



# Quantifluor™ Handheld Fluorometers

Setting Up Your Instrument. It's Easy!

**QUICK START**  
GUIDE

## 1. Connections



For a complete protocol, see the Quantifluor™ Fluorometers Technical Manual TM338 provided on the CD supplied with your instrument. The Technical Manual is also available at: [www.promega.com/tbs](http://www.promega.com/tbs)



Quantifluor™ ST Fluorometer: Use only the AC Adapter provided with your instrument. Powering the Quantifluor™ ST Fluorometer with an unapproved AC adapter may damage the unit.

### Quantifluor™-P Fluorometers

1. Loosen the screw on the back of the instrument and pull the screw up to remove the battery panel.
2. Install the 4 AAA batteries into the appropriate spaces. Replace the battery panel and tighten the screw. The panel has an o-ring, which creates a watertight seal. The battery panel may be difficult to install if there is no lubrication on the o-ring. Use a silicon based grease to lightly lubricate the o-ring, if necessary.

### Quantifluor™-ST

1. Position the instrument on a flat surface.
2. Choose the appropriate AC plug for your wall outlet. Slide the desired AC plug onto the Adapter from the top until it clicks into place.
3. Plug the AC into the Quantifluor™ Fluorometer.

## 2. Power Up

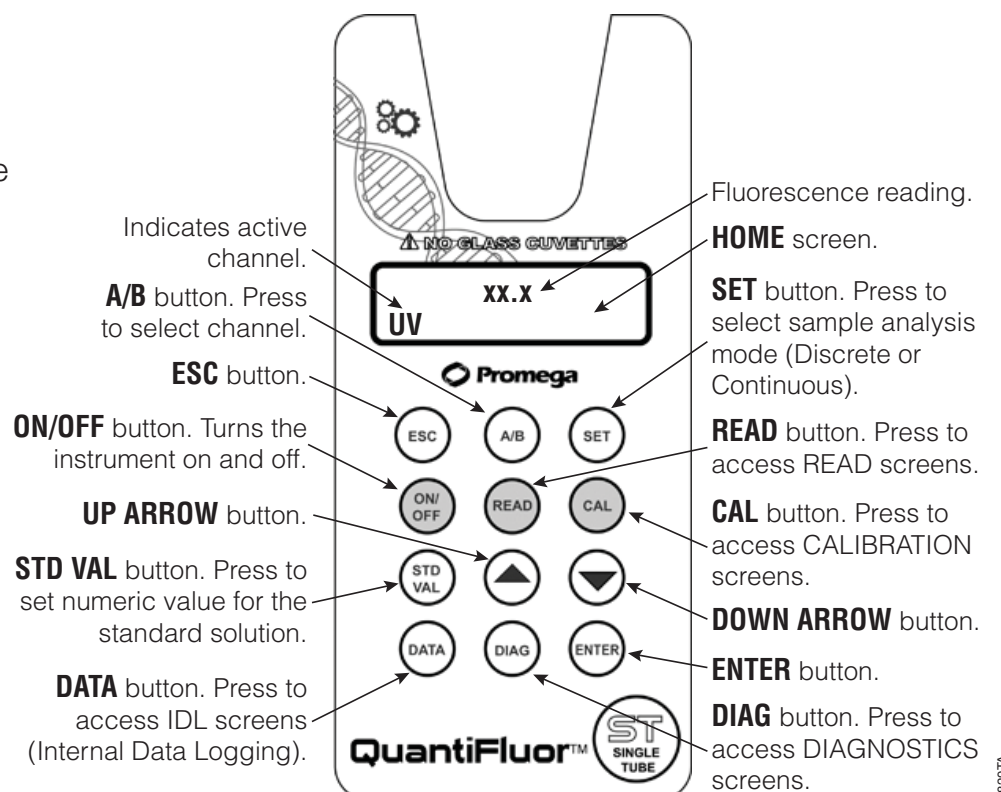
1. To turn on the Quantifluor™ Fluorometer, press the ON/OFF button. The instrument takes five seconds to warm up. After the warm-up, the Quantifluor™ Fluorometer is ready for use. Pressing the **ON/OFF** button again will turn off the unit.

**Quantifluor™-P units only:** If left idle for three minutes, the unit will automatically turn off to save battery power. A set of new batteries can last over 1,000 sample readings. If the batteries have low power or are not positioned properly, the following warning message will appear: "Batt Level < 20% !! Caution !!"

## 3. Calibration

1. Select the right channel by pressing the **A/B** button.
2. Set a standard value by pressing the **STD VAL** button and then the **UP** or **DOWN** arrows.
3. Press **ENTER** when finished.
4. Press the **CAL** button and then **ENTER**.
5. Insert a cuvette with a blank solution and press **ENTER**.
6. Insert a cuvette with the calibration standard and press **ENTER**.
7. Press **ENTER** to accept the calibration.

Note: If the Minicell Adaptor is used, be sure to insert it into the right position in the sample compartment.



Continued on other side.



# Quantifluor™ Handheld Fluorometers

**QUICK START**  
GUIDE

## 4. Sample Analysis

You can read samples in two modes: Discrete or Continuous.

### Discrete Mode

1. Press the **SET** button.
2. Select “Discrete” using the **UP** or **DOWN** arrows and press **ENTER**.
3. Insert the sample and press **READ**.

Note: For low volume samples, use the Minicell Adapter.

### Continuous Mode

1. Press the **SET** button.
2. Select “Continuous” using the **UP** or **DOWN** arrows and press **ENTER**.
3. Select a measure interval using the **UP** or **DOWN** arrows and press **ENTER**.
4. Select a number of measurements you want to obtain using the **UP** or **DOWN** arrows and press **ENTER**.
5. Press the **READ** button and the screen will display the setting selected.
6. Press **ENTER** to accept the settings.
7. Insert the sample and press **ENTER**.



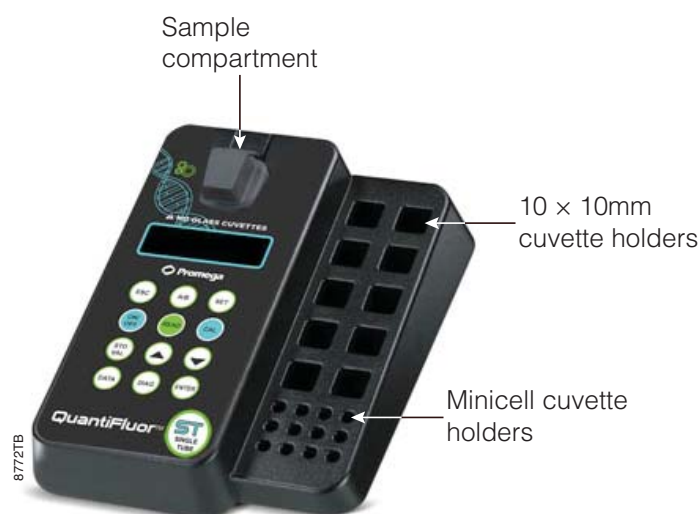
Minicell Adapter.



**Minicell Adapter in BLUE position.**  
With “BLUE” channel selected, insert the Minicell Adapter with the BLUE label facing you.



**Minicell Adapter in UV position.** With “UV” channel selected, insert the Minicell Adapter with the UV label facing you.



## 5. Installation of Spreadsheet Interface Software

1. Exit all Windows programs.
2. Insert the CD and run the Setup program.
3. The Setup Wizard will install the necessary files.
4. When Setup is complete, the “Spreadsheet Interface Software” icon will appear in the “Programs” menu.
5. Restart your computer.
6. Connect the RS-232 cable to Quantifluor™ Fluorometer and computer.
7. Open the software and select the right COM port.
8. Click “Start”.
9. Follow the directions for downloading data provided in the Quantifluor™ Fluorometers Technical Manual #TM338.



AF9FB119 0310FB119

ORDERING/TECHNICAL INFORMATION: Email: [techserv@promega.com](mailto:techserv@promega.com)  
[www.promega.com](http://www.promega.com) • Phone 608-274-4330 or 800-356-9526 • Fax 608-277-2601

©2010 Promega Corporation. All Rights Reserved. Printed in USA 03/10 18277-PM-CR. Part# FB119

  
**Promega**