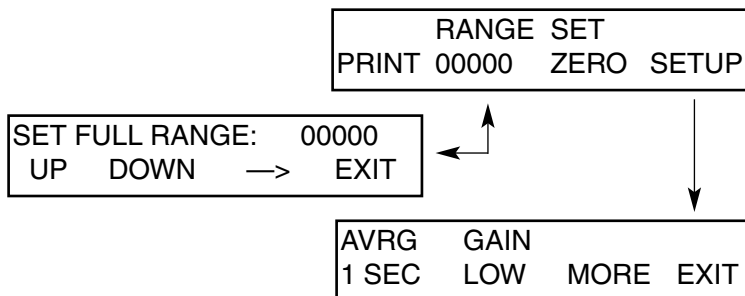




VersaFluor™ Quick Reference Guide

Note: The instrument should be allowed to warm up for at least 20 minutes before taking any readings.

1. **Selection and insertion of two optical filters.**
Insert the emission and excitation filters into the VersaFluor Fluorometer.
2. **Set the gain.**
Press the SETUP button on the main menu and set the gain to LOW, MED, or HIGH.
3. **Zero the instrument.**
Set the range to 00000 by pressing the RANGE button. Adjust the range to 00000. Press EXIT. Place the baseline sample cuvette in the cuvette chamber and press the SET ZERO button. When the blinking stops, the fluorescence display should read 0 ± 5 . If not, re-zero the instrument by pressing the SET ZERO button.
4. **Set the range.**
Set the range by placing a cuvette containing the highest concentration standard in the cuvette holder and closing the sample compartment lid. Wait approximately 10 seconds for the detector to adjust to the light conditions. Press the RANGE button. Adjust the range to the desired setting. The maximum setting is 19,999 RFU. Press EXIT.
5. **Measure the sample.**
Place a sample cuvette in the cuvette holder and close sample compartment lid. Wait approximately 10 seconds for the detector to adjust to the light conditions. Record or print the RFU value in the fluorescence display. Read all remaining samples.



VersaFluor Filter Selection Guide

Application	Excitation Filters	Catalog Number	Emission Filters	Catalog Number
DNA Quant (Hoechst 33258)	EX 360/40 (340–380 nm)	170-2420	EM 460/10 (455–465 nm)	170-2421
β -Galactosidase (MUG)	EX 360/40 (340–380 nm)	170-2420	EM 460/10 (455–465 nm)	170-2421
β -Glucuronidase (MUGluc)	EX 360/40 (340–380 nm)	170-2420	EM 460/10 (455–465 nm)	170-2421
Apopain (Z-DEVD-AFC)	EX 360/40 (340–380 nm)	170-2420	EM 520/10 (515–525 nm)	170-2424
pH indicator (BCECF)	EX 490/10 (485–495 nm)	170-2422	EM 520/10 (515–525 nm)	170-2424
Calcium indicator (Fluo-3)	EX 490/10 (485–495 nm)	170-2422	EM 520/10 (515–525 nm)	170-2424
DNA Quant (PicoGreen™)	EX 480/20 (470–490 nm)	170-2427	EM 520/10 (515–525 nm)	170-2424
Oligo Quant (OliGreen™)	EX 480/20 (470–490 nm)	170-2427	EM 520/10 (515–525 nm)	170-2424
Protein Quant (NanoOrange™)	EX 480/20 (470–490 nm)	170-2427	EM 590/10 (585–595 nm)	170-2425
Fluorescein	EX 490/10 (485–495 nm)	170-2422	EM 520/10 (515–525 nm)	170-2424
Ethidium Bromide	EX 510/10 (505–515 nm)	170-2423	EM 620/10 (615–625 nm)	170-2426
RNA Quant (RiboGreen™)	EX 490/10 (485–495 nm)	170-2422	EM 520/10 (515–525 nm)	170-2424

Helpful Hints

Use filters which are tailored to the sample being measured. Using incorrect filters will greatly degrade the accuracy of the fluorescence measurement.

- Always use calibrated pipets to insure accurate pipetting.
- Mix standards and sample completely by using a disposable transfer pipet.
- Remove any air bubbles in the cuvettes.
- Hold fluorometer cuvettes by the upper edges since the cuvettes have optically clear sides.
- Clean the cuvette sides with a lint-free tissue.
- Read all standards and samples at ambient temperature.
- While taking a reading, keep the samples in the fluorometer only. This helps reduce sample photobleaching.
- Fluorescence measurements can be made continuously. It is not necessary to zero the instrument before each measurement unless the baseline conditions change. Check the zero of the instrument with the baseline fluid (blank) when switching between gain settings.

PicoGreen, OliGreen, NanoOrange, and RiboGreen are trademarks of Molecular Probes Inc.

Bio-Rad Laboratories, 2000 Alfred Nobel Drive, Hercules, CA 94547

4100157 Rev A