



Glowing products for science

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

Fast Probe Master Mix

A qPCR master mix containing Cheetah™ Taq hotstart DNA polymerase, suitable for probe-based real-time PCR applications.



Product attributes

Detection method/readout	PCR/qPCR
Reference dye	No ROX, With ROX
Storage Conditions	Store at -10 to -35 °C

Product Description

Fast Probe Master Mix is designed for fluorescent probe-based PCR applications. It is compatible with probes such as a TaqMan®, BHQ®, and Molecular Beacon. However for most PCR applications we recommend the [Forget-Me-Not™ Universal Probe Master Mix](#), for superior activity.

- 2X probe-based qPCR master mix for fast cycling protocols
- Suitable for all fluorescent probe types, including TaqMan®, BHQ®, and Molecular Beacon
- Contains Cheetah™ Taq hotstart DNA polymerase

Fast Probe Master Mix is available with or without ROX reference dye, for use with high ROX, low ROX, or no ROX instruments.

Cheetah Taq is covered under US and international patents. TaqMan is a registered trademark of Roche Diagnostics.

This datasheet was generated on June 23, 2026 at 06:29:24 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/fast-probe-master-mix/>