## Fast EvaGreen® qPCR Master Mix

A qPCR master mix containing Evagreen® qPCR dye and Cheetah<sup>TM</sup> Taq hotstart DNA polymerase, suitable for qPCR using a fast cycling protocol.



## **Product attributes**

Detection method/readout PCR/qPCR, HRM®

Storage Conditions Store at -10 to -35 °C

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

## **Product Description**

Fast 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 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 Tag are covered under US and international patents

## References

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

Download a list of Selected EvaGreen References.

This datasheet was generated on November 18, 2025 at 07:21:05 AM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/fast-evagreena-apcr-master-mix-200-rxn/">https://biotium.com/product/fast-evagreena-apcr-master-mix-200-rxn/</a>