

## FRS3 Protein

Full-length recombinant protein expressed in Sf9 cells

### Catalog # F80-30G

Lot # I313-2

### Product Description

Recombinant full-length human FRS3 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM\\_006653](#).

### Gene Aliases

SNT2, FRS2B, FRS2beta, MGC17167

### 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.

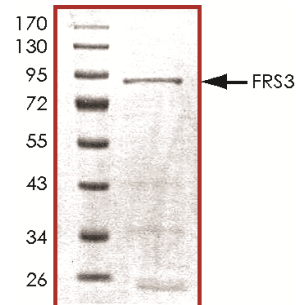
### Scientific Background

FRS3 or fibroblast growth factor receptor substrate 3 is a peripheral plasma membrane protein that is a substrate for the fibroblast growth factor receptor. During FGF or NGF stimulation, FRS3 becomes tyrosine phosphorylated and then serves as a platform for the recruitment of multiple signaling proteins for activation of the Ras-MAP kinase signaling cascade (1). FRS3 is mainly expressed in fibroblast and myoblast cell lines and undergo robust tyrosine phosphorylation in response to several mitogenic ligands. Through interaction with Rho family of small GTPases, FRS3 has been implicated in the reorganization of the actin cytoskeleton and subsequent morphological changes in various cells (2).

### References

- Huang L, et al: SNT-2 interacts with ERK2 and negatively regulates ERK2 signaling in response to EGF stimulation. *Biochem Biophys Res Commun*. 2004 Nov 19;324(3):1011-7.
- Xu H, et al: Novel recognition motif on fibroblast growth factor receptor mediates direct association and activation of SNT adapter proteins. *J. Biol. Chem.* 273: 17987-17990, 1998.

### Purity



The purity of FRS3 was determined to be **>80%** by densitometry. Approx. MW **87kDa**.

## FRS3 Protein

Full-length recombinant protein expressed in Sf9 cells

Catalog Number	F80-30G
Specific Lot Number	I313-2
Purity	>80%
Concentration	0.2 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.**