



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

Forget-Me-Not™ Universal Probe qPCR Master Mix



A high-performance product for fluorescent probe-based PCR applications, including quantitation and SNP genotyping.

Product attributes

Reference dye	No ROX, With Separate ROX
Detection method/readout	PCR/qPCR

Product Description

2X Forget-Me-Not™ Universal qPCR Probe Master Mix is a high-performance product for fluorescent probe-based PCR applications, including quantitation and SNP genotyping.

- Suitable for all fluorescent probe types, including TaqMan®, BHQ®, and molecular beacons
- Sensitive, robust, and reproducible performance at a fraction of the cost of other mixes
- For singleplex or multiplex reactions
- Tracking Buffer helps you catch pipetting mistakes, saving time and precious samples
- Cheetah™ HotStart Taq for fast cycling protocols
- With or without separate ROX for use with any qPCR instrument

The master mix contains [Cheetah™ HotStart Taq DNA Polymerase](#) and dNTPs in a buffer optimized for high qPCR sensitivity and multiplex reactions. Only primers, probe, and template need to be added.

Cheetah Taq is covered under US and international patents. TaqMan is a registered trademark of Roche Diagnostics. CAL Fluor and BHQ are registered trademarks of LGC Biosearch Technologies.

This datasheet was generated on June 3, 2026 at 06:49:26 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/forget-me-not-universal-probe-master-mix/>