

GloMax®-Multi Detection System: Luminometer

INSTRUCTIONS FOR USE OF PRODUCT E7041.

Quick
PROTOCOL

Startup

1. Turn on the GloMax®-Multi Detection System Instrument (the switch is located on the back panel of the instrument).
2. The built-in PC and Windows® CE will load.
3. After warming up, the “Home” screen will appear.
4. Touch “Instrument Control”, and toggle to the Luminescence detection mode.
5. Touch the “PMT Activation” button.

Note: PMT activation takes about 5 minutes. Once activated, the PMT stays activated until the instrument is switched to fluorescence or absorbance mode, or the door is left open for more than one minute.

Select a Protocol

From the “Home” screen:

1. Setup a new protocol using the “New Protocol” wizard.
- OR –
2. Select a saved protocol using the “Select Protocol” wizard.

Injectors (When Needed)

Injectors must be initialized prior to use.

1. Place the reagent bottle in the holder, and place the stainless steel inlet tube into the bottle.
2. Touch the “Setup” button, and follow the step-by-step instructions to Prime, Flush, or Reverse Purge. Injectors must be primed prior to use.

Note: Flush the injectors after use. Check injector tips periodically. Immediately replace bent or damaged tips.

Start a Run

1. Touch the “Plate” button, and highlight the wells to be read (green wells are selected and gray wells are deselected).
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. Do not use force or push on the Microplate Sample Tray Cover to close. 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.

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

ORDERING/TECHNICAL INFORMATION:

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

©2008–2010 Promega Corporation. All Rights Reserved.



Printed in USA. Revised 5/10
Part #9FB089

GloMax®-Multi Detection System: Luminometer

INSTRUCTIONS FOR USE OF PRODUCT E7041.

Quick
PROTOCOL

Viewing and Saving Results

To View Results

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

To Transfer Results File(s)

1. Insert a USB flash drive into the USB port to the left of the touch screen.
2. Touch the “USB” button on the upper right corner of the screen.
3. Touch the “OK” button to export only the selected file, or touch the “Copy All Files” button to export all displayed and undisplayed results files.
4. Move the USB flash drive to a PC or Macintosh® computer to download files.

Note: 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 no loss of data, save the results files to a USB flash drive frequently.

For complete protocol information, see the *GloMax®-Multi Detection System Technical Manual #TM297*, available at: www.promega.com/tbs

ORDERING/TECHNICAL INFORMATION:

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

©2008–2010 Promega Corporation. All Rights Reserved.




Promega

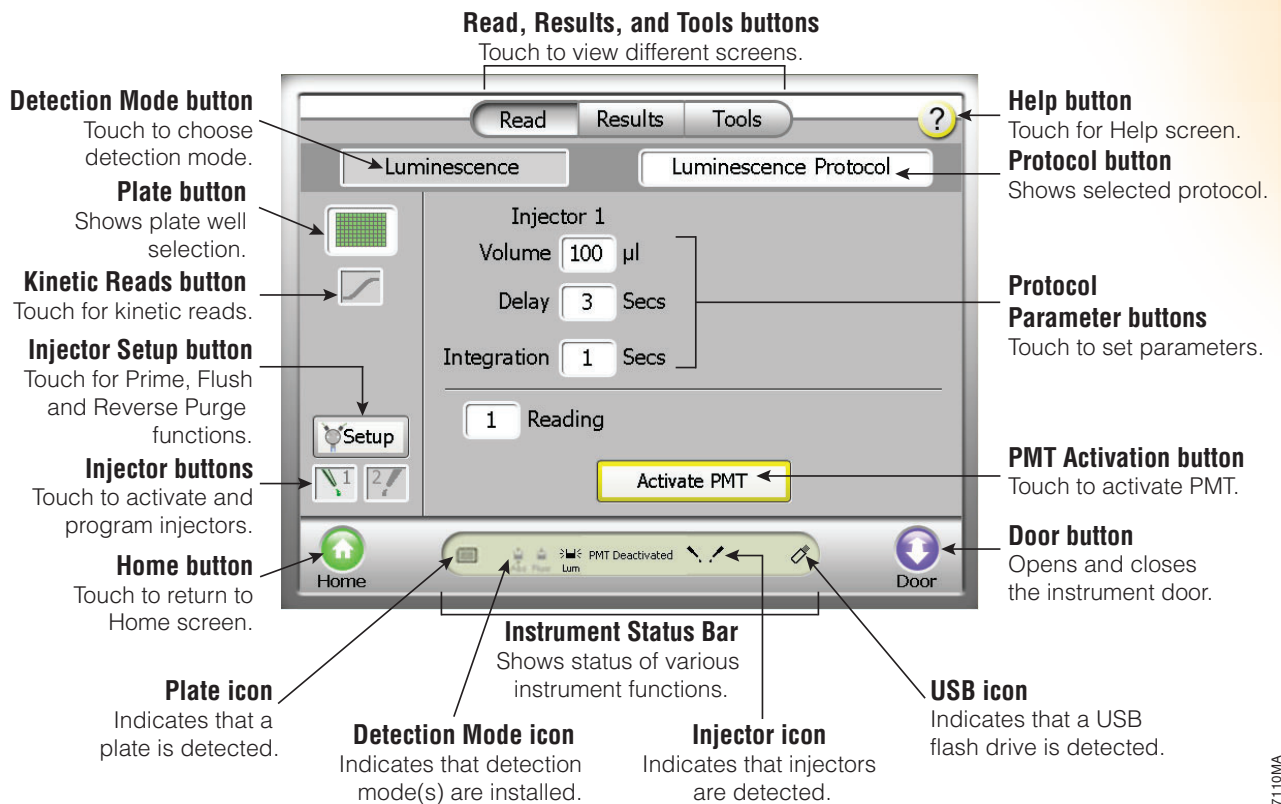
Printed in USA. Revised 5/10
Part #9FB089

GloMax®-Multi Detection System: Luminometer

INSTRUCTIONS FOR USE OF PRODUCT E7041.

Quick
PROTOCOL

“Read” Screen Map



7110MA

ORDERING/TECHNICAL INFORMATION:

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

©2008–2010 Promega Corporation. All Rights Reserved.



Promega

Printed in USA. Revised 5/10
Part #9FB089

GloMax®-Multi Detection System: Luminometer

INSTRUCTIONS FOR USE OF PRODUCT E7041.

Quick
PROTOCOL

“Results” Screen Map

Read, Results, and Tools buttons
Touch to view different screens.

USB button
Touch to transfer results file(s) to USB flash drive.

Help button
Touch for Help screen.

Files button
Touch to view the 50 most recent files. Touch Up and Down arrows to view individual file results.

Plate View button
Toggle to switch between full or partial plate view.

Door button
Opens and closes the instrument door.

Instrument Status Bar
Shows read status and file name.

Black Plate icon
Indicates that a plate read is complete.

Black USB icon
Indicates that files have been saved to USB flash drive.

Ratio button
Appears when in two-injector mode. Toggle to see results in actual reading or in ratio.

Home button
Touch to return to Home screen.

Data is displayed in this area.

Shows results file name.

	1	2	3	4	5	6
A	7.006 e1		4.993 e0			
B						
C		8.365 e0				
D						
E						
F						
G						
H						

Results:Luminescence Protocol
Tuesday, July 03, 2007 9:01 AM

Plate read complete
Results_Luminescence Protocol_7-3-2007_9-01-48 AM

7111MA

ORDERING/TECHNICAL INFORMATION:

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

©2008–2010 Promega Corporation. All Rights Reserved.



Promega

Printed in USA Revised 5/10
Part #9FB089