

QuantiFluor® Single-Tube Fluorometers

The QuantiFluor®-ST and QuantiFluor®-P Fluorometer Instruments have been widely used laboratories worldwide for the past several years, but as of March 1, 2014, we are retiring these instrument models. They continue to offer high fluorescence sensitivity performance for customers worldwide, but we have redesigned and streamlined them into our new product, Quantus™ Fluorometer (Cat.# E6150). All QuantiFluor®-ST and QuantiFluor®-P accessories will continue to be available, as will instrument service, until March 1, 2019.

If you have any questions or would like an assessment based on your...

Expand to Read More »



Learn more about the Helix program



Prices valid for US customers only

PRODUCT	SIZE	CONC.	CATALOG #	*LIST PRICE	ORDER QTY
QuantiFluor®-ST Minicell Adapter Kit (for measuring 50–250µl of sample)	400 each	-	E6112	\$ 619.00	Add to cart
QuantiFluor®-ST Solid Standard	1 each	-	E6113	\$ 167.00	Add to cart
QuantiFluor®-ST AC Adapter Replacement	1 each	-	E6096	\$ 103.00	Add to cart
QuantiFluor®-P Minicell Adapter Kit (for measuring 75–250μl of sample)	400 each	-	E6111	\$ 301.00	Add to cart
PCR Tube Adapter, QuantiFluor® Fluorometers	1 each	-	E6101	\$ 103.00	Add to cart
Minicell Borosilicate Glass Cuvettes	400 each	-	E6091	\$ 189.00	Add to cart
10 × 10mm Square Polystyrene Cuvette (3.5ml capacity)	100 each	-	E6092	\$ 113.00	Add to cart
10 × 10mm Square Methacrylate Cuvette (3.5ml capacity)	100 each	-	E6093	\$ 111.00	Add to cart
Thermal Serial Printer and Universal Power Cable	1 each	-	E2821	\$ 670.00	Add to cart
Thermal Printer Paper	1 each	-	E2851	\$ 26.00	Add to cart
0.5ml PCR Tubes Add to Helix	50 pack	-	E4941	\$ 30.00	Add to cart
0.5ml PCR Tubes	200 pack	-	E4942	\$ 55.00	Add to cart
QuantiFluor® Service Agreement	1 each	-	SA3060	Please Enquire	

You Also May Be Interested In



QuantiFluor® dsDNA System

QuantiFluor® RNA System

Recombinant RNasin® Ribonuclease Inhibitor

GoScript™ Reverse Transcription System



Figure 1. QuantiFluor®-ST and QuantiFluor®-P Handheld Fluorometers.

Articles

Measuring the QuantiFluor™ dsDNA System Using the QuantiFluor™-ST Fluorometer

Measuring the QuantiFluor™ RNA System Using the QuantiFluor™-ST Fluorometer

Related Products

Agarose, LE, Analytical Grade

Promega Barrier Tips

Nuclease-Free Water

Diamond™ Nucleic Acid Dye	
RNA Markers	
FAQs	
QuantiFluor™ Fluorometers FAQ	
Videos	
QuantiFluor™ dsDNA System Video	
QuantiFluor™ RNA System Video	
Guides	
Fluorophores Guide	
Application Notes	
Application Notes: QuantiFluor® ST and P Fluorometers	
E6112, E6113, E6096, E6111, E6101, E6091, E6092, E6093, E2821, E2851, E49 in Diagnostic Procedures.	941, E4942, SA3060 For Research Use Only. Not for Use
View All Reviews »	
Review This Product »	
by	
View All Reviews » Review This Product »	
by	

View All Reviews »			
Review This Product »			
by			
View All Reviews »			
Review This Product »			
by			
Jy			
View All Reviews »			
Review This Product »			
by			
View All Reviews »			
Review This Product »			
by			

View All Reviews » Review This Product »		
by		
View All Reviews » Review This Product »		
by		
View All Reviews » Review This Product »		
by		
View All Reviews » Review This Product »		
by		

View All Reviews »	
VIEW All INEVIEWS //	
Review This Product »	
Review IIIIS Floudct »	
by	
Бу	
Mary All Devieus	
View All Reviews »	
Review This Product »	
Review Inis Product »	
by	
by	
by	
by	
by View All Reviews »	
by View All Reviews »	
by	
by View All Reviews »	
View All Reviews » Review This Product »	
View All Reviews » Review This Product »	
by View All Reviews »	
View All Reviews » Review This Product »	
View All Reviews » Review This Product »	
View All Reviews » Review This Product »	
View All Reviews » Review This Product »	
View All Reviews » Review This Product »	
View All Reviews » Review This Product »	
View All Reviews » Review This Product »	