

UBE2E3

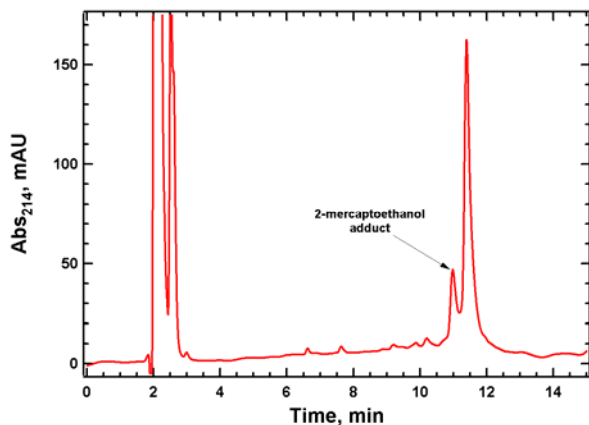
Cat. # UB213

Background: Ubiquitin-conjugating enzyme E2 E3 (UBE2E3 alias UbcH9) is a protein that in humans is encoded by the *UBE2E3* gene. The deduced 207-amino acid protein shares over 94% amino acid identity with the UBC domains of class III E2s, UbcH6, UBE2E2, UbcM2, UbcM3, and UbcD2. But the N-terminal extension exhibited little homology among these, except for UbcM2, which showed 100% identity, and which is thought to be a mouse counterpart. The UBE2E3 form a thioester bond with ubiquitin (Ub) in an E1-dependent manner and mediates the transfer of ubiquitin from a ubiquitin-activating enzyme (E1) to a substrate protein or E3 ligase (2)

Application: Ubiquitin ligation reactions

Product Information

Organism	Human, recombinant; Accession No. Q969T4
Purity:	≥ 95% by RP-HPLC
Molecular Weight:	22,913 Da by MS (calculated 22,912.7)
Tag	none
Physical State:	Liquid, 20 mM Tris, pH 7.4; 150 mM NaCl; 10 mM DTT; 10% glycerol
Quantity:	20 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

- Ito, K., Kato, S., Matsuda, Y., Kimura, M., and Okano, Y. (1999) cDNA cloning, characterization, and chromosome mapping of UBE2E3 (alias UbcH9), encoding an N-terminally extended human ubiquitin-conjugating enzyme. *Cytogenet Cell Genet.* **84**, 99-104.

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