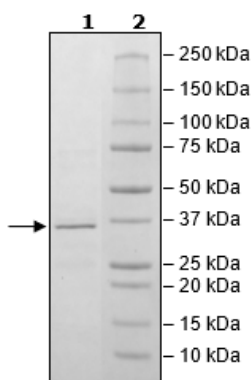


Product Information

Construct:	MCL1 (His-2-329) (Rabbit)
Concentration:	0.40 mg/ml
Species:	Rabbit
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 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:	XP_002715631.2
MW:	36 kDa
Purity:	84%
Assay Conditions:	Assay run according to MCL-1 TR-FRET Assay Kit (BPS Bioscience Kit #79506). Enzyme mixture incubated for 1 hour at room temperature and TR-FRET measured at ex 340 nm / em 620 nm, 665 nm.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



Rabbit MCL-1:MCL-1 Ligand [Biotinylated] Binding Activity

