

GNA13 Protein

Full-length recombinant human protein expressed in E. coli cells

Catalog # G118-30G

Lot # L1885-8

Product Description

Full-length recombinant human GNA13 protein was expressed in E. coli cells using an N-terminal GST tag. The GNA13 gene accession number is [NM_006572](#).

Gene Aliases

G13; G alpha-14

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 150mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 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.

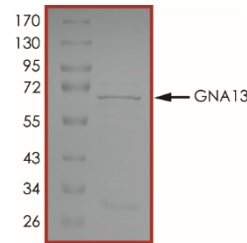
Scientific Background

G-alpha 13 (GNA13) is a guanine nucleotide-binding protein of the G12 class of G-alpha proteins. It activates Rho and regulates cytoplasmic as well as nuclear signaling events such as activation of the Jun N-terminal kinase signaling module, Na^+/H^+ exchangers, focal adhesion assemblies, and transcriptional activation of specific primary response genes. Like GNA12, GNA13 proteins are found up-regulated in several human cancers, and the GNA12/GNA13 signaling plays an important role in cancer cell invasion and metastasis.

References

- Cheong, S. C. et al. Gene expression in human oral squamous cell carcinoma is influenced by risk factor exposure. *Oral Oncol* 45, 712 (2009).
- Kelly, P. A Role for the G12 Family of Heterotrimeric G Proteins in Prostate Cancer Invasion. *J Biol Chem* 281, 26483 (2006).
- <http://www.phosphosite.org/proteinAction?id=2145&showAllSites=true>

Purity



The purity of GNA13 was determined to be **>70%** by densitometry, approx. MW **67 kDa**.

GNA13 Protein

Full-length recombinant human protein expressed in E. coli cells

Catalog #	G118-30G
Lot #	L1885-8
Purity	>70%
Concentration	0.05 µg/µl
Stability	1 yr 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.