

RhoA Mutant (S188A) Protein

Human recombinant protein expressed in E.coli cells

Catalog # R44-32AH

Lot # H2788-1

Product Description

Recombinant human RhoA (S188A) (1-190) was expressed in E. coli cells using an N-terminal His tag. The gene accession number is [NM_001664](#).

Gene Aliases

ARH12; ARHA; RHO12; RHOH12

Formulation

Recombinant protein stored in 50mM sodium phosphate, pH 7.0, 300mM NaCl, 150mM imidazole, 0.1mM PMSF, 0.25mM DTT, and 25% glycerol.

Storage and Stability

Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

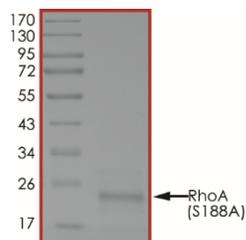
Scientific Background

RhoA is a member of the Rho gene family that interacts with protein kinases and serves as a target of activated GTPase. RhoA play an important role in the initial events of cell protrusion, whereas Rac1 and Cdc42 activate pathways implicated in reinforcement and stabilization of newly expanded protrusions. Rho mediates corneal epithelial migration in response to external stimuli by regulating the organization of the actin cytoskeleton (1). The inhibition of RhoA signaling is required for both vaccinia morphogenesis and virus-induced cell motility. Vaccinia inhibits RhoA signalling using the viral protein F11 and this can enhance the spread of infection (2). It has been reported that cGMP/cGMP-dependent protein kinase can phosphorylate RhoA at Ser188 *in vitro* (3).

References

1. Wu, K. Y.et.al: Local translation of RhoA regulates growth cone collapse. Nature 436: 1020-1024, 2005.
2. Valderrama, F.et.al: Vaccinia virus-induced cell motility requires F11L-mediated inhibition of RhoA signaling. Science 311: 377-381, 2006.
3. Sawada N. et.al: cGMP-dependent protein kinase phosphorylates and inactivates RhoA. Biochem. Biophys. Res. Commun. 280:798-805(2001).

Purity



The purity of RhoA (S188A) was determined to be **>80%** by densitometry. Approx. MW **23kDa**.

RhoA Mutant (S188A) Protein

Human recombinant protein expressed in E. coli cells

Catalog #	R44-32AH
Lot #	H2788-1
Purity	>80%
Concentration	0.1 µg/µl
Stability	1yr at -70°C from date of shipment
Storage & Shipping	Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.

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.