

His₆-UBE2R1

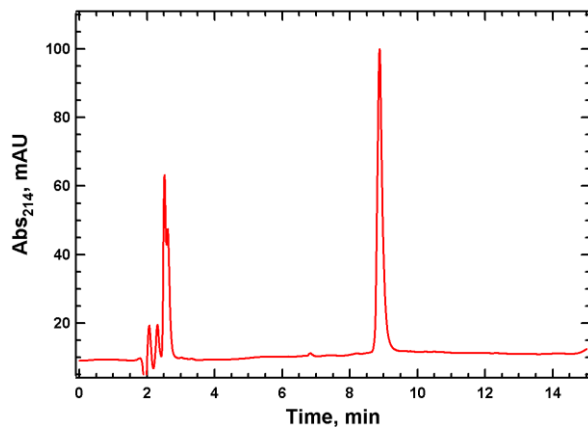
Cat. # UB208H

Background: 236 amino acid product of the CDC34 gene. It controls cellular proliferation through regulation of p27/Kip1 protein levels. Blocking CDC34 activity enhances the anti-myeloma effects of Bortezomib.

Application: Ubiquitin ligation reactions

Product Information

Organism	Human, recombinant; Accession No. P49427
Purity:	≥ 95% by RP-HPLC
Molecular Weight:	27,730.4 Da by MS (calculated 27,730.8)
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**

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