

Certificate of Analysis

PKG Kinase Enzyme System:

Cat.# V4248

Includes:

Part No.	Name	Size
V517A	cGMP-Dependent Protein Kinase (α isozyme)	6,000 units
V745A	cGMP-Dependent Protein Kinase Peptide Substrate (10mg/ml)	1mg
V307C	5X Reaction Buffer A	1.5ml
V641A	cGMP, 1mM	500 μ l
P117A	DTT, 100mM	100 μ l

Description: cGMP-Dependent Protein Kinase (PKG) is purified from bovine lung according to the method of Corbin and Døskeland (1). PKG is a serine/threonine protein kinase present in smooth muscle and a variety of other tissues, including lung, heart and Purkinje cells of the cerebellum (2,3). The gene sequence accession number is NP_776861.1.

Storage Buffer: 10mM potassium phosphate (pH 6.8), 1mM EDTA, 25mM β -mercaptoethanol, 150mM NaCl and 12% sucrose.

Storage Conditions: Avoid multiple freeze-thaw cycles or exposure to frequent temperature changes. These fluctuations can greatly alter product stability. For optimal storage, dispense PKG into smaller aliquots upon receipt and store at $< -65^{\circ}\text{C}$; store the rest of the components at -20°C . The enzyme is stable for 1 month at 4°C . See the expiration date on the Product Information Label.

Unit Definition: One unit is the amount of kinase required to incorporate 1pmol of phosphate into the test heptapeptide, RKRSRAE, per minute at 30°C . See Activity Assay in the Quality Control Assays section for buffer and conditions (4).

Part# 9PIV4248

Revised 4/18



AF9PIV42480418V4248



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

Quality Control Assays

Activity Assay: cGMP-Dependent Protein Kinase activity is determined in an assay buffer containing 97.2 μ g/ml heptapeptide substrate, 40mM Tris-HCl (pH 7.5), 20mM magnesium acetate, 0.2mM ATP, 2.86 μ M cGMP and 30,000cpm/ μ l γ - ^{32}P ATP. The reaction is run for 10 minutes at 30°C and terminated by spotting onto P81 filters.

Molecular Weight: cGMP-Dependent Protein Kinase demonstrates a 78kDa band by SDS-PAGE analysis.

Protein Concentration: Determined by Bradford Assay using BSA as a standard.

Purity: Greater than 90% as determined by SDS-PAGE analysis and Coomassie[®] blue staining.

Stimulation: Kinase activity is stimulated more than threefold by addition of 2.86 μ M cGMP.

References

1. Corbin, J.D. and Døskeland, S.O. (1983) Studies of two different intrachain cGMP-binding sites of cGMP-dependent protein kinase. *J. Biol. Chem.* **258**, 11391-7.
2. Edelman, A.M., Blumenthal, D.K. and Krebs, E.G. (1987) Protein serine/threonine kinases. *Ann. Rev. Biochem.* **56**, 567-613.
3. Beebe, S.J. and Corbin, J.D. (1986) In: *The Enzymes*, Vol. 17, 3rd., Boyer, P.D. and Krebs, E.D., eds.
4. Roskoski, R. (1983) Assays of protein kinase. *Methods Enzymol.* **99**, 3-6.

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.

ADP-Glo is a trademark of Promega Corporation.

© 2012-2018 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.

Signed by:

R. Wheeler, Quality Assurance

Part# 9PIV4248
Printed in USA. Revised 4/18.

Protocol

The ADP-Glo™ Kinase Assay can be used to measure the activity of PKG by quantifying the amount of ADP produced during a kinase reaction. The following is only a brief protocol. For more detailed and complete protocols, see the *ADP-Glo™ Kinase Assay Technical Manual* #TM313 found at: www.promega.com/protocols, and Kinase Enzyme Systems Protocol at: www.promega.com/KESProtocol

1. **Kinase Reaction:** In a 384-well low-volume plate (5µl total reaction volume) or 96-well plate (25µl total reaction volume), perform PKG kinase reaction using 1X Reaction buffer A supplemented with 50µM DTT and containing the following:
 - Kinase (desired amount)
 - 1µg of cGMP-Dependent Protein Kinase Peptide Substrate
 - 2.86µM cGMP
 - ATP (see concentration in Application Note available on the protocol page for Cat.# V4248*)
 - Inhibitor (or DMSO for control)

Incubate at room temperature for 30–60 minutes or any desired time.

2. **ADP-Glo™ Kinase Assay:**
 - Add 5µl of ADP-Glo™ Reagent. (Add 25µl if using a 96-well plate.)
 - Incubate at room temperature for 40 minutes.
 - Add 10µl of Kinase Detection Reagent. (Add 50µl if using a 96-well plate.)
 - Incubate at room temperature for 30–60 minutes.
 - Record luminescence (Integration time: 0.5–1 second).