

Product Contents

Anti-pS⁴⁷³ Akt pAb

Part No.	Size
G744A	40µl

Description: Anti-pS⁴⁷³ Akt pAb is an affinity-purified polyclonal rabbit antibody. The antibody is purified using a phosphorylated peptide that corresponds to the phospho-S⁴⁷³ form of Akt-1.

Formulation: Anti-pS⁴⁷³ Akt pAb is supplied in PBS containing 50µg/ml gentamicin.

Immunogen: Phosphorylated Akt-1-derived peptide.

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 Notes:

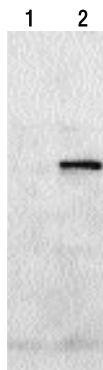
1. Concentration gradients may form in frozen products and should be dispersed upon thawing. Mix well prior to use.
2. The suggested working dilution for Western blotting is a 1:2,500 dilution.

The performance of this product is guaranteed for six months from the date of purchase if stored and handled properly.

Quality Control Assays

Western Blot: Western blot analysis is performed using rabbit Anti-pS⁴⁷³ Akt pAb at a 1:2,500 dilution, along with Donkey Anti-Rabbit, IgG, HRP (H+L) Conjugate (Cat.# V7951) at a 1:10,000 dilution and detected with a chemiluminescence-based detection method (see the protocol on the reverse).

Sensitivity: Western blot analysis shows that the Anti-pS⁴⁷³ Akt antibody detects a phospho-Akt enzyme band at approximately 60kDa in 10µg of activated (PDGF-treated) NIH/3T3 cell extract. No band is detected in 10µg of untreated cell extract.



Detection of phosphorylated Akt by Western blot analysis with Anti-pS⁴⁷³ Akt pAb. NIH/3T3 total cell extracts (10µg per lane) were resolved by polyacrylamide gel electrophoresis and blotted to nitrocellulose. Lane 1: Untreated cells. Lane 2: Cells pretreated with PDGF at 50ng/ml for 20 minutes. Anti-pS⁴⁷³ Akt pAb was used at a 1:2,500 dilution. Detection was performed using chemiluminescence.

Part# 9PIG744

Revised 6/05



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.

© 1999, 2005 Promega Corporation. All Rights Reserved.

Anti-ACTIVE is a registered trademark of Promega Corporation.

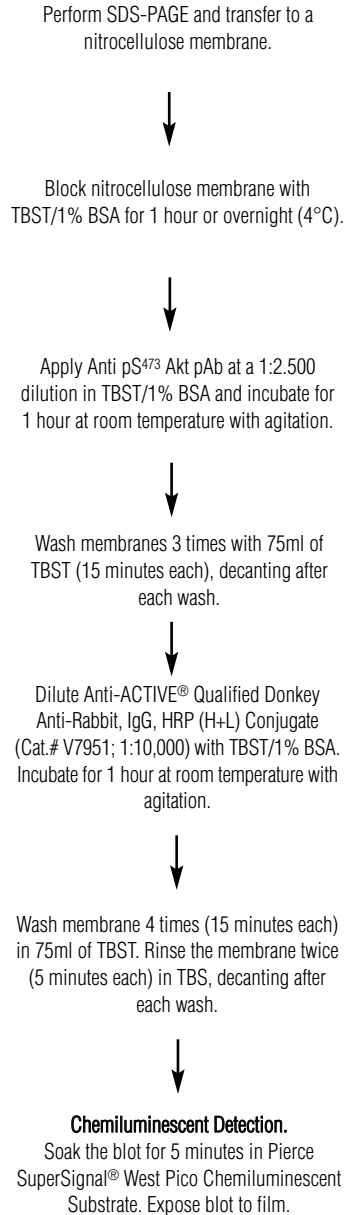
SuperSignal is a registered trademark of Pierce Chemical Company.

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# 9PIG744
Printed in USA. Revised 6/05

Usage Information



Schematic diagram illustrating the use of nitrocellulose membranes in Western blot analysis. Note: It may be necessary to empirically determine the optimal dilutions of primary antibodies for your system. Use of secondary antibodies other than those available from Promega may require additional optimization. TBST = Tris-buffered saline with 0.05% Tween® 20. BSA = Bovine Serum Albumin.