

## Methylamp Taq 5x PCR Mix Ready-to-Load 12.5 mM

(Catalog No. R12022)

### Description

Methylamp Taq 5x PCR Mix Ready-to-Load is a premixed ready-to-use solution containing: all reagents required for PCR (except template, primers and water), additional compound needed for direct loading onto agarose gel and two tracking dyes (blue and yellow) that allow to monitor progress during electrophoresis.

### Composition

- Methylamp Taq DNA polymerase
- 5x 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
- Blue dye: Migration equivalent to 3.5-4.5 kb DNA fragment
- Yellow dye: Migration rate in excess of primers in 1% agarose gel: <35-45 bp
- Compound that increases sample density for direct loading
- 1 ml of mix is sufficient for 250 reactions

### Applications

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

### Recommended PCR reaction mix:

Component	Volume	Final Conc.
Methylamp Taq 5x PCR Mix Ready-to-Load	10 μl	1x
Forward Primer (10 pmol/μl)	0.5-1.5 μl	0.1-0.3 μM
Reverse Primer (10 pmol/μl)	0.5-1.5 μl	0.1-0.3 μM
Template DNA	x μl	5-50 ng/μl
Add H <sub>2</sub> O	Up to 50 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	

### Recommendations

We recommend using Methylamp Taq 5x PCR Mix Ready-to-Load in any PCR applications that will be visualized by agarose gel electrophoresis and ethidium bromide staining. We do not recommend using the Methylamp Taq 5x PCR Mix Ready-to-Load for PCR reactions where detergent free buffer system is required (detergent free 5x PCR Mix available upon request). Methylamp Taq 5x PCR Mix Ready-to-Load is not recommended for use in applications where spectrophotometric measurements (absorbance or fluorescence) are necessary because yellow and blue dyes can interfere with these applications.

### 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 Ready-to-Load.

### Ordering Information

#### Products

Methylamp Taq 5x PCR Mix Ready-to-Load 12.5 mM

#### Size

1 ml

#### Cat. No.

R12022-1

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