UBE2F (NCE2) [6His-tagged]

E2 - NEDD8 Conjugating Enzyme

Alternate Names: NEDD8 conjugating enzyme, MGC18120, NCE2

62-0023-100 Quantity: 100 µg Cat. No. Lot. No. 1372 Storage: -70°C

FOR RESEARCH USE ONLY NOT FOR USE IN HUMANS



CERTIFICATE OF ANALYSIS

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 proteosomal degradation. Three classes of enzymes are involved in the process of NEDDylation; the ubiquitin-like activating enzyme APP-BP1/Uba3 (E1), the ubiquitinlike conjugating enzymes (E2s) and protein ligases (E3s). UBE2F is a member of the ubiquitin-like 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 ubiguitin-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, Ayrault O, Hunt HW, Taherbhoy 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: ~25 kDa

Purity: >98% by InstantBlue™ SDS-PAGE

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

aliquot as required

Protein Sequence:

MGSSHHHHHHSSGLVPRGSHMASMTG GQQMGRGS**M**LTLASKLKRDDGLKGSR TAATASDSTRRVSVRDKLLVKEVAELEAN LPCTCKVHFPDPNKLHCFQLTVTPDEGYYQG GKFQFETEVPDAYNMVPPKVKCLTKIWHP NITETGEICLSLLREHSIDGTGWAPTRTLKDV VWGLNSLFTDLLNFDDPLNIEAAEHHLRDKED **FRNKVDDYIKRYAR**

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

Protease cleavage site: Thrombin (LVPR▼GS) UBE2F (regular text): Start bold italics (amino acid

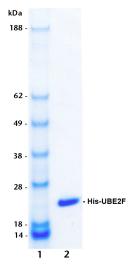
residues 1-185)

Accession number: NP_542409

Quality Assurance

Purity:

4-12% gradient SDS-PAGE InstantBlue™ staining lane 1: MW markers lane 2: 1 µg His-UBE2F



Protein Identification:

Confirmed by mass spectrometry.

E2-NEDD8 Thioester Loading Assay:

The activity of His-UBE2F was validated by loading E1 APP-BP1/Uba3 activated NEDD8 onto the active cysteine of the His-UBE2F E2 enzyme via a transthiolation reaction. Incubation of the APP-BP1/Uba3 and His-UBE2F enzymes in the presence of NEDD8 and ATP at 30°C was compared at two time points, T₀ and T₁₀ minutes. Sensitivity of the NEDD8/ His-UBE2F thioester bond to the reducing agent DTT was confirmed.



Dundee, Scotland, UK

ORDERS / SALES SUPPORT

International: +1-617-245-0003

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 2011**. 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