

# Cul3/Rbx1 [untagged]

## E3 Ligase

Alternate Names: Cul3 = KIAA0617  
Rbx1=HRT1, Regulator of cullins 1, Ring finger protein 75, RNF75, ROC1, ZYP protein

Cat. No. 63-1003-025  
Lot. No. 30206

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). The human Cullin 1-5 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; Skowrya *et al.*, 1997). Cul3 expression in human fibroblasts is induced by phorbol 12-myristate 13-acetate (PMA) and

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:  
Cul3: ~89.0 kDa; Rbx1: ~12.3 kDa

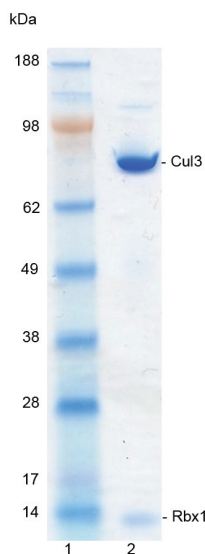
Purity: >95% by InstantBlue™ SDS-PAGE

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

Protein Sequences: Please see page 2

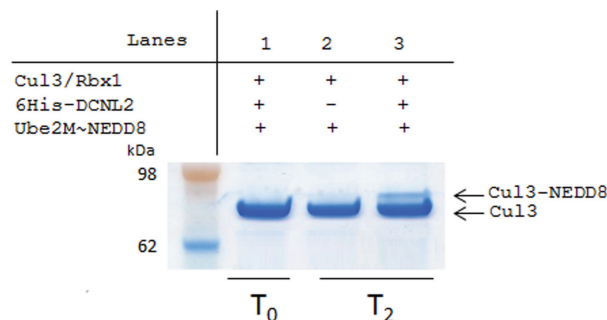
### Quality Assurance

Purity:  
4-12% gradient SDS-PAGE  
InstantBlue™ staining  
Lane 1: MW markers  
Lane 2: 1 µg Cul3/Rbx1



Protein Identification:  
Confirmed by mass spectrometry.

**E3 Ligase Assay:** The activity of Cul3/Rbx1 was validated indirectly through its ability to act as a substrate for neddylation in the presence of the NEDD8 E3 ligase His-DCNL2 and thioester-loaded His-Ube2M~NEDD8. Incubation of Cul3/Rbx1 and thioester loaded His-Ube2M~NEDD8 in the presence or absence of His-DCNL2 at 4°C was compared at two time points T<sub>0</sub> and T<sub>2</sub> minutes. Neddylation of the Cul3 subunit in the presence of His-DCNL2 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: v1.0.1

# Cul3/Rbx1 [untagged]

## E3 Ligase

**Alternate Names:** Cul3 = KIAA0617  
Rbx1=HRT1, Regulator of cullins 1, Ring finger protein 75, RNF75, ROC1, ZYP protein

**Cat. No.** 63-1003-025  
**Lot. No.** 30206

**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

suppressed by salicylate (Du *et al.*, 1998). The Cul3/Kelch like 9 (KLHL9)/Kelch like 13 (KLHL13) complex is an E3 ligase that controls the behavior of Aurora B on mitotic chromosomes and thereby coordinating mitotic progression and completion of cytokinesis (Sumara *et al.*, 2007). Interaction of Cul3 with Kelch like 7 (KLHL7) leads to the ubiquitylation of the dopamine receptor D4 (DRD4) (Rondou *et al.*, 2008).

### References:

Du M, Sansores-Garcia L, Zu Z, Wu KK, (1998) Cloning and expression analysis of a novel salicylate suppressible gene, Hs-CUL-3, a member of cullin/Cdc53 family. *J Biol Chem* **273**, 24289-24292.

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

Kipreos ET, Lander LE, Wing JP, He WW, Hedgecock EM (1996) cul-1 is required for cell cycle exit in *C. elegans* and identifies a novel gene family, *Cell* **85**, 829-839.

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 DCNL3 promotes Cul3 neddylation at membranes, *Proc Natl Acad Sci U S A* **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.

Rondou P, Haegeman G, Vanhoenacker P, Van Craenenbroeck K, (2008) BTB protein KLHL12 targets the dopamine D4 receptor for ubiquitination by a Cul3-based E3 ligase. *J Biol Chem* **283** 11083-11096.

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

Skowyra D, Craig KL, Tyers M, Elledge SJ, Harper JW (1997) F-box proteins are receptors that recruit phosphorylated substrates to the SCF ubiquitin-ligase complex, *Cell* **91**, 209-219.

Sumara I, Quadroni M, Frei C, Olma MH, Sumara G, Ricci R, Peter MA (2007) Cul3-based E3 ligase removes Aurora B from mitotic chromosomes, regulating mitotic progression and completion of cytokinesis in human cells. *Dev Cell* **12**, 887-900.

## Physical Characteristics

Continued from page 1

### Protein Sequence: Cullin 3

GGSMSNLSKGTGSRKDTKMRIRAFPMTM  
DEKYVNSIWDLLKNAIQEIQRKNNSSGLS  
FEELYRNAYTMVLHKKHGEKLYTGLREVVTE  
HLINKVREDVLSLNNNFLTQTLNQAWND  
HQ TAMVMIRDILMYMDRVYVQQNNVENVNL  
GLIIFRDQVVRYGCI RDHLRQTLLEDMI  
ARERKGEVVDRGAIRNACQMLMILGLEGRS  
VYEEDFEAPFLEMSAEFFQMESQKFLAEN  
SASVYIKKVEARINEEIERVMHCLDKSTEEP  
IVKVVVERELISKHMKTIVEMENSGLVHM  
LKNGKTEDLGCYKLF SRVPNGLKTMCECMS  
SYLREQGKALVSEEGEGKNPVDYIQGLDLK  
SRFDRFLLESFNNDRLFQTIAGDFEYFLN  
LNSRSPEYLSLFI DDKLLKKGVKGLTEQE  
VETILDKAMVLF RFMQEKDVFERYYKQHLAR  
RLLTNKSVDSDDEKNNMISKLKTECGC  
QFTSKLEGMFRDMSISNTTMDEFHQHLQAT  
GVSLGGVDLTVRVLT TGWPTQSATPKC  
NIPPAPRHAFEIFRRFYLAHSGRQLTLQH  
HMG SADLNATFYGPVKKEDGSEVGVG  
GAQVTGSNTRKHILQVSTFQMTILMLFN  
NREKYTFEEIQQETDIPERELVRALQS  
LACGKPTQRVLTKEPKSKEIENGHIFT  
VNDQFTSKLHRVKIQTVAAKQGESDPERKET  
RQKVDDDRKHEIEAAIVRIMKSRKMQHNVL  
VAEVTQQLKARFLPSPVVIKKRIEGLIEREY  
LARTPEDRKVYTYVA

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

Cullin3 (regular text): Start **bold italics** (amino acid residues 1-768)

Accession number: NP\_003581.1

Cullin3 [Dac-tagged]/Rbx1 was cleaved with TEV protease [6His-tagged]. The Dac tag and TEV protease [6His-tagged] were removed by capturing on amp sepharose and nickel resin respectively.

### Protein Sequence: Rbx1

MAAAMDVDTPSGTNSGAGKKRFVEVKKW  
NAVALWAWD I VVDNCAICRNHIMDLCEIC  
QANQASATSEECTVAWGVCNHFHFHCISR  
WLKTRQVCPLDNREWEFQKYGH

Rbx1 (regular text): Start **bold italics** (amino acid residues 1-115)

Accession number: NP\_055063.1



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: v1.0.1