

## Recombinant PIK3R1 protein

**Catalog No:** 81097, 81897

**Lot No:** 32317001

**Expressed In:** Baculovirus

**Quantity:** 20, 1000 µg

**Concentration:** 0.4 µg/µl

**Source:** Human

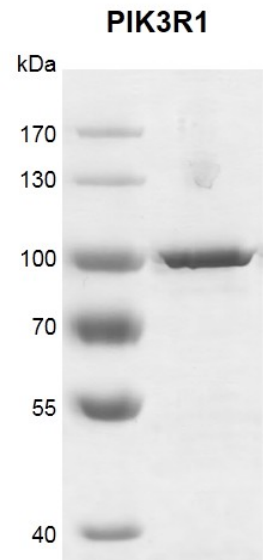
**Buffer Contents:** Recombinant PIK3R1 protein is supplied in 25 mM HEPES-NaOH pH 7.5, 300 mM NaCl, 10% glycerol, 0.04% Triton X-100 and 0.5 mM TCEP.

**Background:** PIK3R1 (Phosphoinositide-3-Kinase Regulatory Subunit 1), also known as GRB1, p85-ALPHA or p85, is a regulatory subunit of phosphatidylinositol 3-kinase. It can bind to activated (phosphorylated) protein-Tyr kinases, through its SH2 domain, and acts as an adapter, mediating the association of the p110 catalytic unit to the plasma membrane. It is necessary for the insulin-stimulated increase in glucose uptake and glycogen synthesis in insulin-sensitive tissues. And it plays an important role in signaling in response to FGFR1, FGFR2, FGFR3, FGFR4, KITLG/SCF, KIT, PDGFRA and PDGFRB, likewise, plays a role in ITGB2 signaling. PIK3R1 also modulates the cellular response to ER stress by promoting nuclear translocation of XBP1 isoform 2 in a ER stress-and/or insulin-dependent manner during metabolic overloading in the liver and hence plays a role in glucose tolerance improvement.

**Protein Details:** Recombinant human PIK3R1 protein was expressed in a baculovirus expression system as the full length protein (accession number NP\_852664.1) with an N-terminal FLAG tag. The molecular weight of the protein is 84.9 kDa.

**Application Notes:** This protein is suitable for use in binding assays, inhibitor screening, and selectivity profiling.

**Storage and Guarantee:** Recombinant proteins in solution are temperature sensitive and must be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is for research use only and is not for use in diagnostic procedures. This product is guaranteed for 6 months from date of arrival.



**Recombinant PIK3R1 protein gel**  
7.5% SDS-PAGE Coomassie staining  
MW: 84.9 kDa  
Purity: >95%