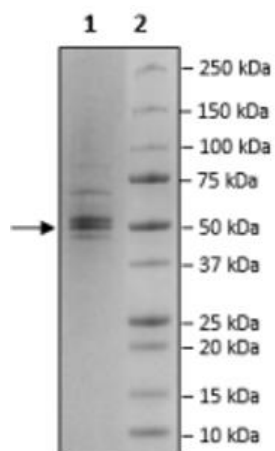


Product Information

Description:	Recombinant human cereblon (CRBN), full length, encompassing amino acids 2-422(end). This construct contains an N-terminal FLAG-tag. The recombinant protein was affinity purified.
Background:	The Cereblon (CRBN) protein via its interaction with the DNA damage-binding protein-1 (DDB1), Cullin 4 (Cul4A or Cul4B), and regulator of Cullins 1 (RoC1) forms a functional E3 Ub ligase complex. In this complex, Cereblon functions as a substrate receptor that mediates the ubiquitination and subsequent proteasomal degradation of target proteins. Like most E3 ligases, Cereblon complex ubiquitinates itself and its auto-ubiquitination promotes its ligase activity. Cereblon complex is involved in many biological processes including cell proliferation and apoptosis, and it has been extensively used in targeted protein degradation (PROTAC®), where CRBN binds to the protein of interest and leads to its degradation. This approach has shown promising results in the treatment of cancer.
Species:	Human
Construct:	Cereblon (FLAG-2-442(end))
Concentration:	0.31 mg/ml
Expression System:	HEK293
Purity:	76%
Format:	Aqueous buffer solution.
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
MW:	51 kDa
Genbank Accession:	NM_016302
Stability:	At least 6 months at -80°C.
Storage:	-80°C
Instructions for Use:	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.
Assay Conditions:	Assay was done according to Cereblon Binding Assay Kit (BPS Bioscience #79899) with various amounts of Cereblon.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



Cereblon Activity

