

Certificate of Analysis

DNA-Dependent Protein Kinase:

Part No.	Size
V581A	2,500 units

Description: DNA-Dependent Protein Kinase (DNA-PK) consists of an approximate 460kDa catalytic subunit and a heterodimeric DNA-binding subunit (Ku) containing a 85kDa and a 70kDa peptide (1). It is purified from HeLa cells.

Storage Buffer: 25mM HEPES (pH 7.5), 50mM KCl, 0.2mM EDTA, 10mM MgCl₂, 1mM DTT, 10% glycerol.

Note: The storage buffer formulation has been changed to remove the IGEPAL CA-630 detergent.

Storage Conditions: See the storage recommendations on the Product Information Label. DNA-PK is stable at 4°C for 1 hour. Avoid multiple freeze-thaw cycles or exposure to frequent temperature changes. These fluctuations can greatly alter product stability. See the expiration date on the Product Information Label.

Unit Definition: One unit is the amount of enzyme required to incorporate 1pmol of phosphate into DNA-PK Peptide Substrate (Cat.# V5671) in one minute at 30°C.

Part# 9PIV581

Revised 8/16



AF9PIV581 0816V581

Quality Control Assays

Activation: The enzyme activity is increased by at least tenfold in the presence of 10µg/ml of linear, double-stranded DNA.

Activity Assay Conditions: 50mM HEPES (pH 7.5), 1mM DTT, 0.1mM EDTA, 0.2mM EGTA, 10mM MgCl₂, 0.1M KCl, 1.14mM DNA-PK Peptide Substrate (Cat.# V5671), 80µg/ml BSA, 0.2mM ATP and 10µg/ml linear double-stranded DNA.

Concentration: See the product information label for batch-specific information.

References

1. Gottlieb T.M. and Jackson S.P. (1993) The DNA-dependent protein kinase: Requirement for DNA ends and association with Ku antigen. *Cell* **72**, 131–42.
2. Carter, T. *et al.* (1990) A DNA-activated protein kinase from HeLa cell nuclei. *Mol. Cell. Biol.* **10**, 6460–71.
3. Smith, G.C.M. and Jackson, S.P. (1999) The DNA-dependent protein kinase. *Genes Dev.* **13**, 916–34.



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.

© 2001, 2005, 2006, 2010, 2011, 2015, 2016 Promega Corporation. All Rights Reserved.

Whatman is a registered trademark of Whatman Paper Company, 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.

Signed by:

R. Wheeler, Quality Assurance

Part# 9PIV581
Printed in USA. Revised 8/16.

Protocol for Use of DNA-Dependent Protein Kinase:

The following assay protocol may be used to verify the activity of purified DNA-PK. It also may be used as a basis for developing an assay for DNA-PK phosphorylation of protein substrates or for DNA-PK activity in cellular extracts. Dilute or dialyze DNA-PK samples to be assayed into DNA-PK dilution buffer.

Materials to Be Supplied by the User

(Solution compositions are provided below.)

- DNA-PK Peptide Substrate (Cat.# V5671)
- ATP, 10mM (Cat.# P1132)
- [γ - 32 P]ATP, 3,000Ci/mmol, 10 μ Ci/ μ l
- acetic acid, 15% and 30%
- Whatman® P-81 phosphocellulose paper
- DNA-PK activation buffer
- 5X DNA-PK reaction buffer
- 10mg/ml BSA

1. Prepare the following reaction as a positive control using a minimum of 10 units of DNA-PK. As additional controls, prepare two reactions lacking either the peptide substrate or the calf thymus DNA.

Component	Volume
5X DNA-PK reaction buffer	10 μ l
DNA-PK activation buffer	5 μ l
ATP, 10mM	1 μ l
DNA-PK Peptide Substrate, 10mg/ml	10 μ l
[γ - 32 P]ATP, 3,000Ci/mmol	0.2 μ l
10mg/ml BSA	0.4 μ l
DNA-PK (added last; see Note)	10–20u
water to final volume of	50 μ l

Before adding DNA-PK, pre-incubate the reaction tubes at 30°C for 3 minutes.

Note: In the presence of reaction buffer, DNA-PK can autophosphorylate and deactivate itself. Therefore, add the DNA-PK sample to the reaction last (2,3).

2. Incubate for 10 minutes at 30°C; then stop the reaction by adding 20 μ l of 30% acetic acid.
3. Spot 35 μ l of the reaction products onto a 2 x 2cm piece of Whatman® P-81 phosphocellulose paper. Allow the reaction products to soak into the paper (approximately 5 seconds).
4. Before the filters dry, wash the filters 5 times for 3–5 minutes each, with swirling, in 15% acetic acid; use 15ml per filter.
5. Using forceps, place the filters on a clean piece of filter paper and allow them to dry completely. Count the samples in a scintillation counter. Reactions using purified DNA-PK should exhibit >tenfold stimulation of 32 P incorporation when double-stranded DNA is added compared to control samples with no activation buffer.

Composition of Buffers and Solutions

5X DNA-PK reaction buffer

250mM	HEPES (pH 7.5)
500mM	KCl
50mM	MgCl ₂
1mM	EGTA
0.5mM	EDTA
5mM	DTT

DNA-PK dilution buffer (1ml)

990 μ l	1X DNA-PK reaction buffer
10 μ l	10mg/ml BSA

DNA-PK activation buffer

100 μ g/ml	calf thymus DNA in 1X TE
----------------	--------------------------