Glowing products for science

## **Fast Probe Master Mix**

A qPCR master mix containing Cheetah<sup>TM</sup> 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

Call us: 800-304-5357 Email: btinfo@biotium.com

## **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 November 4, 2025 at 12:14:43 PM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/fast-probe-master-mix/">https://biotium.com/product/fast-probe-master-mix/</a>