

# Cul5/Rnf7 [untagged]

## E3 Ligase

**Alternate Names:** Cul5 = VACM 1, Vasopressin activated calcium mobilizing receptor  
Rnf7 = Regulator of Cullins 2; ROC2, Rbx2, Sensitive to Apoptosis Gene; SAG

**Cat. No.** 63-1002-025  
**Lot. No.** 30165

**Quantity:** 25 µg  
**Storage:** -70°C

FOR RESEARCH USE ONLY

NOT FOR USE IN HUMANS



CERTIFICATE OF ANALYSIS Page 1 of 2

## Background

The enzymes of the ubiquitylation pathway play a pivotal role in a number of cellular processes including the regulated and targeted proteasome dependent degradation of substrate proteins. Three classes of enzymes are involved in the process of ubiquitylation; activating enzymes (E1s), conjugating enzymes (E2s) and protein ligases (E3s). Cullin-RING-Ligases (CRLs) are one largest class of ubiquitin E3 ligases and the enzymes of the NEDDylation pathway play a pivotal role in the activation of these, akin to ubiquitylation, the E1 activating enzyme (APP-BP1/UBA3 heterodimer) and the E2 conjugating enzymes (UBE2M or UBE2F) are involved in mammalian NEDDylation of the Cullin Ring Ligases (CRLs) (Meyer-Schaller *et al.*, 2009; Huang *et al.*, 2011; Morimoto *et al.*, 2003). Identification of the human Cullin1-5 genes were first described by Kipreos *et al.* (1996). Cullin RING ligases (CRL) comprise the largest sub-family of ubiquitin ligases which are activated by Neddylation. CRLs are involved in cell cycle regulation, DNA replication, DNA damage response (DDR). CRLs contain subunits including, a scaffold protein (cullin family protein), a Ring finger protein either Rbx1 (Cul1-4) or Rbx2 (Cul5) that binds a ubiquitin E2 Ube2M or Ube2F respectively (Sarikas, *et al.*, 2011; Skowyra *et al.*, 1997). Cul-5 has been shown to form a complex with the Ring

Continued on page 2

## Physical Characteristics

**Species:** human

**Source:** insect (Sf21)

**Quantity:** 25 µg

**Concentration:** 0.5 mg/ml

**Formulation:** 50 mM HEPES pH 7.5,  
150 mM sodium chloride,  
2 mM dithiothreitol, 10% glycerol

**Molecular Weight:**

Cul5: ~91.1 kDa; Rnf7: ~12.7 kDa

**Purity:** >85% by InstantBlue™ SDS-PAGE

**Stability/Storage:** 12 months at -70°C;  
aliquot as required

**Protein Sequences:** Please see page 2

## Quality Assurance

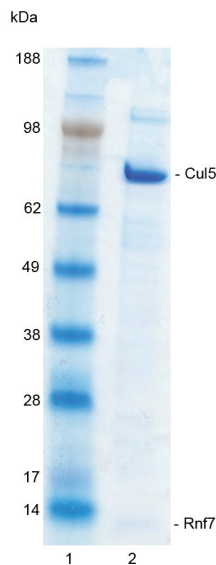
**Purity:**

4-12% gradient SDS-PAGE

InstantBlue™ staining

Lane 1: MW markers

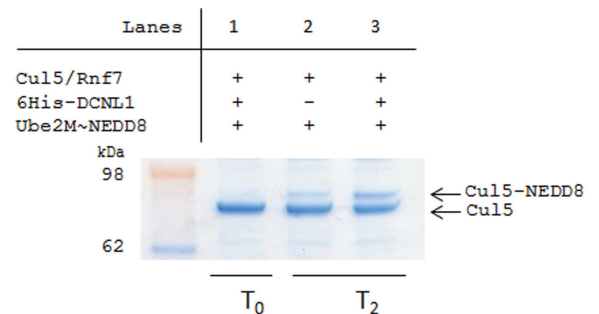
Lane 2: 1 µg Cul5/Rnf7



**Protein Identification:**

Confirmed by mass spectrometry.

**E3 Ligase Assay:** The activity of Cul5/Rnf7 was validated indirectly through its ability to act as a substrate for neddylation in the presence of the NEDD8 E3 ligase His-DCNL1 and thioester loaded His-Ube2M~NEDD8. Incubation of Cul5/Rnf7 and thioester loaded His-Ube2M~NEDD8 in the presence or absence of His-DCNL1 at 4°C was compared at two time points T<sub>0</sub> and T<sub>2</sub> minutes. Increased neddylation of the Cul5 subunit in the presence of His-DCNL1 was demonstrated.



www.ubiquigent.com  
Dundee, Scotland, UK

### ORDERS / SALES SUPPORT

**International:** +1-617-245-0020  
**US Toll-Free:** 1-888-4E1E2E3 (1-888-431-3233)  
**Email:** sales.support@ubiquigent.com

### UK HQ and TECHNICAL SUPPORT

**International:** +44 (0) 1382 381147 (9AM-5PM UTC)  
**US/Canada:** +1-617-245-0020 (9AM-5PM UTC)  
**Email:** tech.support@ubiquigent.com

Email services@ubiquigent.com for enquiries regarding compound profiling and/or custom assay development services.

© Ubiquigent 2013. Unless otherwise noted, Ubiquigent, Ubiquigent logo and all other trademarks are the property of Ubiquigent, Ltd.

**Limited Terms of Use:** For research use only. Not for use in humans or for diagnostics. Not for distribution or resale in any form, modification or derivative OR for use in providing services to a third party (e.g. screening or profiling) without the written permission of Ubiquigent, Ltd.

Lot-specific COA version tracker: v1.0.0

# Cul5/Rnf7 [untagged]

## E3 Ligase

**Alternate Names:** Cul5 = VACM 1, Vasopressin activated calcium mobilizing receptor  
Rnf7 = Regulator of Cullins 2; ROC2, Rbx2, Sensitive to Apoptosis Gene; SAG

**Cat. No.** 63-1002-025

**Lot. No.** 30165

**Quantity:** 25 µg

**Storage:** -70°C

FOR RESEARCH USE ONLY

NOT FOR USE IN HUMANS



CERTIFICATE OF ANALYSIS Page 2 of 2

## Background

Continued from page 1

finger protein Rbx2(Rnf7), the adaptor proteins Elongin B, Elongin C, and the SOCS(suppressors of cytokine signaling) box proteins to form an active CRL-5 E3 ligase (Okumura *et al.*, 2012; Sarikas *et al.*, 2011). Cul-5 also interacts with HSP90 and ErbB2. Cul-5 ubiquitylates ErbB2 – leading to its degradation – in the absence of the traditional adaptors Elongin B/C demonstrating the Elongin B/C independent E3 ligase activity of Cul-5/Rbx2 (Ehrlich *et al.*, 2009).

### References:

Ehrlich ES, Wang T, Luo K, Xiao Z, Niewiadomska AM, Martinez T, Xu W, Neckers L, Yu XF. (2009) Regulation of Hsp90 client proteins by a Cullin5-RING E3 ubiquitin ligase, *PNAS* **106**, 20330-20335.

Huang G, Kaufman AJ, Ramanathan Y, Singh B. (2011) SCCOR (DCUN1D1) promotes nuclear translocation and assembly of the neddylation E3 complex, *J Biol Chem* **286**, 10297-10304.

Meyer-Schaller N, Chou YC, Sumara I, Martin DD, Kurz T, Katheder N, Hofmann K, Berthiaume LG, Sicheri F, Peter M. (2009) The human Dcn1-like protein DCN13 promotes Cul3 neddylation at membranes, *PNAS* **106**, 12365-12370.

Morimoto M, Nishida T, Nagayama Y, Yasuda H. (2003) Neddylation of Cul1 is promoted by Roc1 as a Nedd8-E3 ligase and regulates its stability, *Biochem Biophys Res Commun* **301**, 392-398.

Okumura F, Matsuzaki M, Nakatsukasaka K, Kamura T. (2012) The Role of Elongin BC-Containing Ubiquitin Ligases. *Front Oncol* **2**, 1-13.

Sarikas A, Hartmann, T and Pan, ZQ (2011) The cullin protein family, *Genome Biology* **12**, 220.

Zhou W, Wei, W. and Sun, Y (2013) Genetically engineered mouse models for functional studies of SKP1-CUL1-F-box-protein (SCF) E3 ubiquitin ligases, *Cell Res* **23**, 599-619.

## Physical Characteristics

Continued from page 1

### Protein Sequence: Cullin 5

GGSMATSNNLLKNKGS~~LQ~~FEDKWFMRP I V L K L L  
RQESVTKQQWFDLFS~~VD~~HAVCLWDDKGP A K I  
HQALKEDI~~LE~~F I KQAQARVLSHQDDTALLKAY  
I~~VE~~WRKFF~~TQ~~CDILPKPFCQLEITLMGKQG  
SNKKS~~N~~VEDS I V R K L M L D T W N E S I F S N I K N  
RLQDSAMKLVHAERLGEAFDSQLVIGVRESYVN  
LCSNPEDK~~LQ~~IYRDNFEKAYLDSTERFYRTQA  
P S Y L Q Q N G V Q N Y M K Y A D A K L K E E E K R A L  
R Y L E T R R E C N S V E A L M E C C V N A L V T S F K E T I  
L A E C Q G M I K R N E T E K L H L M F S L M D K V P N G I  
E P M L K D L E E H I I S A G L A D M V A A A E T I T T D  
S E K Y V E Q L L T L F N R F S K L V K E A F Q D D P R F L  
T A R D K A Y K A V V N D A T I F K L E L P L K Q K G V G L K  
T Q P E S K C P E L L A N Y C D M L L R K T P L S K K L T  
S E E I E A K L K E V L L V L K Y V Q N K D V F M R Y H  
K A H L T R R L I L D I S A D S E I E E N M V E W L R E V G  
M P A D Y V N K L A R M F Q D I K V S E D L N Q A F K E M H  
K N N K L A L P A D S V N I K I L N A G A W S R S S E K V F V S  
L P T E L E D L I P E V E E F Y K K N H S G R K L H W H H L M  
S N G I I T F K N E V G Q Y D L E V T T F Q L A V L F A W N  
Q R P R E K I S F E N L K L A T E L P D A E L R R T L W S  
L V A F P K L K R Q V L L Y E P Q V N S P K D F T E G T L F S  
V N Q E F S L I K N A K V Q K R G K I N L I G R L Q L T T E R  
M R E E N E G I V Q L R I L R T Q E A I I Q I M K M R K  
K I S N A Q L O T E L V E I L K N M F L P Q K K M I K E Q  
I E W L I E H K Y I R R D E S D I N T F I Y M A

The residues underlined remain after cleavage and removal of the purification tag.

Cullin 5 (regular text): Start ***bold italics*** (amino acid residues 1-780)

Accession number: NP\_003469.2

Cullin5 [Dac tagged] /Rnf7 was cleaved with TEV protease [6His tagged] (to remove the Dac tag). The Dac tag and TEV protease [6His-tagged] were removed by capturing on amp sepharose and nickel resin respectively.

### Protein Sequence: Rnf7

**MAD**VEDGEETCALASHSGSSGSKSGGDKMF  
SLKKNNAVAMWSWDVECDTCAICRVQVM  
DACLRCAENKQEDCVVVWGECNHSFHNC  
MSLWVKQNNRCP~~LQ~~QDQWVVRIGK

Rnf7 (regular text): Start ***bold italics*** (amino acid residues 1-113)

Accession number: NP\_055060.1



www.ubiquigent.com  
Dundee, Scotland, UK

### ORDERS / SALES SUPPORT

**International:** +1-617-245-0020  
**US Toll-Free:** 1-888-4E1E2E3 (1-888-431-3233)  
**Email:** sales.support@ubiquigent.com

### UK HQ and TECHNICAL SUPPORT

**International:** +44 (0) 1382 381147 (9AM-5PM UTC)  
**US/Canada:** +1-617-245-0020 (9AM-5PM UTC)  
**Email:** tech.support@ubiquigent.com

Email services@ubiquigent.com for enquiries regarding compound profiling and/or custom assay development services.

© Ubiquigent 2013. Unless otherwise noted, Ubiquigent, Ubiquigent logo and all other trademarks are the property of Ubiquigent, Ltd.

Limited Terms of Use: For research use only. Not for use in humans or for diagnostics. Not for distribution or resale in any form, modification or derivative OR for use in providing services to a third party (e.g. screening or profiling) without the written permission of Ubiquigent, Ltd.

Lot-specific COA version tracker: v1.0.0