

Methylamp Hot Taq Probe qPCR Mix (no ROX)

(Catalog No. R12025)

Description

Methylamp Hot Taq Probe qPCR Mix (no ROX) is optimized for real-time quantitative PCR assays and contains all the components necessary to perform qPCR, with the exception of template, primers, and probe. The qPCR Mix contains optimized components and Methylamp Hot Taq DNA polymerase supplied in a proprietary reaction buffer that enables detection of low copy number targets. Methylamp Hot Taq DNA polymerase is activated by a 15 min incubation step at 95°C. This prevents extension of non-specifically annealed primers and primer-dimers formed at low temperatures during qPCR setup.

Methylamp Hot Taq Probe qPCR Mix (no ROX) is compatible for use with any probe system and qPCR cyclers that do not require reference dye, including Bio-Rad iQ™ 5, Opticon™, Opticon™ 2; Chromo 4™, MiniOpticon, CFX96, CFX384; Cepheid SmartCycler™; Corbett Rotor-Gene™ 3000, Rotor-Gene™ 6000; Eppendorf Mastercycler® ep realplex; Roche LightCycler® 480 and Techne Quantica™.

Composition

- Methylamp Hot Taq DNA polymerase
- 5 x qPCR buffer P
- 15 mM MgCl₂: 1 x PCR solution – 3 mM MgCl₂
- dNTPs, including dTTP to improve reaction sensitivity and efficiency compared to dUTP
- No ROX dye
- 1 ml of mix is sufficient for 250 reactions

Applications

- Detection and quantification of DNA and cDNA targets including DNA from bisulfite treatment, ChIP or MeDIP
- Profiling gene expression
- High Resolution Melt (HRM)
- Analysis Microbial detection
- Viral load determination

Recommended PCR reaction mix:

Component	Volume	Final Conc.
Methylamp Hot Taq Probe qPCR Mix	4 µl	1x
Primer Forward (10 pmol/µl)	0.4-0.8 µl	200-400 nM
Primer Reverse (10 pmol/µl)	0.4-0.8 µl	200-400 nM
Probe	1 µl	100-250 nM
DNA Template	1-5 µl	1-50 ng/µl
H ₂ O PCR Grade	Up to 20 µl	
Total	20 µl	

Recommended PCR cycles:

Cycle step	Temp.	Time	Cycles
Initial Denaturation	95°C	15 min	1
Denaturation	95°C	15 s	40
Annealing/ Elongation	60°C	60 s	40

IMPORTANT: To activate the polymerase, include an incubation step **at 95°C for 15 minutes** at the beginning of the qPCR cycle.

Storage Conditions

Routine storage: -20°C. Shipping and temporary storage for up to 1 month at room temperature has no detrimental effects on the quality of Methylamp Hot Taq Probe qPCR Mix (no ROX).

Ordering Information

Products

Methylamp Hot Taq Probe qPCR Mix (no ROX)

Size

1 ml

Cat. No.

R12025-1

This product is for research purposes only. Not intended for use in diagnostic procedures.