

Methylamp Hot Taq Probe qPCR Mix (Capillary)

(Catalog No. R12026)

Description

Methylamp Hot Taq Probe qPCR Mix (Capillary) is optimized for real-time quantitative PCR assays. The ready-to-use mix includes Methylamp Hot Taq DNA polymerase, ultrapure dNTPs and MgCl₂. Only template, primers, probe and water need to be added. 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 (Capillary) can be used with LightCycler® 1.x and LightCycler® 2.0 (Roche Applied Sciences).

Composition

- Methylamp Hot Taq DNA polymerase
- 5x qPCR Buffer P
- 15 mM MgCl₂: 1 x PCR solution – 3 mM MgCl₂
- dNTPs, including dTTP to improve reaction sensitivity and efficiency
- BSA
- 1 ml of mix is sufficient for 250 reactions

Applications

- Detection and quantification of DNA and cDNA targets
- Profiling gene expression
- Microbial detection
- Viral load determination

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 (Capillary).

Ordering Information

Products

Methylamp Hot Taq Probe qPCR Mix (Capillary)

Size

1 ml

Cat. No.

R12026-1

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	15-20 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.

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