

## Certificate of Analysis

### T4 DNA Polymerase:

Part No.	Size (units)
M421A	100
M421F	500

**Enzyme Storage Buffer:** 200mM potassium phosphate (pH 6.5), 2mM DTT and 50% glycerol.

**T4 DNA Polymerase 10X Buffer:** The 10X Buffer supplied with this enzyme has a composition of 250mM Tris-Acetate (pH 7.7), 1M potassium acetate, 100mM magnesium acetate and 10mM DTT.

**Activity of T4 DNA Polymerase in Promega's Restriction Enzyme Buffers:** The activity of T4 DNA Polymerase in restriction enzyme buffers B, C, E and MULTI-CORE™ Buffer is at least 70% of the value obtained under the conditions of the unit activity assay. Activity in other restriction buffers may be 50% or less of the value obtained under the conditions of the unit activity assay.

**Heat-Inactivation:** T4 DNA Polymerase may be inactivated by incubation at 75°C for 10 minutes.

**Inhibitors:** T4 DNA Polymerase is reported to be inhibited by -SH blocking agents (1).

**Molecular Weight:** 112kDa.

**Source:** Purified from an *E. coli* strain expressing a recombinant clone.

**Storage Temperature:** Store at -20°C. Avoid multiple freeze-thaw cycles and exposure to frequent temperature changes. See the expiration date on the Product Information Label.

**Unit Definition:** One unit is defined as the amount of enzyme required to catalyze the incorporation of 55nmol of dTTP into acid-precipitable material in 30 minutes at 39°C using poly(dA):oligo(dT) as a substrate. See the unit concentration on the Product Information Label.

## Quality Control Assays

### Contaminant Activity

**Endonuclease Assay:** To test for endonuclease activity, 1µg of Type I supercoiled plasmid DNA is incubated with 40 units of T4 DNA Polymerase for one hour at 37°C in 1X restriction enzyme Buffer D. Following incubation, the supercoiled DNA is visualized on an ethidium bromide-stained agarose gel to verify the absence of visible nicking or cutting.

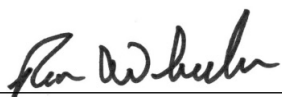
**Single-Strand Endonuclease Assay:** There must be no endonuclease activity detected when 1µg of M13mp18 DNA is incubated with 40 units of T4 DNA Polymerase for 1 hour at 37°C.

**Physical Purity:** The purity is >95% as judged by SDS-polyacrylamide gels with Coomassie® blue staining.

## Reference

1. Kornberg, A. (1992) *DNA Replication*, W.H. Freeman and Company, San Francisco, CA.

Signed by:



R. Wheeler, Quality Assurance

Part# 9PIM421

Revised 10/16



AF9PIM421 1016M421



**Promega**

### Promega Corporation

2800 Woods Hollow Road	
Madison, WI 53711-5399	USA
Telephone	608-274-4330
Toll Free	800-356-9526
Fax	608-277-2516
Internet	<a href="http://www.promega.com">www.promega.com</a>

### PRODUCT USE LIMITATIONS, WARRANTY, DISCLAIMER

Promega manufactures products for a number of intended uses. Please refer to the product label for the intended use statements for specific products. Promega products contain chemicals which may be harmful if misused. Due care should be exercised with all Promega products to prevent direct human contact.

Each Promega product is shipped with documentation stating specifications and other technical information. Promega products are warranted to meet or exceed the stated specifications. Promega's sole obligation and the customer's sole remedy is limited to replacement of products free of charge in the event products fail to perform as warranted. Promega makes no other warranty of any kind whatsoever, and SPECIFICALLY DISCLAIMS AND EXCLUDES ALL OTHER WARRANTIES OF ANY KIND OR NATURE WHATSOEVER, DIRECTLY OR INDIRECTLY, EXPRESS OR IMPLIED, INCLUDING, WITHOUT LIMITATION, AS TO THE SUITABILITY, PRODUCTIVITY, DURABILITY, FITNESS FOR A PARTICULAR PURPOSE OR USE, MERCHANTABILITY, CONDITION, OR ANY OTHER MATTER WITH RESPECT TO PROMEGA PRODUCTS. In no event shall Promega be liable for claims for any other damages, whether direct, incidental, foreseeable, consequential, or special (including but not limited to loss of use, revenue or profit), whether based upon warranty, contract, tort (including negligence) or strict liability arising in connection with the sale or the failure of Promega products to perform in accordance with the stated specifications.

© 1997-2013, 2016 Promega Corporation. All Rights Reserved.

Altered Sites and RiboClone are trademarks of Promega Corporation and are registered with the U.S. Patent and Trademark Office. MULTI-CORE is a trademark of Promega Corporation.

Coomassie is a registered trademark of Imperial Chemical Industries.

All specifications are subject to change without prior notice.

Product claims are subject to change. Please contact Promega Technical Services or access the Promega online catalog for the most up-to-date information on Promega products.

Part# 9PIM421

Printed in USA. Revised 10/16.

## I. Description

T4 DNA Polymerase can be used to fill 5' protruding ends with labeled or unlabeled dNTPs (1) or for the generation of blunt ends from DNA molecules with 3' overhangs (2).

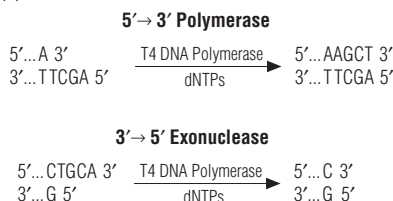


Figure 1. 5'→3' polymerase, and 3'→5' exonuclease activities of T4 DNA Polymerase.

## II. Standard Applications

### A. Filling 5' Overhangs With Unlabeled dNTPs

1. Digest 0.5–2.0µg DNA (in a volume of 50µl) with a restriction enzyme that leaves a 5' overhang.
2. Add 5 units of T4 DNA Polymerase/µg of DNA and 100µM of each dNTP. The recommended reaction buffer for T4 DNA Polymerase is 1X T4 DNA Polymerase Buffer. **Note:** T4 DNA Polymerase also functions well in many restriction enzyme reaction buffers.
3. Incubate at 37°C for 5 minutes. Stop the reaction by adding 2µl of 0.5M EDTA or by heating at 75°C for 10 minutes (3).

### B. Filling 5' Overhangs With Labeled dNTPs

1. After restriction enzyme digestion as described above, resuspend the DNA in restriction enzyme buffer **or** 1X T4 DNA Polymerase Buffer. Add 50µM of each unlabeled dNTP, 2µCi of the labeled [ $\alpha$ -<sup>32</sup>P]dNTP and 2.5 units of T4 DNA Polymerase in a total volume of 20µl.
2. Incubate for 5 minutes at 37°C. Stop the reaction by adding 1µl of 0.5M EDTA or by heating at 75°C for 10 minutes.

### C. Converting a 3' Overhang to a Blunt End

1. Digest 0.5–2.0µg of DNA (in a volume of 50µl) with a restriction enzyme that leaves a 3' overhang. Resuspend the DNA in restriction enzyme buffer **or** 1X T4 DNA Polymerase Buffer. Add 5 units of T4 DNA Polymerase/µg of DNA and 100µM of each dNTP. Incubate at 37°C for 5 minutes.
2. Stop the reaction by adding 2µl of 0.5M EDTA or by heating at 75°C for 10 minutes. With high concentrations of dNTPs (100µM), degradation of the DNA will stop at duplex DNA; however, if the dNTP supply is exhausted, the very active exonuclease activity (200 times more active than that of DNA polymerase I) will degrade the double-stranded DNA (4).

## D. Probe Synthesis

1. In a 20µl reaction volume, resuspend 0.5–2.0µg of linearized DNA in 1X T4 DNA Polymerase Buffer **or** restriction enzyme buffer. Add 1 unit of T4 DNA Polymerase for every microgram of DNA. The rate of exonuclease activity depends on the ratio of enzyme to DNA. After exonuclease digestion, add 1µl of a 2mM solution of each of the 3 unlabeled dNTPs, and add an amount of the [ $\alpha$ -<sup>32</sup>P]dNTP (400Ci/mmol) equivalent to the number of moles excised by the exonuclease.
2. Incubate the reaction at 37°C for 15 minutes. Stop the reaction by adding 1µl of 0.5M EDTA (5).

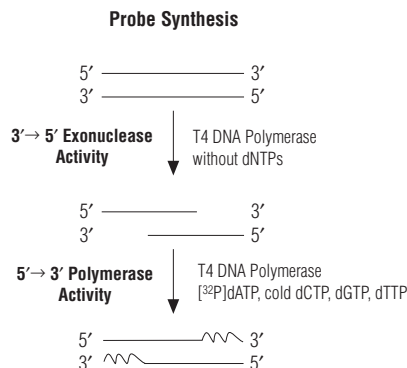


Figure 2. Probe synthesis with T4 DNA Polymerase.

## III. References

1. Challberg, M.D. and Englund, P.T. (1980) Specific labeling of 3' termini with T4 DNA polymerase. *Meth. Enzymol.* **65**, 39–43.
2. Burd, J.F. and Wells, R.D. (1974) Synthesis and characterization of the duplex block polymer d(C15A15)-d(T15G15). *J. Biol. Chem.* **249**, 7094–801.
3. Ausubel, F.M. *et al.* (1993) *Current Protocols in Molecular Biology*, Vol. 2, Greene Publishing Associates, Inc., and John Wiley and Sons, N.Y.
4. Sambrook, J. Fritsch, E.F. and Maniatis, T. (1989) *Molecular Cloning: A Laboratory Manual*, Cold Spring Harbor Laboratory, Cold Spring Harbor, N.Y.
5. O'Farrell, P.H., Kutter, E. and Nakanishi, M. (1980) A restriction map of the bacteriophage T4 genome. *Mol. Gen. Genet.* **179**, 421–35.