

Product Contents

Anti-Human p75 pAb:

Part No.	Size
G323A	200µg

Description: Anti-Human p75 pAb is generated in rabbits (1–3).

Applications: For immunocytochemistry and immunohistochemistry, use antibody at 1–10µg/ml; for immunoprecipitation, use antibody at 1µg/ml.

Formulation: Provided at 1mg/ml as frozen liquid in PBS containing 50µg/ml gentamycin.

Immunogen: Recombinant Human p75 cytoplasmic domain.

Storage Conditions: Store long-term (infrequent use or 1–2 times per month) at –20°C. For daily/weekly use, store at 4°C. Avoid multiple freeze-thaw cycles and exposure to frequent temperature changes. The performance of this product is guaranteed for 6 months from date of purchase if stored and handled properly.

Usage Notes: Anti-Human p75 pAb is raised against a GST fusion protein produced in *E. coli*. This antibody is not suitable for use in Western blot analysis because of the potential for cross-reactivity with the fusion protein.

Quality Control Assays

Activity Assays

Cross-Reactivity: Anti-Human p75 pAb cross-reacts with human, mouse, rat and chicken p75 due to high conservation of the cytoplasmic receptor domain.

Immunocytochemistry: Anti-Human p75 pAb, at a concentration of 2µg/ml (1:500), labels PC12 cells. Pretreatment of the antibody with recombinant rat p75-GST fusion protein at 40µg/ml inhibits this labeling.

Purity: Anti-Human p75 pAb is >90% pure as determined by SDS-PAGE.

References

1. Dobrowsky, R.T. *et al.* (1994) Activation of the sphingomyelin cycle through the low-affinity neurotrophin receptor. *Science* **265**, 1596–9.
2. Chao, M.V. (1994) The p75 neurotrophin receptor. *J. Neurobiol.* **25**, 1373–85.
3. Chao, M.V. and Hempstead, B.L. (1995) p75 and Trk: A two-receptor system. *Trends Neurosci.* **18**, 321–6.

Part# 9PIG323

Revised 8/04



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.

© 1998–2004 Promega Corporation. All Rights Reserved.

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# 9PIG323
Printed in USA. Revised 8/04