

PFKFB2 Protein

Full length recombinant protein expressed in Sf9 cells

Catalog # P322-30G

Lot # Z1335-4

Product Description

Recombinant full-length human PFKFB2 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM_006212](#).

Gene Aliases

PFK-2/FBPase-2

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 50mM 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.

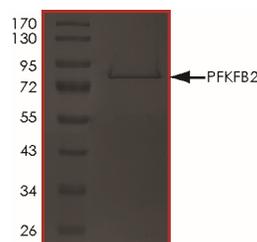
Scientific Background

PFKFB2 is involved in both the synthesis and degradation of fructose-2,6-bisphosphate, a regulatory molecule that controls glycolysis in eukaryotes. PFKFB2 has a 6-phosphofructo-2-kinase activity that catalyzes the synthesis of fructose-2,6-bisphosphate, and a fructose-2,6-bisphosphatase activity that catalyzes the degradation of fructose-2,6-bisphosphate. PFKFB2 regulates fructose-2,6-bisphosphate levels in the heart, while a related enzyme encoded by a different gene regulates fructose-2,6-bisphosphate levels in the liver and muscle (1). PFKFB2 also functions as a homodimer. PFKFB2 play an important role in the regulation of muscle glycolysis during the first moments of exercise (2).

References

- Hilliker, C. E. et.al: Human and rat chromosomal localization of two genes for 6-phosphofructo-2-kinase/fructose-2,6-bisphosphatase by analysis of somatic cell hybrids and in situ hybridization. *Genomics* 10: 867-873, 1991.
- Rovira J. et.al: Upregulation of heart PFK-2/FBPase-2 isozyme in skeletal muscle after persistent contraction. *Pflugers Arch.* 2012 Apr;463(4):603-13. doi: 10.1007/s00424-011-1068-5.

Purity



The purity of PFKFB2 was determined to be **>90%** by densitometry. Approx. MW **84 kDa**.

PFKFB2 Protein

Full length recombinant protein expressed in Sf9 cells

Catalog # P322-30G

Lot # Z1335-4

Purity >90%

Concentration 0.1 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.