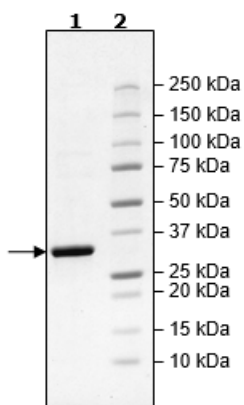


Product Information

Construct:	Abl Kinase (His-248-531)
Concentration:	1.00 mg/ml
Species:	Human
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 3 mM DTT, 130 mM Imidazole, 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_007313
MW:	35 kDa
Purity:	≥90%
Specific Activity:	2,700 pmol/min/μg
Assay Conditions:	Assay was done in a kinase buffer containing 0.2 mM DTT using Abltide substrate (0.2 mg/ml) and 20 μM ATP. Reaction was done at 30°C for 40 min. Amount of ATP transferred was calculated using Kinase-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



Specific Activity

