

UBE2E2

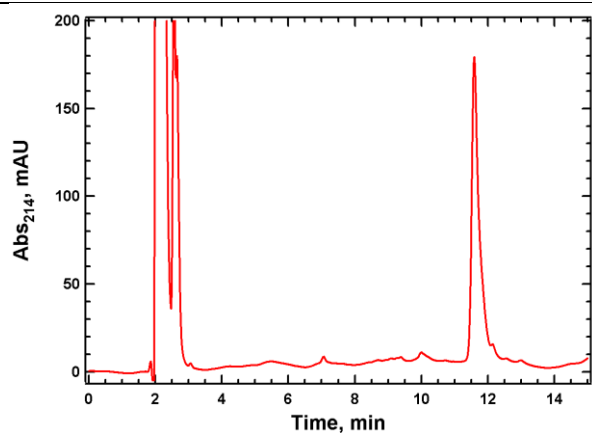
Cat. # UB212

Background: Ubiquitin-conjugating enzyme E2 (UBE2E2) is a protein that in humans is encoded by the *UBE2E2* gene. Amino acid sequence within the UBC domain of UBE2E2 shares over 90% identity with human UbcH6, mouse UbcM2, and *Drosophila* UbcD2, whereas the N-terminal region shows little amino acid sequence similarity with known proteins. The UBE2E2 protein forms a thioester bond with ubiquitin in an E1-dependent manner and mediates the transfer of ubiquitin from a ubiquitin-activating enzyme (E1) to a substrate protein or E3 ligase (1).

Application: Ubiquitin conjugation reactions

Product Information

Organism	Human, recombinant; Accession No. Q96LR5
Purity:	≥ 95% by RP-HPLC
Molecular Weight:	22,255.3 Da by MS (calculated 22,255.1)
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

Kimura, M., Hattori, T., Matsuda, Y., Yoshioka, T., Sumi, N., Umeda, Y., Nakashima, S., and Okano, Y. (1997) cDNA cloning, characterization, and chromosome mapping of UBE2E2 encoding a human ubiquitin-conjugating E2 enzyme. *Cytogenet Cell Genet.* **78**, 107-11.

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