



personal
DETECTION

Now, you can optimize your personal workflow. Promega instruments and reagents integrate easily.



*Fluorometers and Fluorescent
Nucleic Acid Quantitation Dyes*

QuantiFluor™

Nucleic Acid Quantitation Systems

Compact, easy to use, and cost-effective single tube fluorometers for Life Science applications.



INSTRUMENT FEATURES

QUANTIFLUOR™-ST FLUOROMETER

QUANTIFLUOR™-ST FLUOROMETER

The QuantiFluor™-ST is an affordable, sensitive fluorometer designed for quick, easy, and accurate fluorescence measurements. The QuantiFluor™-ST provides high sensitivity for fluorescence measurements and is so easy to use that you can have results within minutes. Single-point calibration saves time and the dual-channel design allows you to switch between two assays with the touch of a button.

UNIVERSAL POWER

The QuantiFluor™-ST is powered using a universal power adapter (included).

SMALL AND SENSIBLY PRICED

The QuantiFluor™-ST is smaller than a standard sized book, saving you valuable bench space. It is sensibly priced to meet your need for accurate fluorescence measurements within a limited budget.

SIMPLE

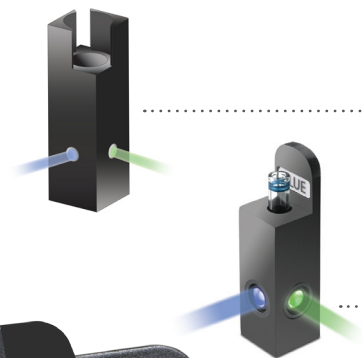
The QuantiFluor™-ST is designed for fast, accurate fluorescence measurements. It is so easy to use that you can have results within minutes without training.

DUAL CHANNEL

Dual UV and Blue Channels allow you to switch between two assays by pressing one button.

INTERNAL DATA LOGGING

An internal datalogging package with software that interfaces with Microsoft Excel can log up to 1,000 data points.



PCR TUBE ADAPTER

An optional PCR Tube Adapter allows measurement of small fluorescence sample volumes (100 µl minimum) in 0.5 mL PCR tubes without sacrificing instrument sensitivity.

MINICELL ADAPTER

An optional Minicell Adapter with two focusing lenses allows you to measure samples in low assay volumes without losing sensitivity. When used with this Minicell Adapter, the QuantiFluor™-ST can detect 50 pg of dsDNA in a 50 µL assay volume.

CALIBRATION

Single-point calibration improves accuracy while saving time.

FLUORESCENCE APPLICATIONS

UV Channel (Ex 365 nm, Em 440 - 470 nm)

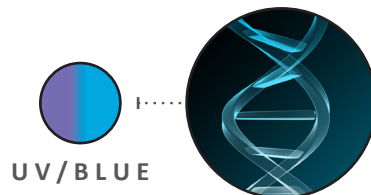
- Hoechst dye 33258 for DN A Quantitation
- 4-methyl-umbelliferone

Blue Channel (Ex 460 nm, Em 515 - 575 nm)

- DNA/RNA Quantitation
- eGFP
- Fluorescein



DUAL CHANNELS



PRODUCT

QuantiFluor™-ST Handheld Fluorometer
with UV/Blue Channels

CAT.#

E6090

INSTRUMENT FEATURES

QUANTIFLUOR™-P FLUOROMETER

QUANTIFLUOR™-P FLUOROMETER

The QuantiFluor™-P is a lightweight, handheld instrument configured for many of the fluorescent probes commonly used in nucleic acid and protein quantitation. While small in size, performance is not compromised. The QuantiFluor™-P has two dedicated channels that offer the ability to quickly analyze a sample for two different fluorescent signals.

BATTERY POWERED

The QuantiFluor™-P is powered by four size AAA batteries.

EASY TO USE

No special training is needed so you can focus on your data, not on how to use the software.

INEXPENSIVE AND COMPACT

The small, lightweight QuantiFluor™-P can be stored in a drawer or on a shelf. With its low price, you can buy several to easily manage the demands of a busy lab.

DEDICATED OPTICAL CHANNELS

Two optical channels allow you to measure two different fluorophores in the same assay.

SENSITIVE

Four logs of dynamic range.

INTERNAL DATALOGGING

An internal datalogging package with software that interfaces with Microsoft Excel can log up to 1,000 data points.



PCR TUBE ADAPTER

An optional PCR Tube Adapter allows measurement of small fluorescence sample volumes (100 µl minimum) in 0.5 mL PCR tubes without sacrificing instrument sensitivity.

MINICELL ADAPTER

The optional Minicell Adapter allows you to measure sample in low volumes between 75 - 250 µl.

FLUORESCENCE APPLICATIONS

UV/Blue Channel

- UV Channel (Ex 365 nm, Em 440 - 470 nm)**
- Hoechst dye 33258 for DNA Quantitation
 - 4-methyl-umbelliferone

Blue Channel (Ex 460 nm, Em 515 - 575 nm)

- DNA/RNA Quantitation
- eGFP
- Fluorescein

Green/Blue Channel

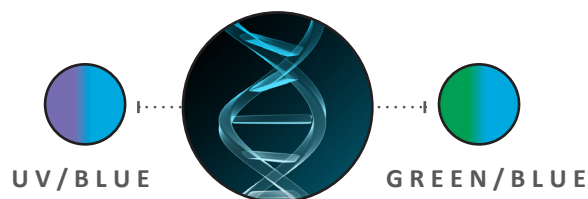
Green Channel (Ex 525 nm, Em >570 nm)

- Nucleotide or Protein Labeling
- Enzyme Activity
- Rhodamine
- Cy3

Blue Channel (Ex 460 nm, Em 515 - 575 nm)

- DNA/RNA Quantitation
- eGFP
- Fluorescein

DEDICATED OPTICAL CHANNELS



PRODUCT

CAT.

QuantiFluor™-P Handheld Fluorometer
with Green/Blue Channels

E6100

QuantiFluor™-P Handheld Fluorometer
with UV/Blue Channels

E6105

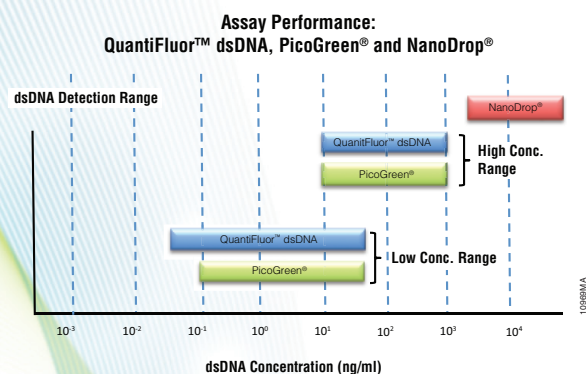
QUANTIFLUOR™ NUCLEIC ACID QUANTITATION

HIGHLY SENSITIVE, COST EFFECTIVE, AND EASY TO USE DYES FOR DSDNA, SSDNA, AND RNA QUANTITATION

These kits include all the necessary reagents to quickly set up and quantitate your nucleic acid samples. They are easy to set up on both the QuantiFluor™ and GloMax®-Multi instruments, but can be read on any fluorometer with the appropriate optical channel.

QUANTIFLUOR™ DSDNA SYSTEM

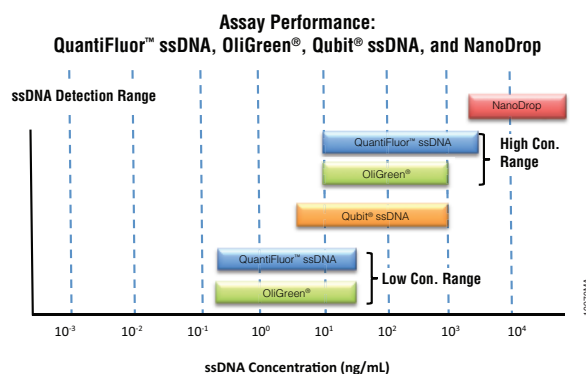
- **Specificity:** Highly specific to dsDNA, minimal binding to ssDNA, RNA, protein and interfering compounds.
- **Sensitivity:** 40,000 times more sensitive over absorbance at 260nm (NanoDrop spectrophotometer) for low-concentration samples. Performs better or equal to PicoGreen® dye, and can detect as little as 50pg/mL.



The QuantiFluor™ dsDNA System is 40,000 times more sensitive than NanoDrop® with better sensitivity than Quanti-IT™ PicoGreen®.

QUANTIFLUOR™ SSDNA SYSTEM

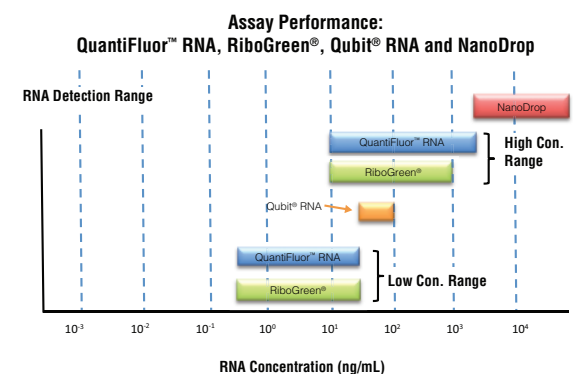
- **Increase your Sensitivity:** 10,000 times more sensitive over absorbance at 260nm over absorbance at 260nm (NanoDrop®) for those samples that are low in concentration; five times more sensitive than Qubit® ssDNA.
- **Save Precious Sample for Downstream Assays:** Less template ssDNA required than spectrophotometry.



The QuantiFluor™ ssDNA System is 10,000 times more sensitive than NanoDrop®, five times more sensitive than Qubit® ssDNA, and two times higher dynamic range than Quanti-IT™ OliGreen®.

QUANTIFLUOR™ RNA SYSTEM

- **Increase your Sensitivity:** 20,000 times more sensitive over absorbance at 260nm (NanoDrop®) for those samples that are low in concentration; 1,000 times more sensitive than Qubit® RNA.
- **Save Precious Sample for Downstream Assays:** Less template RNA required than spectrophotometry.



The QuantiFluor™ RNA System is 20,000 times more sensitive than NanoDrop®, 1,000 times more sensitive than Qubit® RNA, and two times higher dynamic range than Quanti-IT™ RiboGreen®.



PRODUCT

QuantiFluor™ dsDNA System

QuantiFluor™ ssDNA System

QuantiFluor™ RNA System

CAT.

E2670

E3190

E3110

INSTRUMENT SPECIFICATIONS

QUANTIFLUOR™ - ST

Available Detection Modes	Fluorescence
Light Source	Wavelength-matched LED
Detector	Silicon photodiode
Wavelength Selection	2 modes
Wavelengths	UV (Ex 365 nm, Em 440 - 470 nm) Blue (Ex 460 nm, Em 515 - 575 nm)
Detection Limit	1 ng/ml dsDNA with DNA Quantitation Dye, 1 ng/ml RNA using RNA Quantitation Dye, 10 ng/ml dsDNA using Hoechst Dye 33258, 100 ng/ml protein using Protein Quantitation Dye
Linear Dynamic Range	4 logs, assay dependent
Read Out	Direct concentration
Calibration	Single-point calibration
Discrete Sample Average	Sample readings averaged over 5 seconds to improve accuracy
Read Type	Discrete or Continuous
Sample Format	10 x 10 mm plastic cuvette, 0.5 mL PCR tube (with optional PCR tube adapter, 100 µl minimum), or 50 - 250 µl minicell vial (with optional adapter)
User Interface	Menu-driven keypad and LCD display
Data Output	ASCII format through a 9-pin RS-232 serial cable at 9600 baud rate (optional). Optional thermal printer available.
External PC Requirements (optional)	Windows XP SP2 or later
Computer Interface	RS-232 port
Power	Instrument: 9V DC, 0.3A AC Adapter: 100-240V AC, 50/60 Hz, 0.25A
Alarms	High blank
Dimensions	7.25" D x 5.5" W x 2.7" H (18.42 cm D x 14 cm W x 6.9 cm H)
Weight	28.8 oz (0.67 kg)
Operating Temperature	59 - 86 °F (15 - 30 °C)
Warranty	One year
Approvals	CE, UL-listed power supply

For research use only. Not for use in diagnostic procedures. QuantiFluor is a trademark of Promega Corporation. All other trademarks are the sole property of their respective owners. For the most up-to-date specifications, visit www.promega.com.

QUANTIFLUOR™ - P

Available Detection Modes	Fluorescence
Light Source	Wavelength-matched LED
Detector	Silicon photodiode
Wavelength Selection	2 modes
UV/Blue QuantiFluor™-P Wavelengths	UV (Ex 365 nm, Em 440 - 470 nm), Blue (Ex 460 nm, Em 515 - 575 nm)
Green/Blue QuantiFluor™-P Wavelengths	Green (Ex 525 nm, Em >570 nm) Blue (Ex 460 nm, Em 515 - 575 nm)
Detection Limit	1 ng/ml dsDNA with DNA Quantitation Dye, 1 ng/ml RNA using RNA Quantitation Dye, 10 ng/ml dsDNA using Hoechst Dye 33258, 100 ng/ml protein using Protein Quantitation Dye
Linear Dynamic Range	4 logs, assay dependent
Read Out	Direct concentration
Calibration	Single-point calibration
Discrete Sample Average	Sample readings averaged over 5 seconds to improve accuracy
Read Type	Discrete or Continuous
Sample Format	10 x 10 mm plastic cuvette, 0.5 mL PCR tube (with optional PCR tube adapter, 100 µl minimum), or 75 - 250 µl minicell vial (with optional adapter)
User Interface	Menu-driven keypad and LCD display
Data Output	ASCII format through a 9-pin RS-232 serial cable at 9600 baud rate (optional). Optional thermal printer available.
External PC Requirements (optional)	Windows XP SP2 or later
Computer Interface	RS-232 port
Power	4 AAA Alkaline batteries
Alarms	Low battery and high blank
Auto Shutoff	After 3 minutes of inactive use
Dimensions	7.25" D x 3.5" W x 1.75" H (18.4 cm D x 8.9 cm W x 4.45 cm H)
Weight	14 oz (0.4 kg)
Operating Temperature	59 - 86 °F (15 - 30 °C)
Warranty	One year
Approvals	CE

