

Product Contents

Anti-Luciferase pAb:

Part No.	Size
G745A	200µg

Description: Anti-Luciferase pAb is a purified polyclonal antibody raised in goats against recombinant firefly luciferase (*Photinus pyralis*).

Formulation: Anti-Luciferase pAb is supplied in PBS containing 50µg/ml gentamicin.

Protein Concentration: 1.0mg/ml.

Storage Conditions: Store at 4°C for daily/weekly use, or aliquot and store at -20°C for long-term storage. Avoid multiple freeze-thaw cycles and exposure to frequent temperature changes.

Usage Note: Concentration gradients may form in frozen products and should be dispersed upon thawing. Mix well prior to use.

The performance of this product is guaranteed for six months from the date of purchase if stored and handled properly.

Part# 9PIG745

Revised 1/05



Quality Control Assays

Purity: This antibody is greater than 90% pure IgG as determined by SDS-PAGE followed by Coomassie® blue staining and visual inspection.

Immunocytochemistry: The antibody labels cultured HeLa cells that have been transfected with recombinant firefly luciferase (*Photinus pyralis*) when used at a concentration of 20µg/ml in conjunction with an anti-goat FITC conjugate as a detection system. Nontransfected cells show less labeling.

Western Blot: Western blot analysis is performed using the Anti-Luciferase pAb at a 1:1,000 dilution along with a donkey anti-goat IgG AP conjugate at a 1:5,000 dilution and Western Blue® Stabilized Substrate for Alkaline Phosphatase (Cat.# S3841). Western blot analysis shows that the antibody detects 2ng of recombinant firefly luciferase (*Photinus pyralis*).



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.

© 1999, 2005 Promega Corporation. All Rights Reserved.

Western Blue is a trademark of Promega Corporation and is registered with the U.S. Patent and Trademark Office.

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# 9PIG745
Printed in USA. Revised 1/05