

## Background

Cereblon (CRBN) complex is an E3 Ligase that mediates ubiquitination and proteasomal degradation of target proteins. CRBN functions as a substrate adaptor, providing substrate specificity without possessing inherent enzymatic activity. It is linked to the scaffolding protein Cullin 4A (Cul4A) and its regulator, the RING-box protein (RBX1), via DNA damage-binding protein 1 (DDB1). The ligase activity of the complex is determined by the Cullin–RBX1 module, which catalyzes the transfer of ubiquitin from the RBX1-bound E2 enzyme to target substrates.

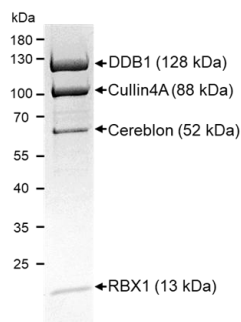
## Application(s)

- Protein degradation
- PROTAC and Molecular Glue discovery
- Selectivity Profiling

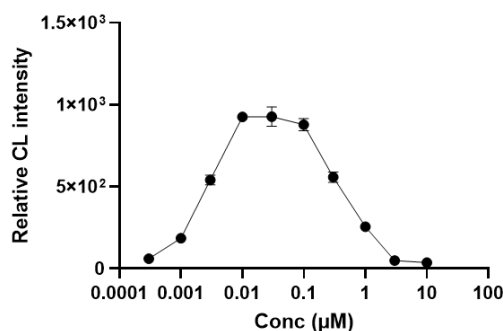
## Product Specifications

<b>Affinity Tag</b>	None
<b>Purity</b>	≥ 90% estimated by SDS-PAGE
<b>Molecular Weight</b>	CRBN: 51 kDa, DDB1: 128 kDa, Cul4A: 88 kDa, Rbx1: 13 kDa
<b>Quantity</b>	10 µg, 50 µg
<b>Species</b>	Human. Genbank Accession No.: CRBN, NM_016302; DDB1, NM_001923; Cul4A, NM_003589; Rbx1, NM_014248
<b>Expression System</b>	HEK293
<b>Physical State</b>	Liquid
<b>Buffer</b>	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 20% glycerol
<b>Stability &amp; Storage</b>	1 year at -80°C. Avoid repeated freeze/thaw cycles

## Product QC



**SDS-PAGE analysis of purified CRBN complex.** Twenty µg of the protein was loaded on a 4-20% SDS-PPAGE gel and stained with Coomassie brilliant blue



**In vitro ubiquitination assay to test the activity of the CRBN complex.** In vitro ubiquitination reaction was performed in the presence of various doses of LC2, a VHL degrader of KRAS G12C. Ubiquitinated KRAS G12C was captured on the microtiter plate coated with TUBEs and detected using anti-KRAS antibody. Chemiluminescence intensities were plotted against PROTAC doses to evaluate the extent of ubiquitination.

## References

1. Meyers M, et al., *ACS Chem Biol.* 2024;19(1):58-68.
2. Barankiewicz J, et al., *Cancers (Basel).* 2022;14(18):4492.
3. Gang, Lu., et al., *Science.* 2014; 343(6168): 305-309.
4. Zhu, Y.X., et al., *Blood.* 2011; 118: 4771-4779.