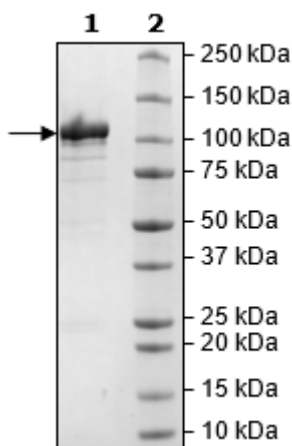


Product Information

Description:	Recombinant human IKK β (inhibitor of nuclear factor kappa B kinase subunit beta), full length encompassing amino acids 2-756 (end). This construct contains an N-terminal GST-tag. A short peptide linker placed between GST and the protein includes a PreScission Protease Cleavage Site to facilitate removal of the GST portion. The recombinant protein was affinity purified.
Species:	Human
Construct:	IKK β (GST-2-756(end))
Concentration:	0.17 mg/ml
Expression System:	Sf9
Purity:	$\geq 90\%$
Format:	Aqueous buffer solution.
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 20% glycerol, and 3 mM DTT
MW:	113 kDa
Genbank Accession:	NM_001556
Stability:	At least 6 months at -80°C .
Storage:	-80°C
Instructions for Use:	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.
Specific Activity:	≥ 40 pmol/min/ μg
Assay Conditions:	Assay was done using the IKK β Kinase Assay Kit (BPS Bioscience #78507) with various amounts of IKK β .
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



IKK β Activity

