

PRODUCT INFORMATION



CRBN^{mid} (human, recombinant)

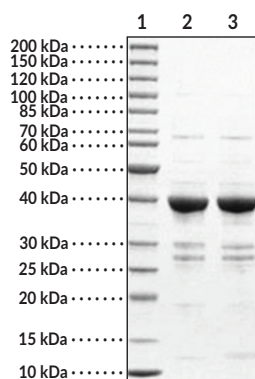
Item No. 45179

Overview and Properties

Synonym:	Cereblon ^{mid}
Source:	Recombinant human N-terminal His-tagged CRBN ^{mid} expressed in <i>E. coli</i>
Amino Acids:	41-187-GSG-249-426
Uniprot No.:	Q965W2
Molecular Weight:	39.8 kDa
Storage:	-80°C (as supplied)
Stability:	≥1 year
Purity:	≥90% estimated by SDS-PAGE
Supplied in:	50 mM HEPES, pH 7.5, with 1 mM DTT, 200 mM sodium chloride, and 20% glycerol
Endotoxin Testing:	<1.0 EU/μg, determined by the LAL endotoxin assay
Protein Concentration:	<i>batch specific</i> mg/ml

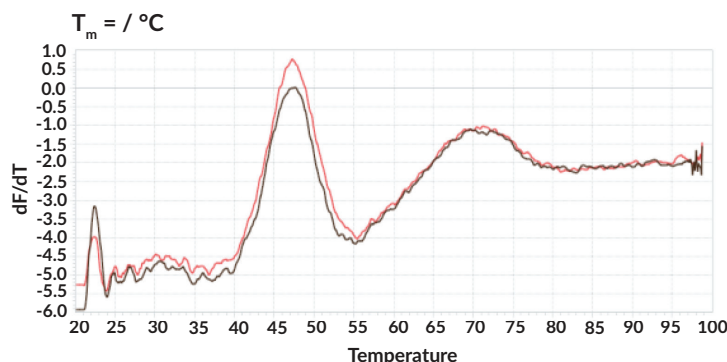
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Images



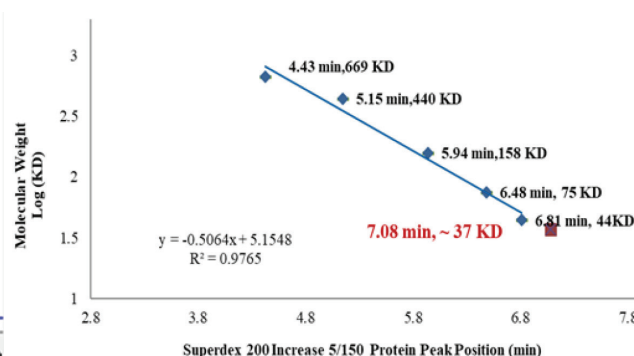
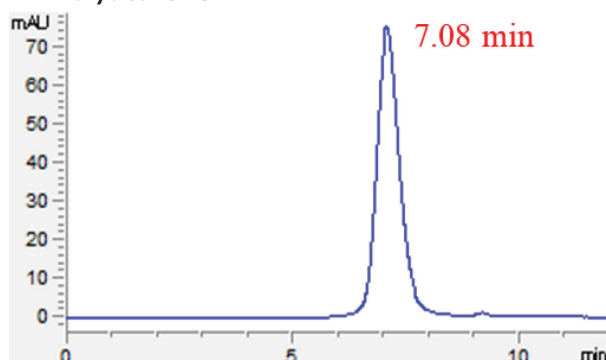
Lane 1: MW Markers
Lane 2: CRBN^{mid} (2 μg) (reduced)
Lane 3: CRBN^{mid} (2 μg) (non-reduced)

SDS-PAGE Analysis of CRBN^{mid}. This protein has a calculated molecular weight of 39.8 kDa.



Melting Profile of Purified CRBN^{mid}

Analytical SEC



WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 02/24/2026

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

Cereblon (CRBN) is the substrate receptor component of the cullin 4 RING ligase (CRL4) complex.^{1,2} It is composed of an N-terminal unstructured region, a Lon protease-like domain, a helical bundle, and a thalidomide-binding domain.^{1,3} CRBN is ubiquitously expressed and localizes to the nucleus, cytoplasm, and plasma membrane.¹ As part of the CRL4 complex with cullin 4, RING box protein 1 (RBX1), and DNA damage-binding protein 1 (DDB1), it facilitates the ubiquitination of substrates for proteasomal degradation. CRBN is involved in the regulation of ion channel expression, AMPK activation, energy metabolism, and endoplasmic reticulum (ER) stress.^{1,2,4} Mutations in CRBN are associated with autosomal recessive non-syndromic intellectual disability, and CRBN is also the target of thalidomide-induced teratogenicity.^{1,2,5} CRBN-targeting ligands, including thalidomide, pomalidomide, and lenalidomide, have been widely used to induce protein degradation using proteolysis-targeting chimera (PROTAC) technology.⁶ CRBN^{midi} is an engineered construct that contains the Lon protease-like domain, a Gly-Ser-Gly linker, part of the helical bundle, and the thalidomide-binding domain with C78I, I92V, K116N, Q134E, R283W, C287N, V293S, G302D, L342R, C343E, T359I, and L423I stabilizing mutations.³ These modifications enable the expression and purification of CRBN without co-expression of DDB1 in *E. coli* and the formation of binary and ternary complexes with degraders using crystallography. Cayman's CRBN^{midi} (human, recombinant) protein has a calculated molecular weight of 39.8 kDa.

References

1. Kim, H.K., Ko, T.H., Nyamaa, B., *et al.* Cereblon in health and disease. *Pflugers Arch* **468(8)**, 1299-1309 (2016).
2. Kim, H.K., Seol, J.E., Ahn, S.W., *et al.* Cereblon: Promise and challenges for combating human diseases. *Pflugers Arch* **473(11)**, 1695-1711 (2021).
3. Kroupova, A., Spiteri, V.A., Rutter, Z.J., *et al.* Design of a Cereblon construct for crystallographic and biophysical studies of protein degraders. *Nat. Commun.* **15(1)**, 8885 (2024).
4. Park, N., Marquez, J., Pham, T.K., *et al.* Cereblon contributes to cardiac dysfunction by degrading Ca_v1.2α. *Eur. Heart J.* **43(20)**, 1973-1989 (2022).
5. Ito, T., Ando, H., Suzuki, T., *et al.* Identification of a primary target of thalidomide teratogenicity. *Science* **327(5971)**, 1345-1350 (2010).
6. Lai, A.C. and Crews, C.M. Induced protein degradation: An emerging drug discovery paradigm. *Nat. Rev. Drug Discov.* **16**, 101-114 (2017).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM