

## Product Contents

### Anti-GFAP pAb:

Part No.	Size
G560A	100µg

**Description:** Anti-GFAP (Glial Fibrillary Acidic Protein) pAb is raised in rabbits against bovine GFAP and purified using Protein A.

**Formulation:** Anti-GFAP 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# 9PIG560

Revised 6/05



AF9PI G560 0605 G560

## Quality Control Assays

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

**Immunohistochemistry:** The antibody, when used at a 1:1,000 dilution and detected with a donkey anti-rabbit Cy<sup>3</sup> antibody conjugate, labels astrocytes in rat spinal cord sections. Preincubation of the antibody with a ten-fold excess of purified GFAP abolishes astrocyte labeling.

**Western Blot:** Western blot analysis, performed using the Anti-GFAP pAb at a 1:1,000 dilution along with Anti-Rabbit IgG (Fc), AP Conjugate (Cat.# S3731) at a 1:2,000 dilution and Western Blue® Stabilized Substrate for Alkaline Phosphatase (Cat.# S3841), shows that the antibody detects 10ng of purified GFAP.



**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.

© 1999, 2005 Promega Corporation. All Rights Reserved.

Western Blue is a registered trademark of Promega Corporation.

Coomassie is a registered trademark of Imperial Chemical Industries, Ltd. Cy is a registered trademark of Amersham Biosciences, 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# 9PIG560  
Printed in USA. Revised 6/05