

RGS1 Protein

Full-length recombinant protein expressed in E. coli cells

Catalog # R39-30H

Lot # M035-4

Product Description

Recombinant full-length human RGS1 Protein was expressed in E. coli cells using an N-terminal His tag. The gene accession number is [BC015510](#).

Gene Aliases

IR20, IER1, 1R20, BL34

Formulation

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

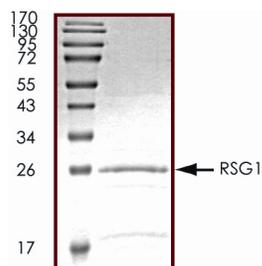
Scientific Background

RGS1 is a member of the regulator of G-protein signalling family and attenuates the signalling activity of G-proteins. RGS1 binds to activated GTP-bound G alpha subunit and acts as a GTPase activating protein (GAP) thereby increasing the rate of conversion of GTP to GDP and terminating the signal (1). RGS1 is extensively up-regulated in renal cell carcinoma (RCC) tissues and melanoma. In melanoma, RGS1 expression is significantly correlated with increased tumor thickness, mitotic rate and presence of vascular involvement. Furthermore, there is significant association between increasing RGS1 expression and reduced relapse-free survival as well as disease-specific survival (DSS) survival (2).

References

1. Sierra D A, et al: Evolution of the regulators of G-protein signaling multigene family in mouse and human. *Genomics* 79: 177-185, 2002.
2. Rangel J, et al: Novel role for RGS1 in melanoma progression. *Am J Surg Pathol.* 2008 Aug;32(8):1207-12

Purity



The purity of RGS1 was determined to be **>90%** by densitometry. Approx. MW **26kDa**.

RGS1 Protein

Full-length recombinant protein expressed in E. coli cells

Catalog Number	R39-30H
Specific Lot Number	M035-4
Purity	>90%
Concentration	0.2 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.