

## T7 RNA Polymerase

**Cat. No. 18033-019**

**Conc.: 50 U/μl**

**Size: 2,500 units**

**Store at -20°C (not frost-free).**

### Description:

T7 RNA Polymerase is a DNA-dependent RNA polymerase which has been isolated from *E. coli* expressing the T7 RNA polymerase gene on a plasmid (1). The enzyme has an extremely high specificity for T7 promoter sequences (2) and will synthesize large quantities of RNA from a DNA fragment inserted downstream from a promoter. A strong class III promoter (3) has been used to construct various cloning vectors, and inserts into the multiple cloning site of these vectors can be transcribed to generate discrete RNA's.

### Components:

18033-019 T7 RNA Polymerase  
Y90108 5X T3/T7 Buffer  
Y00147 0.1 M DTT

### Unit Definition:

One unit incorporates 1 nmol of labeled nucleotide into acid-precipitable material in 1 hour at 37°C.

### Storage Buffer:

20 mM Tris-HCl (pH 7.5)  
0.1 M NaCl  
0.1 mM EDTA  
1 mM DTT  
50% (v/v) glycerol  
0.01% (w/v) Triton® X-100

### 5X T3/T7 Buffer:

0.2 M Tris-HCl (pH 8.0)  
40 mM MgCl<sub>2</sub>  
10 mM spermidine-(HCl)<sub>3</sub>  
125 mM NaCl  
Refer to Functional Assay  
Conditions on reverse side for  
further details.

Doc. Rev.: 050602

This product is distributed for laboratory research only. CAUTION: Not for diagnostic use. The safety and efficacy of this product in diagnostic or other clinical uses has not been established.

Quality Control:

This product has passed the following quality control assays: functional absence of exonuclease, endo-ribonuclease and DNA nicking activities; performance in a transcription reaction.

The enclosed buffers were assayed with the enzyme and met quality control specifications.

Functional Assay Conditions:

2  $\mu$ l 5X T3/T7 Buffer  
70  $\mu$ M [ $\alpha$ -<sup>32</sup>P]UTP (280  $\mu$ Ci of 400 Ci/mmol)  
0.4 mM each ATP, CTP, GTP  
5 mM DTT  
0.1  $\mu$ g linearized template DNA  
50 units T7 RNA Polymerase  
Reaction Volume: 10  $\mu$ l  
Incubation: 10 minutes at 37°C

NOTE: The reaction is not set up on ice due to potential precipitation of DNA in the presence of spermidine.

References:

1. Davanloo, P., Rosenberg, A. H., Dunn, J. J., and Studier, F. W. (1984) *Proc. Natl. Acad. Sci. USA* 81, 2035.
2. Chamberlin, M., McGrath, J., and Waskell, L. (1970) *Nature* 228, 227.
3. Studier, F. W., and Dunn, J. J. (1983) *Cold Spring Harbor Symposia on Quantitative Biology XLVII*, 999.

Triton® is a registered trademark of Rohm & Haas, Co.