

PRODUCT INFORMATION



Cullin 2/RBX1 Complex (human, recombinant)

Item No. 45183

Overview and Properties

Synonyms: Cullin 2: CUL2
RBX1: E3 Ubiquitin-protein Ligase RBX1, E3 Ubiquitin-protein Transferase RBX1, Protein ZYP, Regulator of Cullins-1, RING Box Protein 1, RING Finger Protein 75, RNF75, ROC1

Source: Active recombinant human N-terminal His-Strep II-tagged cullin 2/RBX1 expressed in insect cells

Amino Acids: 2-745 (cullin 2) (full length)

Uniprot Nos.: Q13617/P62877

Molecular Weight: 0.1 (cullin 2) and 12.3 (RBX1) kDa

Storage: -80°C (as supplied)

Stability: 1 year

Purity: ≥90% estimated by SDS-PAGE

Supplied in: 50 mM HEPES, pH 7.5, with 200 mM sodium chloride, 20% glycerol, and 1 mM DTT

Endotoxin Testing: <1.0 EU/μg, determined by the LAL endotoxin assay

Protein

Concentration: *batch specific* mg/ml

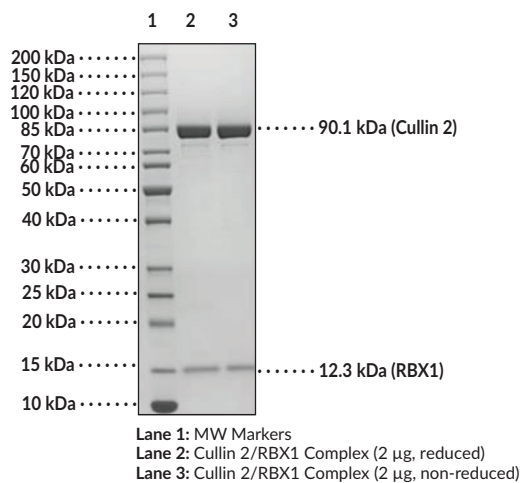
Activity: *batch specific* U/ml

Specific Activity: *batch specific* U/mg

Special Conditions: Rapid thaw with running water. Avoid repeated freeze-thaw cycles.

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

Image



SDS-PAGE Analysis of Cullin 2/RBX1 Complex.

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, 03/13/2026

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

The cullin 2 RING ligase (CRL2) complex is an E3 ubiquitin ligase complex.^{1,2} It is composed of a von Hippel-Lindau (VHL) E3 ligase, a substrate recognition receptor that interacts with elongin B and -C, cullin 2 as the scaffold, and RING box protein 1 (RBX1), which binds to ubiquitin-loaded E2.¹ NEDDylation of cullin 2 alters the conformation of cullin 2 and RBX1 to promote ubiquitin transfer to the CRL2 substrate.² The CRL2 complex is involved in germline development and viral infection.¹ A SNP in *CUL2*, the gene encoding cullin 2, is associated with an increased risk for Crohn's disease and ulcerative colitis.¹ Cayman's Cullin 2/RBX1 Complex (human, recombinant) can be used for enzyme activity assays.

References

1. Cai, W. and Yang, H. The structure and regulation of Cullin 2 based E3 ubiquitin ligases and their biological functions. *Cell Div.* **11**, 7 (2016).
2. Boh, B.K., Smith, P.G., and Hagen, T. Neddylation-induced conformational control regulates cullin RING ligase activity *in vivo*. *J. Mol. Biol.* **409**(2), 136-145 (2011).
3. Rivas, M.A., Beaudoin, M., Gardet, A., *et al.* Deep resequencing of GWAS loci identifies independent rare variants associated with inflammatory bowel disease. *Nat. Genet.* **43**(11), 1066-1073 (2011).

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