

GNAZ Protein

Recombinant human protein expressed in E. coli cells

Catalog # G131-30G

Lot # Q2270-6

Product Description

Recombinant human GNAZ (2-end) protein was expressed in E. coli cells using an N-terminal GST tag. The GNAZ gene accession number is [NM_002073](#).

Gene Aliases

GZ-alpha

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.

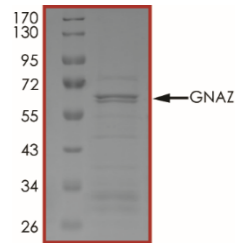
Scientific Background

G-alpha (z) (GNAZ) is a guanine nucleotide-binding protein that belongs to the G-alpha family, G(i/o/t/z) subfamily.

References

- <http://www.phosphosite.org/proteinAction.action?id=4926&showAllSites=true>

Purity



The purity of GNAZ was determined to be **>70%** by densitometry, approx. MW **65 kDa**.

GNAZ Protein

Recombinant human protein expressed in E. coli cells

Catalog #	G131-30G
Lot #	Q2270-6
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.