

## Product Information

<b>Construct:</b>	MCL1 (His-2-327)
<b>Concentration:</b>	0.81 mg/ml
<b>Species:</b>	Human
<b>Formulated In:</b>	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 3 mM DTT, 20% glycerol, and 200 mM Imidazole
<b>Expression System:</b>	<i>E. coli</i>
<b>Format:</b>	Aqueous buffer solution
<b>Stability:</b>	At least 6 months at -80°C. Avoid freeze/thaw cycles.
<b>Storage:</b>	-80°C
<b>Genbank Accession:</b>	NM_021960
<b>MW:</b>	36 kDa
<b>Purity:</b>	≥90%
<b>Assay Conditions:</b>	Figure 1: Inhibition of MCL-1 by A-1210477, was measured using the MCL-1 TR-FRET Assay Kit (BPS Bioscience #79506). Figure 2: Assay run according to MCL-1 TR-FRET Assay Kit (BPS Bioscience #79506).
<b>Applications:</b>	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

## Quality Control Data

## 4-20% SDS-PAGE Coomassie Staining

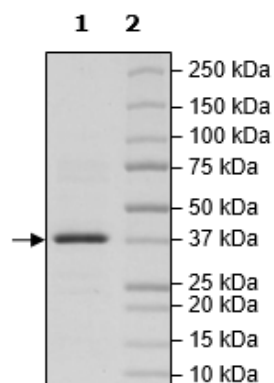


Figure 1: Inhibition of MCL1 Binding

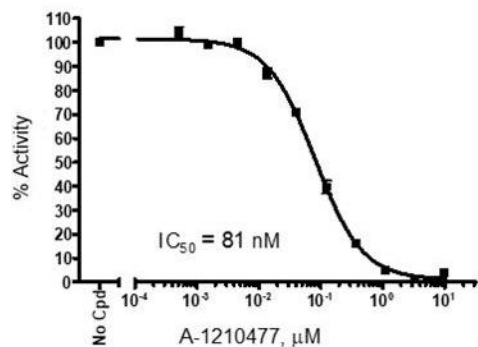


Figure 2: Human MCL-1:MCL-1 Ligand [Biotinylated] Binding

