thermo scientific

T7 RNA Polymerase

Catalog Number EP0111, EP0112, EP0113

Pub. No. MAN0016017 Rev. B.00

WARNING! Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. Safety Data Sheets (SDSs) are available from thermofisher.com/support.

Product description

Bacteriophage T7 RNA Polymerase is a DNAdependent RNA polymerase with strict specificity for its respective double-stranded promoter. The enzyme catalyzes the $5' \rightarrow 3'$ synthesis of RNA on either single-stranded DNA or double-stranded DNA downstream from the promoter. T7 RNA Polymerase accepts modified nucleotides (e.g., biotin-, digoxigenin-, fluorescein-labeled nucleotides) as substrates for RNA synthesis.

Contents and storage

Cat. No.	Components	Amount	Source	Molecular Weight	Storage
EP0111	T7 RNA Polymerase	5000 U, 20 U/µL		99 kDa monomer	-25 °C to -15 °C
	5X Transcription Buffer	1.25 mL	<i>E.coli</i> cells with a cloned gene encoding this		
EP0112	T7 RNA Polymerase	5 x 5000 U, 20 U/µL			
	5X Transcription Buffer	5 x 1.25 mL			
EP0113	T7 RNA Polymerase	25000 U, 200 U/µL HC	enzyme		
EFUIIS	5X Transcription Buffer	5 x 1.25 mL			

BSA included

Applications

Synthesis of unlabeled and labeled RNA that can be used:

- for hybridization (1), in vitro RNA translation (2);
- as aRNA (3), siRNA (4), substrate in RNase protection assays (5), template for genomic DNA sequencing (6);
- in studies of RNA secondary structure and RNA-protein interactions (7), RNA splicing (8).

Definition of Activity Unit

One unit of the enzyme incorporates 1 nmol of AMP into a polynucleotide fraction in 60 minutes at 37 °C.

Storage Buffer

The enzyme is supplied in: 50 mM Tris-HCI (pH 8.0), 150 mM NaCl, 5 mM DTT, 0.1 mg/mL BSA, 0.03% (v/v) ELUGENT Detergent and 50% (v/v) glycerol.

5X Transcription Buffer

200 mM Tris-HCI (pH 7.9 at 25 °C), 30 mM MgCl2, 50 mM DTT, 50 mM NaCl and 10 mM spermidine.

Inhibition and Inactivation

- Inhibitors: metal chelators, enzyme activity is reduced by 50% at NaCl or KCl concentration above 150 mM.
- Inactivated by heating at 70 °C for 10 min or by addition of EDTA.



For Research Use Only. Not for use in diagnostic procedures.

Protocol for in vitro transcription

 Linearize template DNA with a restriction enzyme. Extract DNA with phenol/chloroform, then with chloroform/isoamyl alcohol, and precipitate with ethanol. Dissolve DNA in DEPC-treated Water (#R0601).
Prenare the following reaction mixture:

5X Transcription buffer	10 μL			
ATP/GTP/CTP/UTP Mix, 10 mM each	10 μL (2 mM final concentration)			
Linear template DNA	1 µg			
Thermo Scientific RiboLock RNase Inhibitor (#EO0381)	1.25 μL (50 U)			
T7 RNA Polymerase	30 U			
DEPC-treated water (#R0601)	to 50 μL			

3. Incubate at 37 °C for 2 hours.

- 4. Optional: To remove template DNA add 2 µL (2 U) of DNase I, RNase-free (#EN0521), mix and incubate at 37 °C for 15 min.
- 5. Inactivate DNase I by phenol/chloroform extraction.

Note

- The transcription reaction should be performed under conditions that exclude contamination with RNases. The tips, tubes and water should be nuclease free. All the solutions should be made up in nuclease free water. Wearing gloves is advisable.
- The reaction mixture should be prepared at room temperature, since DNA may precipitate in the presence of spermidine at 4°C.
- Under the conditions described above, more than 10 µg RNA per 1 µg template DNA is obtained.
- The yield of proper length transcripts decreases if the template DNA is incompletely linearized due to a read-through reaction and accumulation of longer transcripts of a variable length.
- The reaction mixture can be scaled up or down.

Protocol for synthesis of radiolabeled RNA probes of high specific activity

1. Linearize template DNA with a restriction enzyme. Extract DNA with phenol/chloroform, then with chloroform/isoamyl alcohol, and precipitate with ethanol. Dissolve DNA in DEPC-treated Water (#R0601).

2. Prepare the following reaction mixture:

5X Transcription buffer	4 µL
3 NTP mix, 10 mM each	1 µL (0.5 mM final concentration)
100 μM CTP (#R0451)	2.4 µL (12 µM final concentration)
[α- ³² P]-CTP, ~30 TBq/mmol (800 Ci/mmol)	1.85 MBq (50 μCi)
Linear template DNA	0.2-1.0 µg
RiboLock™ RNase Inhibitor (#EO0381)	0.4 µL (20 U)
T7 RNA Polymerase	20 U
DEPC-treated water (#R0601)	to 20 µL

3. Incubate at 37 °C for 2 hours.

4. Stop the reaction by cooling at -20 °C.

5. Determine the percentage of label incorporated into RNA.

Note

• RNA synthesized under the conditions described above usually has a specific activity of 3-5 x10⁸dpm/µg.

• RNA can be radiolabeled with [³²P], [³⁵S] or [³H]-ribonucleotides. The use of 1.85 MBq (50 μCi) of 5'-[α-³²P]-CTP,

~30 TBq/mmol (800 Ci/mmol), 11.1 MBq (300μ Ci) of 5'-[α -³⁵S]-UTP, >37 TBq/mmol (>1000 Ci/mmol), 0.925 MBq (25 μ Ci) of 5,6-[³H]-UTP, 1.1-2.2 TBq/mmol (30-60 Ci/mmol) for 20 μ L reaction mixture is recommended.

• The yield of full-length transcripts is reduced when the final concentration of labeled NTP is below 12 µM.

References

1. Melton, D.A., et al., Efficient *in vitro* synthesis of biologically active RNA and RNA hybridization probes from plasmids containing a bacteriophage SP6 promoter, Nucleic Acids Res., 12, 7035-7056, 1984.

2. Krieg, P.A., Melton, D.A., Functional messenger RNAs are produced by SP6 *in vitro* transcription of cloned cDNAs, Nucleic Acids Res., 12, 7057-7070, 1984.

3. Melton, D.A., Injected antisense RNAs specifically block messenger RNA translation *in vivo*, Proc. Natl. Acad. Sci. USA, 82, 144-148, 1985.

4. Bernstein, E., et al., Role for bidentate ribonuclease in the initiation step of RNA interference, Nature, 409, 363-366, 2001.

5. Peebles, C.L., et al., A self-splicing RNA excises an intron lariat, Cell, 44, 213-223, 1986.

6. Church, G.M., Gilbert, W., Genomic sequencing, Proc. Natl. Acad. Sci. USA, 81, 1991-1995, 1984.

7. Witherell, G.W., et al., Cooperative binding of R17 coat protein to RNA, Biochemistry, 29, 11051-11057, 1990.

8. Krainer, A.R., et al., Normal and mutant human β-globin pre-mRNAs are faithfully and efficiently spliced *in vitro*, Cell, 36, 993-1005, 1984.

9. Jorgensen, E.D., et al., Specific contacts between the bacteriophage T3, T7, and SP6 RNA polymerases and their promoters, J. Biol. Chem., 266, 645-651, 1991.

Limited product warranty

Life Technologies Corporation and/or it affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale at www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have any questions, please contact Life Technologies at www.thermofisher.com/support.



Thermo Fisher Scientific Baltics UAB | V.A. Graiciuno 8, LT-02241 Vilnius, Lithuania

For descriptions of symbols on product labels or product documents, go to thermofisher.com/symbols-definition.

The information in this guide is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Important Licensing Information: These products may be covered by one or more Limited Use Label Licenses. By use of this product, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2019 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

