Fast Plus EvaGreen® qPCR Master Mix

A qPCR master mix containing Evagreen® qPCR dye and CheetahTM Taq hotstart DNA polymerase, suitable for qPCR using a fast cycling protocol.



Product attributes

 Detection method/readout
 PCR/qPCR, HRM®

 Reference dye
 No ROX, Low ROX, High ROX

 Storage Conditions
 Store at -10 to -35 °C

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

Product Description

Fast Plus EvaGreen® qPCR Master Mix contains Evagreen® qPCR dye and Cheetah™ Taq hotstart DNA polymerase. It is suitable for qPCR using a fast cycling protocol. However for most PCR applications we recommend the Forget-Me-Not™ series of master mixes, for superior activity.

- 2X qPCR master mix for fast cycling protocols
- EvaGreen® dye-based
- Contains Cheetah™ Taq hotstart DNA polymerase

Fast Plus Evagreen® qPCR Master Mix is a 2X master mix, and is supplied with a separate vial of ROX reference dye, for use with high ROX, low ROX, or no ROX instruments.

EvaGreen Dye and Cheetah Taq are covered under US and international patents

References

Mao, et al. Characterization of EvaGreen Dye and the implication of its physicochemical properties for gPCR applications. *BMC Biotechnology* **7**, 76 (2007).

Download a list of Selected EvaGreen References.

This datasheet was generated on November 7, 2025 at 07:20:09 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/fast-plus-evagreen-apcr-master-mix/