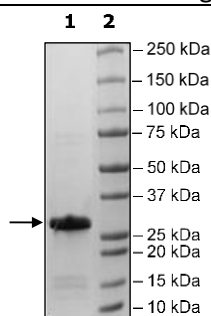


Product Information

Construct:	BCL-XL (1-212-His)
Concentration:	0.97 mg/ml
Species:	Human
Formulated In:	45 mM Tris-HCl, pH 8.0, 124 mM NaCl, 2.4 mM KCl, 10% glycerol, 225 mM imidazole and 3 mM DTT
Expression System:	<i>E. coli</i>
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	Z23115
MW:	25 kDa
Purity:	≥90%
Assay Conditions:	Assay was performed according to the BCL-XL TR-FRET Assay Kit (BPS Bioscience, #50223) with BCL-XL titrated from 42 ng/rxn – 0 ng/rxn.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Control Data

4-20% SDS-Page Coomassie Staining



BCL-XL Activity

