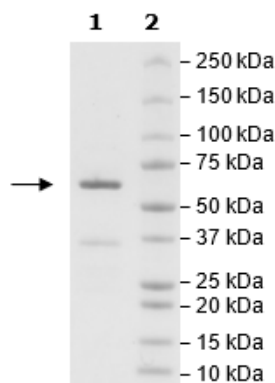


Product Information

Construct:	Aurora Kinase B (GST-2-344(end)-His)
Concentration:	0.98 mg/ml
Species:	Human
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 20% glycerol, and 3 mM DTT
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_004217
MW:	66 kDa
Purity:	80%
Specific Activity:	≥2.3 pmol/min/μg
Assay Conditions:	Assay was done in Kinase Buffer 1 (BPS Bioscience #79334) using 50 μM DTT, 0.1 mg/ml MBP as a substrate with 20 μM ATP incubated at Room Temperature for 60 min. The amount of ATP transferred was calculated using ADP-Glo reagent (Promega).
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



Aurora B Protein Kinase Activity

