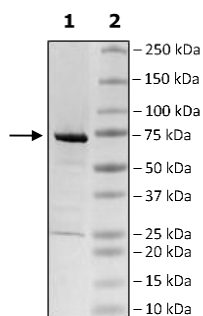


Product Information

Construct:	GSK3 β (GST-2-433(end)-His)
Concentration:	0.32 mg/ml
Species:	Human
Formulated In:	40 mM Tris-HCl pH 8.0, 110 mM NaCl, 2.2 mM KCl, 3 mM DTT, 250 mM Imidazole and 20% glycerol.
Expression System:	Sf9
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	NM_002093
MW:	75 kDa
Purity:	$\geq 90\%$
Specific Activity:	≥ 300 pmol/min/ μ g
Assay Conditions:	Assay was done according to the GSK3 β Assay Kit (BPS Bioscience #79700) with GSK3 β titrated from 48 ng/rxn – 0.75 ng/rxn.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



Specific Activity

