

Product Information

Description:	Complex of recombinant full length human Cereblon, amino acids 1-442(end), and full length DDB1 (damage specific DNA binding protein 1), amino acids 1-1140(end). The cereblon construct includes an N-terminal FLAG-tag, while the DDB1 construct contains a C-terminal Strep-tag. The proteins were co-expressed and purified as a complex.
Background:	Cereblon (CRBN) is the substrate-binding component of the E3 protein ligase complex DDB1-CUL4A-RBX1, which is involved in the ubiquitination and proteasomal degradation of target proteins. Binding of CRBN to a substrate engages its E3 ligase activity and results in the ubiquitination and ultimate degradation of the substrate. Many proteins are known targets of CRBN, including several transcription factors, growth factors and kinases. CRBN has become a target of choice for the development of many PROTACs of therapeutical interest.
Species:	Human
Construct:	Cereblon (FLAG-1-442(end)) / DDB1 (1-1140(end)-Strep)
Concentration:	0.13 mg/ml
Expression System:	HEK293
Purity:	77%
Format:	Aqueous buffer solution.
Formulated In:	45 mM Tris-HCl, pH 8.0, 124 mM NaCl, 2.4 mM KCl, and 10% glycerol
MW:	Cereblon: 53 kDa; DDB1: 131 kDa
Genbank Accession:	Cereblon: NM_016302; DDB1: NM_001923
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.
Applications:	Useful for SDS-PAGE.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining

