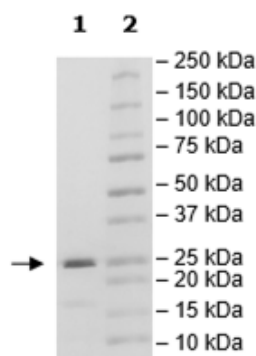


Product Information

Construct:	BCL-2 (1-211-His)
Concentration:	0.78 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:	<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:	NM_000633
MW:	25 kDa
Purity:	≥90%
Assay Conditions:	Assay was done according to BCL-2 TR-FRET Assay Kit (BPS Bioscience #50222) with BCL2 titrated from 12 ng/rxn-0 ng/rxn.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



BCL-2 Activity

