

## GNA11 Protein

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

**Catalog # G116-30G**

Lot # L2232-5

### Product Description

Full-length recombinant human GNA11 protein was expressed in E. coli cells using an N-terminal GST tag. The GNA11 gene accession number is [NM\\_002067](#).

### Gene Aliases

FBH; FBH2; FHH2; GNA-11; HHC2; HYPOC2

### 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  $-70^{\circ}\text{C}$ . 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.

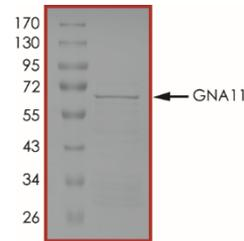
### Scientific Background

Guanine nucleotide-binding proteins (G proteins) are heterotrimeric signal-transducing molecules consisting of alpha, beta, and gamma subunits, and function as transducers downstream of G protein-coupled receptors (GPCRs) in numerous signaling cascades. The alpha subunit binds guanine nucleotide, can hydrolyze GTP, and can interact with other proteins. Guanine nucleotide-binding protein subunit alpha-11 (GNA11) acts as an activator of phospholipase C. Involved in signaling of gonadotropin-releasing hormone receptor, negatively regulating cell growth. Down-regulation may be involved in human breast cancers.

### References

- <http://www.phosphosite.org/proteinAction?id=1101&showAllSites=true>
- <http://www.uniprot.org/uniprot/P29992>

### Purity



The purity of GNA11 was determined to be **>70%** by densitometry, approx. MW **66 kDa**.

## GNA11 Protein

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

Catalog #	G116-30G
Lot #	L2232-5
Purity	>70%
Concentration	0.05 µg/µl
Stability	1yr at $-70^{\circ}\text{C}$ from date of shipment
Storage & Shipping	Store product at $-70^{\circ}\text{C}$ . 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](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

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