

## Methylamp Taq 5x PCR Mix 12.5 mM

(Catalog No. R12021)

### Description

Methylamp Taq 5x PCR Mix is a premixed ready-to-use solution containing all reagents required for PCR (except template, primers and water).

### Composition

- Methylamp Taq DNA polymerase
- 5 x Reaction Buffer 1: 0.4 M Tris-HCl, 0.1 M (NH<sub>4</sub>)<sub>2</sub>SO<sub>4</sub>, 0.1% w/v Tween-20
- 12.5 mM MgCl<sub>2</sub>: 1x PCR solution – 2.5 mM MgCl<sub>2</sub>
- 1 mM dNTPs of each: 1x PCR solution – 200 μM dATP, 200 μM dCTP, 200 μM dGTP and 200 μM dTTP
- 1 ml of mix is sufficient for 250 reactions

### Applications

- Suited for a wide range of PCR assays
- TA cloning

### Recommendations

We recommend using Methylamp 5x PCR Mix in any PCR applications that will be visualized by agarose gel electrophoresis and ethidium bromide staining. We do not recommend using the 5x PCR Mix for PCR reactions where detergent free buffer system is required (detergent free 5x PCR Mix available upon request).

### Safety Warnings and Precautions

This product is designed for research purposes and in vitro use only. According to common laboratory safety practice, it is recommended to wear protective clothing, gloves and safety glasses. Please refer to [www.epigentek.com](http://www.epigentek.com) for Material Safety Data Sheet of the product. Some applications this product is used in may require a license which is not provided by the purchase of this product. Users should obtain the license if required.

### Storage Conditions

Routine storage: -20°C. Shipping and temporary storage for up to 1 month at room temperature or storage for up to 6 months at 2-8°C has no detrimental effects on the quality of Methylamp Taq 5x PCR Mix.

### Ordering Information

#### Products

Methylamp Taq 5x PCR Mix 12.5 mM

#### Size

1 ml

#### Cat. No.

R12021-1

### Recommended PCR reaction mix:

Component	Volume	Final Conc.
Methylamp Taq 5x PCR Mix	4 μl	1x
Forward Primer (10 pmol/μl)	0.2-0.6 μl	0.1-0.3 μM
Reverse Primer (10 pmol/μl)	0.2-0.6 μl	0.1-0.3 μM
Template DNA	x μl	5-50 ng/μl
Add H <sub>2</sub> O	Up to 20 μl	

### Recommended PCR cycles:

Cycle step	Temp.	Time	Cycles
Initial Denaturation	95°C	3-5 min	1
Denaturation	95°C	20-40 s	25-30
Annealing	54-66°C	30-60 s	25-30
Elongation	72°C	40 s-4 min	25-30
Final Elongation	72°C	5-10 min	

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