

GRB7 Protein

Full length human recombinant protein expressed in E.coli

Catalog # G45-30GG

Lot # 1052-3

Product Description

Recombinant full-length human GRB7 was expressed by E.coli using an N-terminal GST tag. The gene accession number is [NM_001030002](#).

Gene Aliases

B47

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.

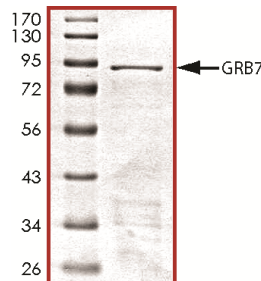
Scientific Background

GRB7 or growth factor receptor-bound protein 7 belongs to a small family of adapter proteins which interact with a number of receptor tyrosine kinases and signaling molecules. GRB7 encodes a growth factor receptor-binding protein that interacts with epidermal growth factor receptor (EGFR) and ephrin receptors. GRB7 is highly expressed in liver and kidney (1). GRB7 isoforms are involved in cell invasion and metastatic progression of human esophageal carcinomas (2). GRB7 plays an important role in the integrin signaling pathway and cell migration by binding with focal adhesion kinase (FAK).

References

- Skolnik, E. Y. et.al: Cloning of PI3 kinase- associated p85 utilizing a novel method for expression/cloning of target proteins for receptor tyrosine kinases. Cell 65: 83-90, 1991.
- Tanaka, S. et.al: A novel variant of human Grb7 is associated with invasive esophageal carcinoma. J. Clin. Invest. 102: 821-827, 1998.

Purity



The purity of GRB7 was determined to be **>90%** by densitometry. Approx. MW **91 kDa**.

GRB7 Protein

Full length human recombinant protein expressed in E.coli

| | |
|---------------------|---|
| Catalog Number | G45-30GG |
| Specific Lot Number | 1052-3 |
| 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.