files (see Technical Manual TM340 for more information about viewing the results).



Important: Only the 50 most recent results files are displayed on the touch screen. All results files are stored on the internal memory of the instrument. To ensure that data is saved, save the results files to a USB flash drive frequently.

8. Helpful resources

Detailed instructions on setting up and operating your GloMax[®]-Multi+ Detection System can be found in Technical Manual TM340, included on the CD that shipped with your instrument and available at:

www.promega.com/protocols Helpful sections of TM340 include:

- | | |
|---|--|
| • Touch Screen Basics | • Fluorometer Operation |
| • GloMax [®] -Multi+ Detection System with Shaking | • Switching Fluorescent Optical Kits |
| • GloMax [®] -Multi+ Detection System with Temperature Control | • Installing Injectors |
| • Luminometer Operation | • Removing and Installing the Microplate Sample Tray Cover |
| • Visible Absorbance Photometer Operation | • Troubleshooting |

