

qPCR Master Mixes

EvaGreen® dye-based or probe mixes

Forget-Me-Not™ EvaGreen® qPCR Master Mix

Forget-Me-Not™ qPCR Master Mix is a hot-start EvaGreen® dye-based master mix for use in real-time PCR applications and DNA melt curve analysis. Forget-Me-Not™ features a unique combination of a master mix containing a low concentration of blue dye, plus a DNA template buffer containing a higher concentration of blue dye. When you add the 2X Forget-Me-Not™ Master Mix to your reaction tube, it appears light blue. Then, when you add template containing Forget-Me-Not™ Template Buffer to the reaction, the color turns dark blue. Forget-Me-Not™ allows you to see at a glance whether you forgot to add master mix or template to any of your reactions, so you can catch pipetting mistakes and avoid wasting time, reagents, and your precious DNA samples. The 2X Forget-Me-Not™ Master Mix also can be used without the Template Buffer if you prefer.

Unrivaled Real-time PCR Performance

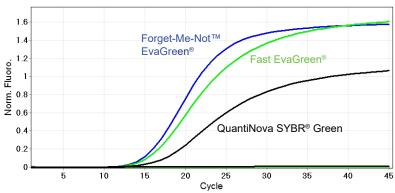


Figure 1. Real-time PCR data comparing Forget-Me-Not™ (blue line) with Biotium's Fast EvaGreen® (green line) and Qiagen's QuantiNova® SYBR® Green (black line) master mixes. Amplification curves on linear scale. EvaGreen® dye-based master mixes yield higher signal compared to the SYBR® Greenbased mix. Forget-Me-Not™ performs as well or better than the other master mixes.

Forget-Me-Not™ EvaGreen® Advantages

- Unbeatable prices! As low as 20¢/rxn
- Contains EvaGreen[®] dye for superior qPCR performance and melt curve analysis
- Two-color tracking to streamline sample set-up and minimize errors
- PCR products can be directly visualized on a gel
- Hot-start Cheetah[™] Taq is ready after just 2 minutes at 95°C



Two Color Tracking to Minimize Errors



Figure 2. PCR tubes containing Forget-Me-Not™ qPCR Master Mix (1X) on left and Forget-Me-Not qPCR Master Mix (1X) plus DNA template in Template Buffer on the right.

EvaGreen® PCR dye

- Bright and sensitive DNA-binding dye
- Detected in the SYBR® Green channel
- Less PCR inhibition than SYBR[®] Green
- Directly visualize PCR products in gels
- Environmentally friendly
- The only qPCR dye used in droplet digital PCR (ddPCR)
- The best dye for high resolution melt curve analysis (HRM)



Forget-Me-Not™ Universal Probe Master Mix

Forget-Me-Not™ Universal Probe Master Mix is a high-performance product for fluorescent probe-based PCR applications, including quantitation and SNP genotyping. This kit is suitable for all fluorescent probe-based technologies, including hydrolysis probes (such as TaqMan® and dual-labeled BHQ® probes) and displacement probes (like Molecular Beacons). Forget-Me-Not™ Universal Probe Master Mix shows excellent concordance of results in singleplex and multiplex reactions, has broad instrument compatibility, and can be used in both standard and fast protocols.

Forget-Me-Not™ Universal Probe Master Mix comes with an optional blue tracking buffer containing an inert blue dye. You have the choice of adding Tracking Buffer to the master mix, to the DNA template, or not to use the tracking buffer in your reactions.

Ideal for multiplex probe reactions

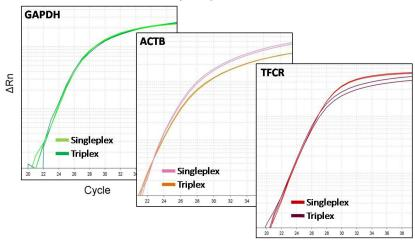


Figure 3. Singleplex and triplex PCR of GAPDH, ACTB and TFCR from human genomic DNA. Reactions contained 20 ng of human genomic DNA, 400 nM of each primer, and 200 nM of each hydrolysis probe (GAPDH-FAM/BHQ®-1, ACTB-CAL Fluor® Orange 560/BHQ®-1, TFCR-CAL Fluor® Red 610/ BHQ®-2).

Cheetah™ Taq

- Proprietary chemically-modified hot-start Taq polymerase
- · Completely inhibited on the bench, but fully active after just 2 minutes of heating
- Ideal for fast PCR reactions
- Used in all of Biotium's qPCR Master Mixes

Ordering Information

Cat. #	Product name	Unit size
31041-T	Forget-Me-Not™ EvaGreen® qPCR Master Mix	1 mL, trial size
31041-1		5 x 1 mL
31041-20mL		2 x 10 mL
31042-T	Forget-Me-Not™ EvaGreen® qPCR Master Mix with ROX	1 mL, trial size
31042-1		5 x 1 mL
31042-20mL		2 x 10 mL
31043-T	Forget-Me-Not™ Universal Probe Master Mix	1 mL, trial size
31043-1		5 x 1 mL
31044-T	Forget-Me-Not™ Universal Probe Master Mix with ROX	1 mL, trial size
31044-1		5 x 1 mL
31000-T	EvaGreen® Dye, 20X in Water	1 mL, trial size
31000		5 x 1 mL
31019	EvaGreen® Dye, 2000X in DMSO	50 uL
29050	Cheetah™ Taq	500 U
29052	ROX Passive Reference Dye, 25 uM in TE	5 x 1 mL
40052	dNTP Set, 100 mM Each	Set of 1 mL each
40054	dNTP Mix, 10 mM Each	5 x 1 mL

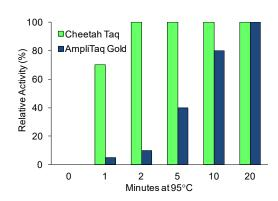


Figure 4. Cheetah™ Taq requires only 2 minutes of hot-start for full recovery of activity. Comparison of hot-start recovery of polymerase activity for Cheetah™ Taq and AmpliTaq Gold® following incubation at 95°C in 50 mM pH 8.0 Tris.

