

Anti-Phosphoserine (Anti-pS)

Rabbit Polyclonal Antibody

Catalog # S95-65R

Lot # E360-3

Cited Applications

For WB, ELISA, IP, IHC

Ideal working dilutions for each application should be empirically determined by the investigator.

Specificity

Recognizes peptides and proteins phosphorylated on serine residues.

Cross Reactivity

Pan-specific antibody. No reaction to phosphotyrosine peptides and proteins.

Host/Isotype/Clone#

Rabbit

Immunogen

Phosphoserine peptide-KLH conjugates

Formulation

PBS, 50% glycerol, pH7.

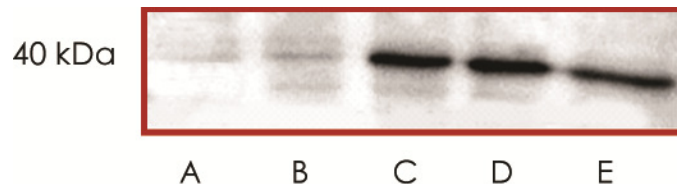
Stability

Store at 4°C (add 0.1% NaN₃) for several months, and at -20°C for longer periods. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles.

Scientific Background

Post-translational modification of proteins and peptides is a robust way to regulate function of existing proteins or peptides. Protein phosphorylation represents one of the most abundant and important post-translational modifications of proteins. Phosphorylation on serine residues in proteins is performed by a variety of serine specific protein kinases in the cell which are involved in a wide variety of signaling networks. The Phospho-Serine antibody detects phosphorylation on serine residues in proteins and peptides.

Sample Data



Representative western blot of the phosphoserine protein profile with Anti-Phosphoserine (1:250) using lysates of human melanoma cells (MMRU) (A), increased UV-treated MMRU (B, C, D) and UV-treated MMRU in phosphatase inhibitor condition (E).

Anti-Phosphoserine (Anti-pS)

Rabbit Polyclonal Antibody

Catalog Number

S95-65R

Specific Lot Number

E360-3

Purification

Affinity chromatography

Concentration

0.25 µg/µl

Stability

1 yr at -20°C from date of shipment

Storage & Shipping

Store product at -20°C. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on ice packs.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.