

UBE2T

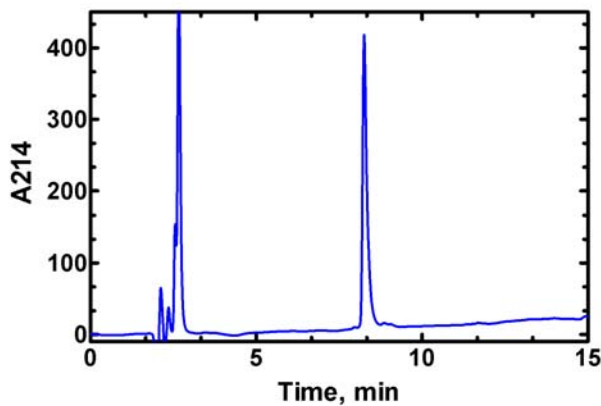
Cat. # UB219

Background: UBE2T can mediate the transfer of ubiquitin from ubiquitin-activating enzyme E1 to a substrate protein or E3 ligase. This E2 enzyme has been shown to be essential for the Fanconi anemia pathway. UBE2T conjugating activity is involved with the monoubiquitylation of FANCD2, co-functioning with the E3 ligase FANCL.

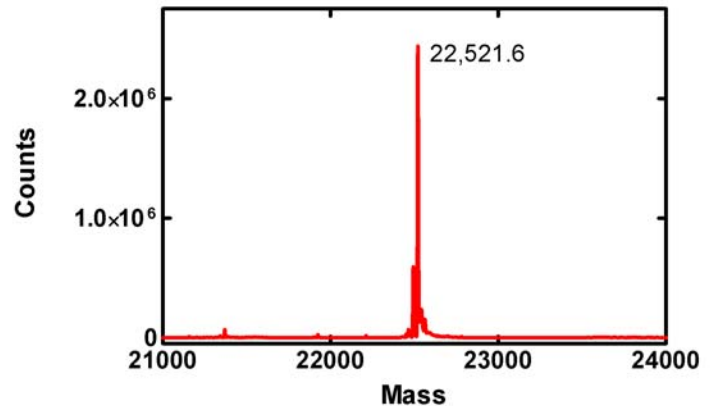
Application: Ubiquitin ligation reactions

Product Information

Organism	Human, Accession No. Q9NPD8
Source	Recombinant, <i>E. coli</i>
Purity:	≥ 80% by RP-HPLC
Molecular Weight:	22,520.9 Da
Tag	none
Physical State:	Liquid, 25 mM Tris, pH 7.4; 150 mM NaCl; 10 mM Dtt; 10% glycerol
Quantity:	3 nmol (75 μL of a 40 μM solution)
Solubility:	>3 mg/mL
Storage:	-80° C. Avoid repeated freeze/thaw cycles



RP-HPLC



Deconvoluted mass spectrum

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