



Glowing products for science

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

## Fast Plus EvaGreen® qPCR Master Mix

A qPCR master mix containing Evagreen® qPCR dye and Cheetah™ 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

## 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 qPCR applications](#). *BMC Biotechnology* 7, 76 (2007).

Download a list of [Selected EvaGreen References](#).

This datasheet was generated on June 4, 2026 at 08:53:01 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/fast-plus-evagreen-qpcr-master-mix/>