

Product Contents

Anti-TrkB In pAb:

Part No.	Size
G156A	100µg

Description: Anti-TrkB In pAb is a polyclonal antibody raised in chickens against a fusion protein containing amino acids 482–501 (HISNGSNTSPSSSEGGPDAVI) of the carboxy terminus of the mouse trkB intracellular domain. The immunogen is injected into chickens, and IgY is isolated from egg yolk using the EGGstract® IgY Purification System (Cat.# G2610).

Formulation: Anti-TrkB In pAb is supplied in 10mM potassium phosphate (pH 7.4) containing 100mM sodium chloride. There are no preservatives added.

Immunogen: The immunogen is a fusion protein generated in *E. coli* using a proprietary technology containing the HISNGSNTSPSSSEGGPDAVI sequence expressed with a novel synthetic carrier protein.

Protein Concentration: 0.5mg/ml.

Storage Conditions: Store in aliquots at –20°C. For daily/weekly use store at 4°C. Avoid multiple freeze-thaw cycles.

Part# 9PIG156

Revised 5/05



Quality Control Assays

Immunohistochemistry: The antibody labels the cell bodies of Purkinje cells in fixed/frozen rat cerebellum sections when used at 5µg/ml in conjunction with a peroxidase/diaminobenzidine detection system.

Sensitivity: A 96-well plate is coated with immunogen at a concentration of 1µg/ml. Anti-TrkB In pAb is added to the first row of wells at 10µg/ml and is then two-fold serially diluted to 0.01µg/ml and detected using Anti-Chicken IgY HRP Conjugate (Cat.# G1351) and TMB substrate. The antibody must demonstrate an OD of twice background at <0.4µg/ml.



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.

© 2000, 2005 Promega Corporation. All Rights Reserved.

Products may be covered by pending or issued patents or may have certain limitations. Please visit our Web site for more information.

EGGstract is a registered trademark of Promega Corporation.

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# 9PIG156
Printed in USA. Revised 5/05