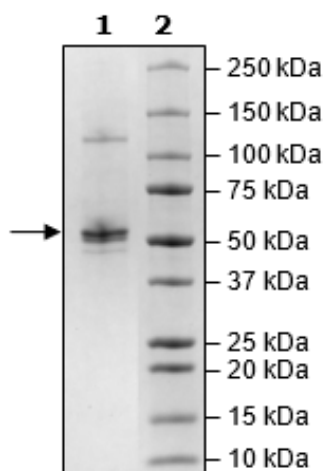


Product Information

Construct:	Cereblon (FLAG-2-442(end))
Concentration:	0.13 mg/ml
Species:	Human
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 20% glycerol, and 100 µg/ml FLAG peptide
Expression System:	HEK293
Format:	Aqueous buffer solution.
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	NM_016302
MW:	52 kDa + glycans
Glycosylation:	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
Purity:	84%
Assay Conditions:	Assay was performed according to Cereblon Binding Assay Kit (BPS Bioscience #79899) with CRBN titrated from 2400 ng/rxn – 0 ng/rxn incubated for 90 minutes at room temperature. Fluorescence polarization measured at ex635 nm, em682 nm.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



Cereblon Activity

