

## Product Contents

### Anti-ACTIVE® MAPK pAb, Rabbit (pTEpY):

Part No.	Size
V803A	40µl

**Description:** Anti-ACTIVE® MAPK pAb is a polyclonal rabbit antibody. The antibody is affinity purified using a dually phosphorylated peptide that corresponds to the active form of the mitogen-activated protein (MAP) kinase enzymes.

**Immunogen:** Dually phosphorylated Thr/Glu region (pTEpY) derived from the active form of MAP kinase enzymes.

**Storage Buffer:** Anti-ACTIVE® MAPK pAb is supplied in PBS.

**Storage Conditions:** Store at -20°C. Avoid multiple freeze-thaw cycles and exposure to frequent temperature changes. If product is to be used in the short term, store at 4°C after adding sodium azide or thimerosal as a preservative. The performance of this product is guaranteed for at least six months from the date of purchase if stored and handled properly.

**Usage Note:** For Western blotting, the suggested working dilution of Anti-ACTIVE® MAPK pAb is 1:5,000.

Part# 9PIV803

Revised 3/05



## Quality Control Assays

**Western Blotting:** In all cases, Western blot analysis is performed using Anti-ACTIVE® MAPK pAb at a 1:5,000 dilution, along with Donkey Anti-Rabbit IgG (H+L), AP, Anti-ACTIVE® Qualified (Cat.# V7971) at a 1:5,000 dilution and Western Blue® Stabilized Substrate for Alkaline Phosphatase (Cat.# S3841).

**Sensitivity:** Using Western blot analysis, the antibody detects MAPK enzyme bands at 42/44kDa in 1µg of NGF-treated PC12 cell extract.

**Specificity:** Using Western blot analysis, the intensity of the MAPK enzyme bands at 42/44kDa in 10µg of quiescent (-NGF) PC12 cell extract is less than their intensity in 1µg of NGF-treated PC12 cell extract.

**Immunocytochemistry:** When used at 1:500 dilution along with a donkey anti-rabbit Cy™3 detection system, the Anti-ACTIVE® MAPK pAb strongly labels NGF-treated PC12 cells, while only background levels of staining are seen with untreated PC12 cells. The protocol used can be found in Promega Technical Bulletin #TB262.



**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.

© 1998-2005 Promega Corporation. All Rights Reserved.

Anti-ACTIVE and Western Blue are trademarks of Promega Corporation and are registered with the U.S. Patent and Trademark Office.

Cy is a trademark of Amersham Pharmacia Biotech 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# 9PIV803  
Printed in USA. Revised 3/05