

Background

UBE2T mediates the transfer of ubiquitin from a ubiquitin-activating enzyme (E1) to a substrate protein or an E3 ligase. This E2 enzyme is essential for the Fanconi anemia pathway. UBE2T's conjugating activity is specifically involved in the monoubiquitination of FANCD2, in cooperation with the E3 ligase FANCL.

Alternate Names

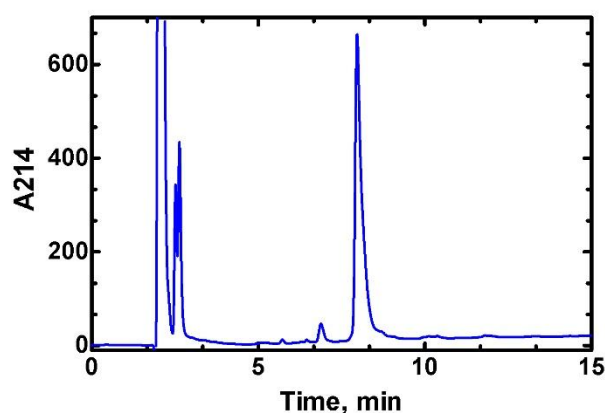
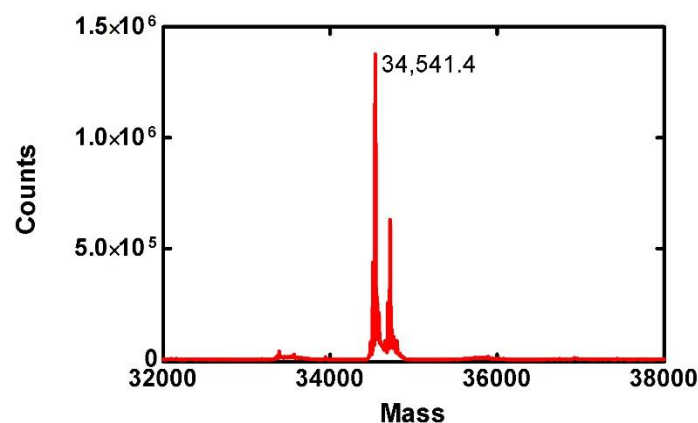
Ubiquitin-Conjugating Enzyme E2 T, Cell Proliferation-Inducing Gene 50 Protein, HSPC150, FANCT, PIG50

Application(s)

Ubiquitin ligation reactions

Product Specifications

Tag	His ₆ -SUMO
Purity	> 80% by RP-HPLC
Molecular Weight	34,540.3 Da
Quantity	20 µl or 75 µl of a 40 µM solution (0.8 and 3 nmoles, respectively)
Species	Human, Accession No. Q9NPD8
Expression System	<i>E. Coli</i>
Physical State	Liquid
Buffer	25 mM Tris, pH 7.4; 150 mM NaCl; 10 mM DTT; 10% glycerol
Solubility	> 3 mg/ml
Stability & Storage	1 year at -80° C. Avoid repeated freeze/thaw cycles

Product QC**RP-HPLC****Deconvoluted mass spectrum****References**

1. Machida YJ., et al., Mol Cell. 2006;23(4):589-96.
2. Rickman KA., et al., Cell Rep. 2015;12(1):35-41.

All products are for research use only • Not intended for human or animal diagnostic or therapeutic uses
Copyright © 2025 LifeSensors, Inc. All Rights Reserved