

Certificate of Analysis

DNA Polymerase I:

Part No.	Size (units)
M205A	500
M205B	2,500

DNA Polymerase I 10X Reaction Buffer (M195A): The DNA Polymerase 10X Reaction Buffer supplied with this enzyme has a composition of 500mM Tris-HCl (pH 7.2 at 25°C), 100mM MgSO₄ and 1mM DTT.

Enzyme Storage Buffer: DNA Polymerase I is supplied in 50mM Tris-HCl (pH 7.5 at 25°C), 1mM DTT, 0.1mM EDTA and 50% (v/v) glycerol.

Source: Purified from an *E. coli* strain expressing a recombinant clone (1).

Storage Temperature: See the Product Information Label for storage recommendations. 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 10nmol of deoxyribonucleotides into trichloroacetic acid insoluble form in 30 minutes at 37°C in 67mM potassium phosphate (pH 7.4), 6.7mM MgCl₂, 1mM DTT, 50µg/ml activated calf thymus DNA and 33µM each dNTP. See the unit concentration on the Product Information Label.

Quality Control Assays

Activity Assay

Unit Activity Assay: See unit definition above.

Contaminant Activity

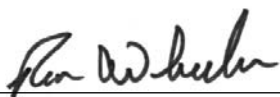
Endonuclease Assay: To test for endonuclease activity, 1µg of Type I supercoiled plasmid DNA is incubated in 25 units of DNA Polymerase I for 5 hours at 37°C. Following incubation, the supercoiled DNA is visualized on an ethidium bromide-stained agarose gel to verify the absence of visible nicking or cutting.

Physical Purity: The purity is ≥90% as judged by SDS-polyacrylamide gels with Coomassie® blue staining.

Reference

1. Kelley, W.S., Chalmers, K. and Murray, N.E. (1977) Isolation and characterization of a lambdapolA transducing phage. *Proc. Natl. Acad. Sci. USA* **74**, 5632-6.

Signed by:



R. Wheeler, Quality Assurance

Part# 9PIM205

Revised 4/18



AF9PIM205 0418M205



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	www.promega.com

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.

© 1997-2018 Promega Corporation. All Rights Reserved.

Coomassie is a registered trademark of Imperial Chemical Industries, Ltd.

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# 9PIM205
Printed in USA. Revised 4/18.

