

Product Contents

Anti-ACTIVE® JNK pAb, Rabbit, (pTPpY):

Part No.	Size
V793A	40µl
V793B	120µl

Description: Anti-ACTIVE® JNK pAb is a polyclonal antibody from rabbit serum. The antibody is affinity purified using a dually phosphorylated peptide that corresponds to the active form of the JNK enzymes.

Immunogen: Dually phosphorylated JNK-derived peptide.

Storage Buffer: 10mM sodium phosphate (pH 7.4) containing 20mM sodium chloride and no preservative.

Storage Conditions: Store at -20°C for long-term storage. For daily or weekly use store at 4°C after adding preservative. Avoid multiple freeze-thaw cycles. The performance of this product is guaranteed for six months from the date of purchase if stored and handled properly.

Part# 9PIV793

Revised 6/05



Quality Control Assays

Western Blotting: Western blot analysis is performed using the Anti-ACTIVE® JNK pAb, Rabbit, at a 1:5,000 dilution, followed by detection with a 1:10,000 dilution of Donkey Anti-Rabbit IgG AP Conjugate (Cat.# V7971) and a colorimetric detection system. See *Anti-ACTIVE® MAPK, p38 and JNK Polyclonal Antibodies and Anti-ACTIVE® Qualified Secondary Antibody Conjugates Technical Bulletin #TB262* for more details.

Sensitivity: Using Western blot analysis, the antibody can detect 2ng of purified recombinant active JNK enzyme.

Specificity: The antibody can detect activated JNK in 20µg of sorbitol-treated PC12 cell extracts. The signal from sorbitol-treated PC12 cells is greater than with 20µg of cell extract from quiescent (untreated) PC12 cells.

Immunocytochemistry: When used at a 1:1,000 dilution along with an anti-rabbit Cy®3 conjugate detection system, the Anti-ACTIVE® JNK pAb, Rabbit, strongly labels sorbitol-treated PC12 cells and weakly labels untreated cells except for an occasional positive cell in the untreated cell population. See *Anti-ACTIVE® MAPK, p38 and JNK Polyclonal Antibodies and Anti-ACTIVE® Qualified Secondary Antibody Conjugates Technical Bulletin #TB262* for more details.



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.

© 2001, 2005 Promega Corporation. All Rights Reserved.

Anti-ACTIVE is a registered trademark of Promega Corporation.

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# 9PIV793
Printed in USA. Revised 6/05.