

# DCNL2 [6His-tagged]

## E3 Ligase

Alternate Names: C13orf17 protein, FLJ10704, FLJ20092, DCUN1D2

Cat. No. **63-2004-025**

Lot. No. **30161**

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 NEDDylation pathway play a pivotal role in the activation of the largest class of ubiquitin E3 ligases called Cullin-RING-Ligases (CRLs). Akin to ubiquitylation three classes of enzymes are involved in the process of mammalian NEDDylation; E1 activating enzyme (APP-BP1/UBA3 heterodimer), E2 conjugating enzymes (UBE2M or UBE2F) and E3 ligases the defective in Cul neddylation 1 domain-containing proteins (DCUN1D1-5) (Meyer-Schaller *et al.*, 2009; Huang *et al.*, 2011). There are 5 human DCUN1D1-5 proteins are also named defective in Cul neddylation 1 like proteins (DCNL1-5) (Meyer-Schaller *et al.*, 2009). Cloning of DCNL2 was first described by Kurz *et al.* (2005) and Lamesch *et al.* (2007). The DCNLs have distinct amino-terminal domains, but share a conserved C-terminal potentiating neddylation (PONY) domain (Kurz *et al.*, 2008). It has been determined that the interaction between the DCNLs and Cul1 occurs through the PONY domain and the Winged Helix DNA binding domain (WHB) respectively (Kurz *et al.*, 2008; Scott *et al.*, 2011). Pairwise analysis of 30 combinations of the five DCNL PONY domains and six cullin WHB subdomains by isothermal titration calorimetry have all shown interaction albeit with differing affinities (Monda *et al.*, 2013).

Continued on page 2

### Physical Characteristics

**Species:** human

**Source:** *E. coli*

**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:** ~33.3 kDa

**Purity:** >98% by InstantBlue™ SDS-PAGE

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

### Protein Sequence:

**MGSSHHHHHSSGLEVLFGQPGSPPEFPGVDMH**  
KLKSSQKDKVRQFMACTQAGERTAIYCLTQNE  
WRLDEATDSFFQNPDSLHRESMRNAVVDKK  
KLERLYGRYKDPQDENKIGVDGIQQFCFCD  
DLSLDPASISVLVIAWKFRRAATQCEFSRKE  
FLDGMTELGCDSEKLLKALLPRLEQELKDTAK  
FKDFYQFTTFAKNPQGKGLDLEMAVAYWKL  
LSGRFKFLDLWNTFLMEHHKRSIPRDTWNL  
DFGNMIADDMSNYDEEGAWPVLIDDFVEYAR  
PVVTGGKRSFLF

Tag (**bold text**): N-terminal His

Protease cleavage site: PreScission™ (LEVLFGQ▼GP)

DCNL2 (regular text): Start **bold italics** (amino acid residues 1-259)

Accession number: NP\_001014305.1

### Quality Assurance

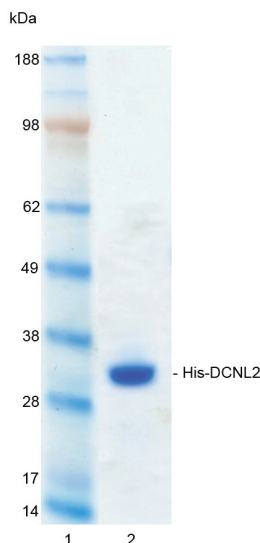
#### Purity:

4-12% gradient SDS-PAGE

InstantBlue™ staining

Lane 1: MW markers

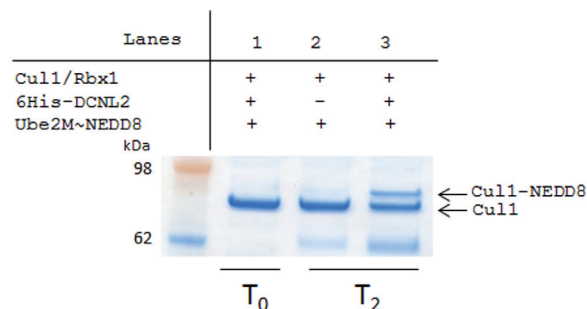
Lane 2: 1 µg His-DCNL2



#### Protein Identification:

Confirmed by mass spectrometry.

**E3 Ligase Assay:** The activity of His-DCNL2 was validated through its ability to enhance the neddylation of Cul1/Rbx1 acting as a substrate in the presence of the thioester-loaded His-Ube2M~NEDD8. Incubation of Cul1/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. Increased neddylation of the Cul1 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: v1.0.0

# DCNL2 [6His-tagged]

## E3 Ligase

Alternate Names: C13orf17 protein, FLJ10704, FLJ20092, DCUN1D2

Cat. No. **63-2004-025**

Lot. No. **30161**

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

### References:

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

Kurz T, Chou YC, Willems AR, Meyer-Schaller N, Hecht ML, Tyers M, Peter M, Sicheri F. (2008) Dcn1 functions as a scaffold-type E3 ligase for cullin neddylation, *Mol Cell* **29**, 23-35.

Kurz T, Ozlü N, Rudolf F, O'Rourke SM, Luke B, Hofmann K, Hyman AA, Bowerman B, Peter M. (2005) The conserved protein DCN-1/Dcn1p is required for cullin neddylation in *C. elegans* and *S. cerevisiae*, *Nature* **435**, 1257-1261.

Lamesch P, Li N, Milstein S, Fan C, Hao T, Szabo G, Hu Z, Venkatesan K, Bethel G, Martin P, Rogers J, Lawlor S, McLaren S, Dricot A, Borick H, Cusick ME, Vandenhaute J, Dunham I, Hill DE, Vidal M. (2007) hORFeome v3.1: a resource of human open reading frames representing over 10,000 human genes, *Genomics* **89**, 307-315.

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.

Monda J,K,Scott DC, Miller DJ, Lydeard J, King D, Harper JW, Bennett EJ, Schulman BA. (2013) Structural Conservation of Distinctive N-terminal Acetylation-Dependent Interactions across a Family of Mammalian NEDD8 Ligation Enzymes, *Structure* **21**, 42-53.

Scott D.C, Monda JK, Bennett EJ, Harper JW, Schulman B.A. (2011) N-terminal acetylation acts as an avidity enhancer within an interconnected multiprotein complex, *Science* **334**, 674-678.



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