



EAGLE
BIOSCIENCES

**Polyclonal Protein Kinase
Consensus Sequence
Antibody**

Product Number: PSA00-A1000

Aliquot: 1.0 mL

Storage: -20°C

PRODUCT SPECIFICATION SHEET

**Immunoglobulin
type:**

Polyclonal rabbit IgG

Specificity:

This antibody was generated against the G-T-P-E-Y-L-A-P-E peptide and is excellent for Western blot detection of protein kinases containing this sequence.

Applications:

A dilution of 1: 50 is recommended for Western blotting using alkaline phosphatase. A further dilution may be possible with more sensitive detection systems.

Note:

Use of 1% normalized goat serum added to a usual blocking agent is recommended to reduce non-specific interactions.

Contents/vial:

1.0 ml of rabbit antiserum in PBS

Storage:

Stable for several weeks at 4°C, several months at -20°C, and indefinitely at -70°C. Do not refreeze this product.

Product Developed and Manufactured in the USA

EAGLE BIOSCIENCES, INC.
20A NW Blvd., Suite 112
Nashua, NH 03063

Phone: 617-419-2019
Fax: 617-419-1110
Email: info@EagleBio.com
www.EagleBio.com