

## UBE2Q2

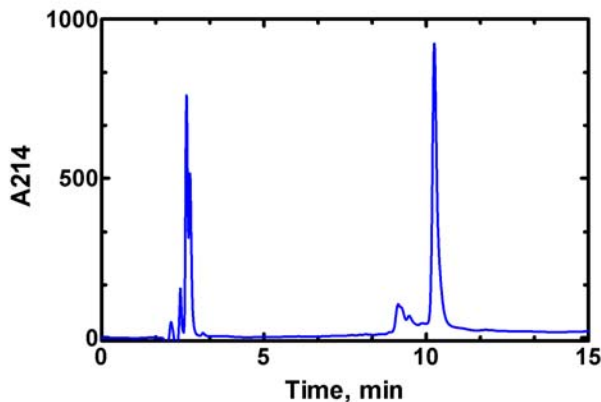
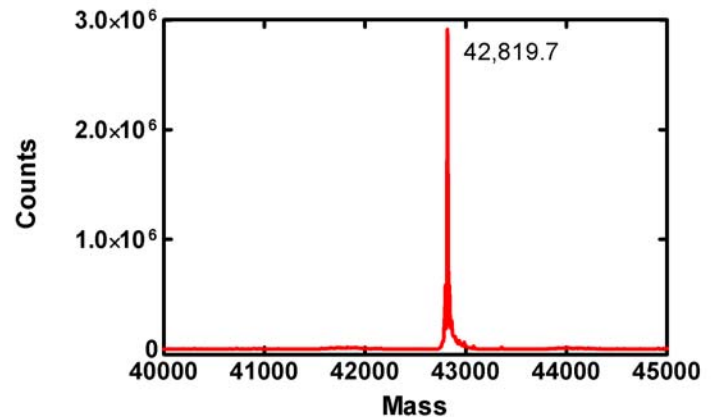
Cat. # UB225

**Background:** UBE2Q2, a member of the Ubc2 family of ubiquitin-conjugating E2 enzymes, was originally discovered as a differentially expressed gene in the endometrium of rabbits<sup>1</sup>. Further studies have shown that the enzyme is involved in cell cycle regulation and inhibition of its activity leads to mitotic arrest in the presence of microtubule inhibiting agents<sup>2</sup>. UBE2Q2 message and protein levels are elevated in human head and neck squamous cell carcinomas<sup>3</sup>.

**Application:** Ubiquitin ligation reactions

### Product Information

<b>Organism</b>	Human, Accession No. Q8WVN8
<b>Source</b>	Recombinant, <i>E. coli</i>
<b>Purity:</b>	≥ 80% by RP-HPLC
<b>Molecular Weight:</b>	42,818.4 Da
<b>Tag</b>	none
<b>Physical State:</b>	Liquid, TBS, 10% glycerol
<b>Quantity:</b>	3 nmol (75 µL @ 40 µM)
<b>Solubility:</b>	>3 mg/mL
<b>Storage:</b>	-80° C. Avoid repeated freeze/thaw cycles

**RP-HPLC****Deconvoluted mass spectrum**

### References

1. Melner, M.H., Haas, A.L., Klein, J.M., Brash, A.R., Boeglin, W.E., Nagdas, S.K., Winfrey, V.P., and Olson, G.E. Demonstration of ubiquitin thiolester formation of UBE2Q2 (UBCi), a novel ubiquitin-conjugating enzyme with implantation site-specific expression. *Biol Reprod.* **75**:395-406 (2006).
2. Banerjee, S., Brooks, W.S., and Crawford, D.F. Inactivation of the ubiquitin conjugatin enzyme UBE2Q2 causes a prophase arrest and enhanced apoptosis in response to microtubule inhibiting agents. *Oncogene* **26**:6509-6517 (2007).
3. Maeda, H., Miyajima, N., Kano, S., Tsukiyama, T., Okumura, F., Fukuda, S., and Hatakeyama, S. Ubiquitin-conjugating enzyme UBE2Q2 suppresses cell proliferation and is down-regulated in recurrent head and neck cancer. *Mol. Cancer Res.* **7**:1553-1562 (2009).

All products are for research use only • not intended for human or animal diagnostic or therapeutic uses

Copyright © 2009 LifeSensors, Inc. All Rights Reserved