UBE2F (NCE2) [untagged]

E2 - NEDD8 Conjugating Enzyme

Alternate Names: NEDD8 conjugating enzyme, MGC18120, NCE2

Cat. No.	62-0025-100	Quantity:
Lot. No.	30120	Storage:
FOR RESEARCH USE ONLY		

100 µg -70°C

NOT FOR USE IN HUMANS

CERTIFICATE OF ANALYSIS

UBIQUIGENT

Background

The enzymes of the NEDDylation pathway play a pivotal role in a number of cellular processes including the indirect regulation and targeting of substrate proteins for proteasomal degradation. Three classes of enzymes are involved in the process of NEDDylation; the ubiquitin-like activating enzyme APP-BP1/Uba3 (E1), the ubiquitin-like conjugating enzymes (E2s) and protein ligases (E3s). UBE2F is a member of the E2 conjugating enzyme family and the human gene was first described by Huang et al. (2009). UBE2F acts as a NEDD8 conjugating enzyme both in vitro and in vivo. UBE2F accepts the ubiquitin-like protein NEDD8 from the Uba3-NAE1 (APP-BP1/Uba3) E1 complex and catalyzes its covalent attachment to other proteins. The specific interaction of UBE2F with the E3 ubiquitin ligase RBX2, but not RBX1, suggests that the RBX2-UBE2F complex NEDDylates specific target proteins such as CUL5, a component of one of the many Cullin Ring Ligases (CRLs) (Huang et al., 2009).

Reference:

Huang DT, Avrault O, Hunt HW, Taherbhov AM, Duda DM, Scott DC, Borg LA, Neale G, Murray PJ, Roussel MF, Schulman BA (2009) E2-RING expansion of the NEDD8 cascade confers specificity to cullin modification. Mol Cell 33, 483-95.

Physical Characteristics

Species: human

Source: E. coli expression

Quantity: 100 µg

Concentration: 1 mg/ml

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

Molecular Weight: ~23 kDa

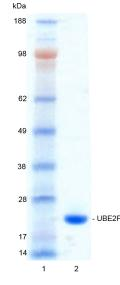
Purity: >95% by InstantBlue™ SDS-PAGE

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

Quality Assurance

Purity:

4-12% gradient SDS-PAGE InstantBlue[™] staining Lane 1: MW markers Lane 2: 1 µg UBE2F



Protein Sequence:

GSHMASMTGGQQMGRGS**M**LTLASKLKRDDG LKGSRTAATASDSTRRVSVRDKLLVKEVAEL EANLPCTCKVHFPDPNKLHCFOLTVTPDEGYY QGGKFQFETEVPDAYNMVPPKVKCLTKIWHP NITETGEICLSLLREHSIDGTGWAPTRTLKDV VWGLNSLFTDLLNFDDPLNIEAAEHHLRDKED FRNKVDDYIKRYAR

The residues underlined remain after cleavage and removal of the purification tag. UBE2F (regular text): Start **bold italics** (amino acid residues 1-185)

Accession number: NP 542409

Protein Identification: Confirmed by mass spectrometry.

E2-NEDD8 Thioester Loading Assay:

The activity of UBE2F was validated by loading E1 (APP-BP1/Uba3) activated NEDD8 onto the active cysteine of the UBE2F E2 enzyme via a transthiolationreaction. Incubation of the APP-BP1/ Uba3 and UBE2F enzymes in the presence of NEDD8 and ATP at 30°C was compared at two time points, T_0 and T_{10} minutes. The sensitivity of the NEDD8/ UBE2F thioester bond to the reducing agent DTT was confirmed.



ORDERS / SALES SUPPORT US Toll-Free: 1-888-4E1E2E3 (1-888-431-3233) International: +1-617-245-0003 Email: sales.support@ubiquigent.com

UK HQ and TECHNICAL SUPPORT International: +1-617-245-0020

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