

## PFKFB1 Protein

Full length recombinant protein expressed in Sf9 cells

**Catalog # P321-30G**

Lot # J617-2

### Product Description

Recombinant full-length human PFKFB1 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM\\_002625](#).

### Gene Aliases

F6PK; HL2K; PFRX

### 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.

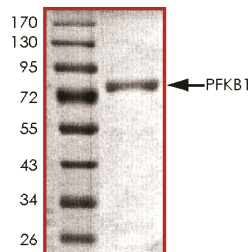
### Scientific Background

PFKFB1 is a member of the family of bifunctional 6-phosphofructo-2-kinase: fructose-2, 6-biphosphatase enzymes which forms a homodimer that catalyzes both the synthesis and degradation of fructose-2,6-biphosphate using independent catalytic domains (1). Fructose-2,6-biphosphate is an activator of the glycolysis pathway and an inhibitor of the gluconeogenesis pathway. Regulation of fructose-2,6-biphosphate levels through the activity of this enzyme is thought to regulate glucose homeostasis (2).

### References

- Algaier, J.et.al: Molecular cloning, sequence analysis, and expression of a human liver cDNA coding for fructose-6,2-kinase:fructose-2,6-bisphosphatase. *Biochem. Biophys. Res. Commun.* 153: 328-333, 1988.
- Darville, M. I. et.al: Complete nucleotide sequence coding for rat liver 6-phosphofructo-2-kinase/fructose-2,6-bisphosphatase derived from a cDNA clone. *FEBS Lett.* 224: 317-321, 1987.

### Purity



The purity of PFKFB1 was determined to be **>95%** by densitometry. Approx. MW **80 kDa**.

## PFKFB1 Protein

Full length recombinant protein expressed in Sf9 cells

Catalog Number	P321-30G
Specific Lot Number	J617-2
Purity	>95%
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](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.**