

GloMax[®] 20/20 Luminometer

INSTRUCTIONS FOR USE OF PRODUCTS E5311, E5321 AND E5331

Quick
PROTOCOL

Quick Start Guide

Powering the Luminometer

1. Connect the power supply to the GloMax[®] 20/20 Luminometer.
2. Turn the GloMax[®] 20/20 Luminometer "ON".
3. (optional) Connect the GloMax[®] 20/20 Luminometer to a PC using the 9-pin serial cable.
4. (optional) Install the GloMax[®] Spreadsheet Interface Software to export data from the luminometer to an Excel spreadsheet.

Choose a Protocol

5. Touch "Protocol" on the touchscreen to run a Promega protocol, create a new protocol or access a saved protocol. You may also choose to run the GloMax[™] Luminometer default protocol. The default protocol provides a simple method to measure samples without injection for an integration time of 1 second.
6. To run a Promega assay, touch "Run Promega Protocol," then select the new protocol that corresponds to your assay. The next screen will allow you to view and adjust the settings; however, the optimal settings are preset for Promega protocols. Touch the name of the setting once to highlight the setting you wish to change; then use the arrow buttons to increase or decrease the value. When you are satisfied, touch "OK" to go to the Home Screen and begin measuring your samples or "Save As" to save the protocol. Touch "Cancel" to exit from the protocol setup.
7. Touch "Create New Protocol" to set up your own protocol with or without injectors. The Measurement Parameter Screen will list the different options available. Follow the directions outlined in Step 6 to adjust the settings.

Prepare Injectors

8. Prime injectors before starting any auto-injection protocol. Touch the "Inj Func" key to access the injector functions. Select the prime option that suits your needs.
9. Each prime cycle pushes 1,000 μ l of reagent through the injection system for each injector. Place a waste container underneath the injector tips to capture the volume dispensed during priming.

For additional information see the GloMax[®] 20/20 Luminometer Technical Manual #TM276, available online at: www.promega.com.

ORDERING/TECHNICAL INFORMATION:

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




Promega

Printed in USA 3/09
Part# 9FB081

© 2006, 2009 Promega Corporation. All Rights Reserved.

GloMax[®] 20/20 Luminometer

INSTRUCTIONS FOR USE OF PRODUCTS E5311, E5321 AND E5331

Quick
PROTOCOL

Quick Start Guide

Measure Samples

10. Insert the sample into the luminometer. For injection protocols, use a microcentrifuge tube and the Microcentrifuge Tube Holder. For noninjection protocols, use a petri dish or microcentrifuge tube.
11. Close the instrument lid and touch the "Measure Luminescence" bar at the top of the touchscreen.
Note: Touching any part of the touchscreen during a protocol run will abort sample measurement.
12. The resulting data will appear on the luminometer screen. Use the arrow keys to scroll through data.
13. If you wish to use the Sample ID feature, touch the "Sample ID" key. You may name samples individually using the touchscreen keypad.
14. Touch the "Parameters" key to adjust your protocol parameters. When you are finished, touch "OK" to return to the Home Screen.
15. To save a protocol for future use, touch the "Parameters" key, then the "Save As" key. Use the keypad to name your protocol and store it for future use.

Flush Injectors

16. Clean the injection system after use. Touch the "Inj Func" key to access the injector functions. Flush the injectors that were used by touching the flush option that best suits your needs.
17. Place a waste container to capture the flushed liquid. Each flush cycle pushes 3ml of liquid through the injection system. Repeat the flush cycle several times. For best cleaning, use 1 cycle of deionized water, 1 cycle of 70% ethanol and 1 cycle of deionized water.

Touchscreen

18. Save the protocol in use. Record all of the data collected.
19. Turn the luminometer "OFF".
20. Gently wipe the touchscreen with a Kimwipes[®] dampened with 70% ethanol. Repeat if necessary.
21. Turn the luminometer "ON" to resume measurements.

ORDERING/TECHNICAL INFORMATION:

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

© 2006, 2009 Promega Corporation. All Rights Reserved.
Prices and specifications subject to change without prior notice.




Promega

Printed in USA 3/09
Part #9FB081