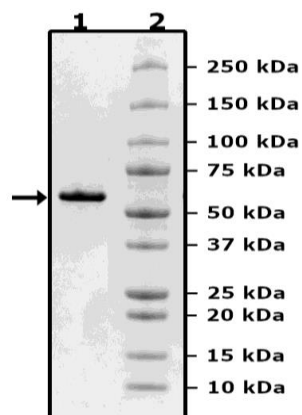


Product Information

Construct:	PKM2 (PKM2 Variant 1) (His-2-531(end))
Concentration:	0.73 mg/ml
Species:	Human
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 87 mM Imidazole, 20% glycerol, and 3 mM DTT
Expression System:	<i>E. coli</i>
Format:	Aqueous buffer solution.
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	NM_002654
MW:	59 kDa
Purity:	≥90%
Specific Activity:	≥60 pmol/min/μg
Assay Conditions:	Assay was done in buffer containing 50 mM Tris (pH 7.4), 10 mM MgCl ₂ , 50 mM KCl, 2 mM ADP, 10 mM phosphoenolpyruvate (PEP) and PKM2 (BPS Bioscience #50295) titrated at various concentrations. Reactions were done in 50 μl duplicates and incubated at room temperature for 15 minutes. ATP production was detected using Kinase-Glo [®] Max (Promega).
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



PKM2 Activity

