

His₆-UBE2M

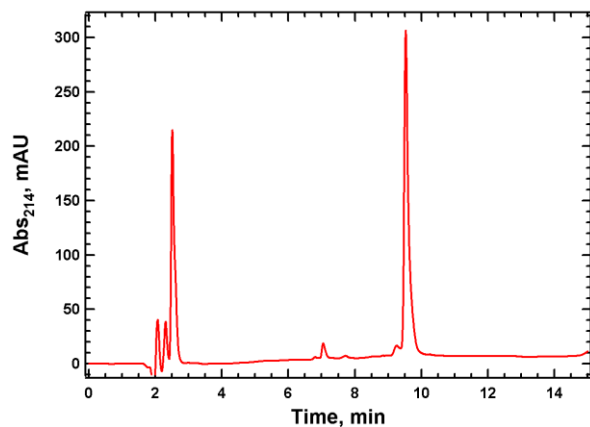
Cat. # UB217H

Background: UBE2M (Ubc12) is an E2 that is involved in the conjugation of the ubiquitin-like protein (Ubl) NEDD8 to target substrates (NEDDylation) along with the APPBP1/Uba3 activating enzyme. NEDDylation is required for cell cycle progression. In addition, NEDD8 is conjugated to cullin family members which are a part of SCF ubiquitin E3 ligase complexes, thereby activating them.

Application: Ubiquitin ligation reactions

Product Information

Organism	Human, recombinant; Accession No. P61081
Purity:	≥ 90% by RP-HPLC
Molecular Weight:	21894.2 Da by MS (calculated 21894)
Tag	His ₆
Physical State:	Liquid, 50 mM Tris, pH 7.5; 150 mM NaCl; 10 mM DTT; 10% glycerol
Quantity:	20 μL or 75 μL of a 40 μM solution (0.8 and 3 nmoles, respectively)
Solubility:	>3 mg/mL
Storage:	-80° C. Avoid repeated freeze/thaw cycles



RP-HPLC

References

- Huang, D.T. et al. E2-RING expansion of the NEDD8 cascade confers specificity to cullin modification. *Mol Cell* **33**, 483-95 (2009).
- Huang, D.T. et al. Structural basis for recruitment of Ubc12 by an E2 binding domain in NEDD8's E1. *Mol Cell* **17**, 341-50 (2005).
- Wada, H., Yeh, E.T. & Kamitani, T. Identification of NEDD8-conjugation site in human cullin-2. *Biochem Biophys Res Commun* **257**, 100-5 (1999).
- Wada, H., Yeh, E.T. & Kamitani, T. A dominant-negative UBC12 mutant sequesters NEDD8 and inhibits NEDD8 conjugation in vivo. *J Biol Chem* **275**, 17008-15 (2000).

All products are for research use only • not intended for human or animal diagnostic or therapeutic uses
Copyright © 2009 LifeSensors, Inc. All Rights Reserved