



Which real-time instruments are compatible with GoTaq® qPCR Master Mix?

GoTaq® qPCR Master Mix (Cat.# A6001) can be used with any real-time instrument capable of detecting SYBR® Green I or FAM™ dye. The GoTaq® qPCR Master Mix is optimized for use on the Applied Biosystems 7500 and 7500 FAST Real-Time PCR Systems and is compatible with many other instruments. Some real-time instruments require a higher level of reference dye than what is included in the GoTaq® qPCR Master Mix. A separate tube of 100X CXR Reference Dye is provided with the GoTaq® qPCR Master Mix for use with these instruments.

The following real-time instruments have been tested and found to be compatible with the GoTaq® qPCR Master Mix:

Instruments that do not require addition of 100X CXR Reference Dye

- Applied Biosystems 7500 and 7500 FAST Real-Time PCR Systems
- Stratagene Mx3000P® and Mx3005P® qPCR Systems
- Roche LightCycler® 480 and LightCycler® 1.5 and 2.0
- BioRad Chromo4™ Real-Time Detector
- Eppendorf Mastercycler® ep *realplex4* and *realplex4 S*
- BioRad iCycler iQ™5 and MyCycler™ **Note:** May require addition of fluorescein reference dye(1)

Instruments that require addition of 100X CXR Reference Dye to 1X per reaction

- Applied Biosystems 7000 Sequence Detection System
- Applied Biosystems 7300 Real-Time PCR System
- Applied Biosystems 7700 Sequence Detection System
- Applied Biosystems 7900HT Real-Time PCR System
- Applied Biosystems StepOne™ and StepOnePlus™ Real-Time PCR Systems

Reference

1. Reece, K.L. *et al.* (2009) [Introducing GoTaq® qPCR Master Mix: The bright choice for dye-based qPCR. Promega PubHub.](#)

GoTaq is a registered trademark of Promega Corporation.

FAM, StepOne and StepOnePlus are trademarks of Applied Biosystems Corporation. iCycler is a registered trademark and Chromo4, iQ and MyCycler are trademarks of Bio-Rad Laboratories, Inc. LightCycler is a registered trademark of Roche Diagnostics, GmbH. Mastercycler is a registered trademark of Eppendorf-Netheler-Hinz GmbH. Mx3000P and Mx3005P are registered trademarks of Stratagene. SYBR is a registered trademark of Molecular Probes, Inc.

Products may be covered by pending or issued patents or may have certain limitations. Please visit our web site for [more information](#).

All prices and specifications are subject to change without prior notice.

Product claims are subject to change. Please contact Promega Technical Services or access the Promega online catalog for the most up-to-date information on Promega products.