



GoTaq® Real-Time qPCR and RT-qPCR Systems for Dye-Based Detection



The **GoTaq® qPCR Master Mix** is a ready-to-use 2X master mix for use in real-time quantitative PCR (qPCR and RT-qPCR). The system contains BRYT Green® dye, a novel fluorescent DNA-binding dye with minimal PCR inhibition for maximum PCR efficiency and greater fluorescence enhancement upon binding to double-stranded DNA (dsDNA) than SYBR® Green I. Containing the GoTaq® Hot Start Polymerase, optimized buffer and proprietary dye, the GoTaq® qPCR Master Mix provides robust real-time PCR with earlier quantification cycle values and broad range detection for increased reliability, reproducibility and s...

[Expand to Read More »](#)



[Learn more about the Helix program](#)



Prices valid for US customers only

PRODUCT	SIZE	CONC.	CATALOG #	*LIST PRICE	QTY	
GoTaq® qPCR Master Mix	200 × 50µl reactions	-	A6001	\$ 345.00	<input type="text"/>	Add to cart
Protocol MSDS Components Add to Helix						
GoTaq® qPCR Master Mix	1,000 × 50µl reactions	-	A6002	\$ 1,383.00	<input type="text"/>	Add to cart
Protocol MSDS Components Add to Helix						
GoTaq® 2-Step RT-qPCR System	50 × 20µl RT reactions + 200 × 50µl qPCR reactions	-	A6010	\$ 498.00	<input type="text"/>	Add to cart
Protocol MSDS Components Add to Helix						
GoTaq® 1-Step RT-qPCR System	200 × 50µl reactions	-	A6020	\$ 605.00	<input type="text"/>	Add to cart
Protocol MSDS Components Add to Helix						

Available Separately

PRODUCT	SIZE	CONC.	CATALOG #	*LIST PRICE	QTY	
CXR Reference Dye	100µl	-	C5411	\$ 23.00	<input type="text"/>	Add to cart
MSDS Components Add to Helix						

Storage Conditions

Upon arrival, store all components at -30 to -10°C , protected from light. For immediate use, components may be stored at $2-8^{\circ}\text{C}$, protected from light, for up to 3 months.

For product intended use please see Patents & Disclaimers tab.

You Also May Be Interested In



[GoTaq® Hot Start Polymerase](#)

[GoScript™ Reverse Transcription System](#)

[Maxwell® 16 LEV simplyRNA Purification Kits](#)

[QuantiFluor® RNA System](#)