

Data Sheet

PFKFB4

Human, recombinant, N-terminal His-tag

Catalog #: 71199

Lot #: 141209

Conc.: 1.77 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 200 mM imidazole and 20% glycerol.

Stability: At least 6 months at -80°C . Avoid freeze/thaw cycles. Storing diluted enzyme is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).

References:

1. Minchenko, O.H., *et al.*, *Acta Biochimica Polonica*. 2005; **52(4)**: 881-888.
2. Goidts, V., *et al.*, *Oncogene*. 2012; **31**: 3235-3243.

Description: Human recombinant 6-phosphofructo-2-kinase/fructose-2,6-bisphosphatase 4 (PFKFB4), also known as 6PF-2-K/Fru-2,6-P2ase 4, and PFK/FBpase 4, a.a. 2-469 with an N-terminal His-tag, GenBank Accession No. NM_004567/NM_004558, MW = 55 kDa, expressed in an *E. coli* cell expression system.

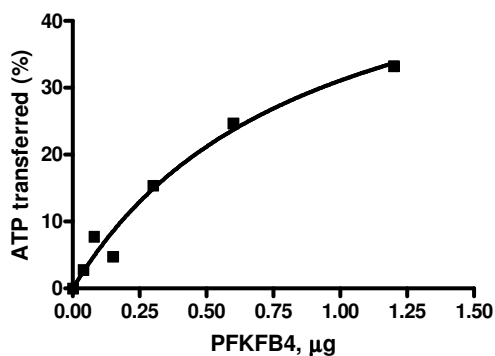
Specific Activity: 16 pmol/min/ μg

Assay Conditions: The reaction mixture contained 50 mM Tris-HCl, pH 7.5, 12.5 mM MgCl_2 , 1 mM β -D-fructose-6-phosphate, 20 μM ATP, 2 mM DTT, 5 mM KH_2PO_4 , and various amounts of PFKFB4 in a final volume of 50 μL . The reaction was incubated at 30°C for 30 min, the amount of ATP consumed was calculated using kinase-glo plus reagent (Promega).

Application: Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance

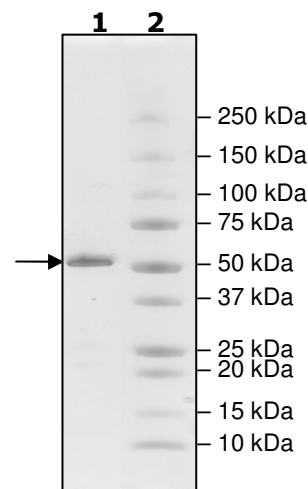
Specific Activity



**4-20% SDS-PAGE
Coomassie staining**

Lane 1:
2 μg PFKFB4
Lane 2:
Protein Marker

MW: 55 kDa
Purity: $\geq 90\%$



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com
 Please visit our website at: www.bpsbioscience.com