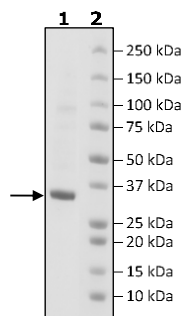


Product Information

Construct:	ABL (T315I) (His-229-512)
Concentration:	2.91 mg/ml
Species:	Human
Formulated In:	25 mM Tris-HCl, pH 8.0, 100 mM NaCl, 0.05% Tween-20, 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_005157
MW:	35 kDa
Purity:	86%
Specific Activity:	2500 pmol/min/μg
Assay Conditions:	ABL (T315I) activity was confirmed using the ABL (T315I) Assay Kit (BPS Bioscience, #79576) and following the protocol provided with the kit.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



Specific Activity

