

## UBE2C

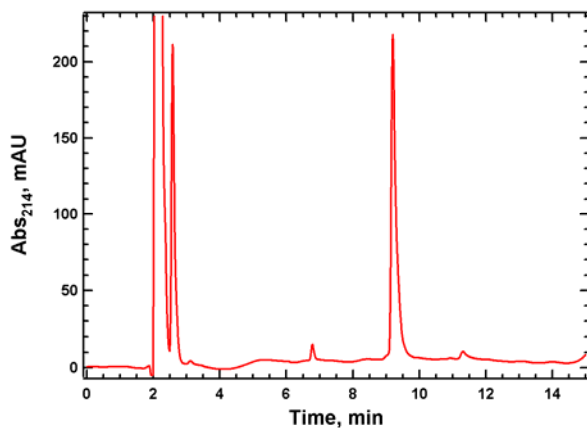
Cat. # UB222

**Background:** Ube2C (UbcH10) is an E2 conjugating enzyme and mediates the transfer of ubiquitin from an ubiquitin-activating enzyme (E1) to a substrate protein or E3 ligase. Ube2C interacts with the Anaphase Promoting Complex (APC) and as a member of this complex functions to ubiquitylate substrates and target these substrates, which include several cyclins, for degradation<sup>1</sup>. As such UbcH10 plays a critical role in cell cycle progression. Recently UbcH10 has been shown to play a key role in tumorigenesis<sup>2,3</sup>.

**Application:** Ubiquitin ligation reactions

### Product Information

<b>Organism</b>	Human, recombinant; Accession No. Q00762
<b>Purity:</b>	≥ 95% by RP-HPLC
<b>Molecular Weight:</b>	19,652.6 Da by MS (calculated 19,652.3)
<b>Tag</b>	none
<b>Physical State:</b>	Liquid, 25 mM Tris, pH 7.4; 150 mM NaCl; 10 mM DTT; 10% glycerol
<b>Quantity:</b>	20 or 75 µL of a 40 µM solution (0.8 and 3 nmoles, respectively)
<b>Solubility:</b>	>3 mg/mL
<b>Storage:</b>	-80° C. Avoid repeated freeze/thaw cycles



**RP-HPLC**

### References

1. Townsley FM, Aristarkhov A, Beck S, Hershko A, Ruderman JV. Dominant-negative cyclin-selective ubiquitin carrier protein E2-C/UbcH10 blocks cells in metaphase. *Proc Natl Acad Sci U S A*. **94**,2362-67 (1997).
2. van Ree JH, Jeganathan KB, Malureanu L, van Deursen JM. Overexpression of the E2 ubiquitin-conjugating enzyme UbcH10 causes chromosome missegregation and tumor formation. *J Exp Med*. **207**,83-100 (2010).
3. Chen SM, Jiang CY, Wu JY, Liu B, Chen YJ, Hu CJ, Liu XX. RNA interference-mediated silencing of UBCH10 gene inhibits colorectal cancer cell growth in vitro and in vivo. *Clin Exp Pharmacol Physiol*. DOI 10.1111/j.1440-1681.2009.05348.x (2009).

All products are for research use only • not intended for human or animal diagnostic or therapeutic uses  
Copyright © 2010 LifeSensors, Inc. All Rights Reserved