

## Ubiquitin, K48R

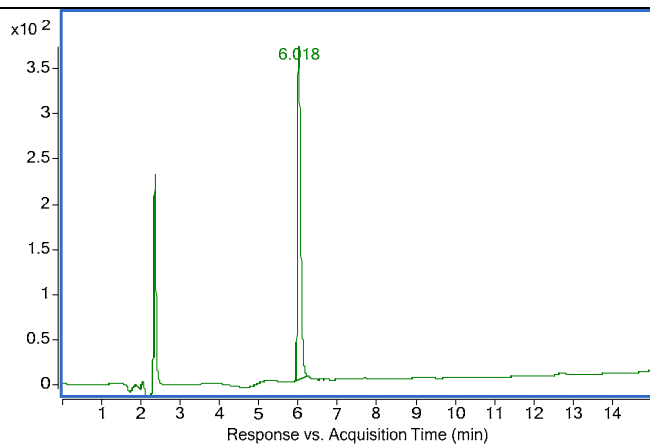
Cat. # SI217

**Background:** This tag-free recombinant form of human ubiquitin is engineered to have a single substitution of arginine for lysine at position 48. This mutant ubiquitin cannot form polyubiquitin chains through the lysine 48 linkage.

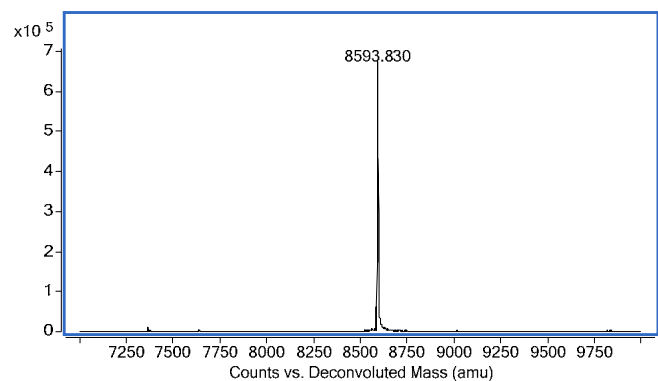
- Application:**
- Synthesis of polyubiquitin chains that lack the lysine 48-linkage for structural or DUB linkage specificity studies
  - Determination if polyubiquitin chain formation is through the lysine 48-linkage

### Product Information

<b>Purity:</b>	≥ 95% by RP-HPLC and SDS-PAGE
<b>Molecular Weight:</b>	8,592.9 Da
<b>Physical State:</b>	Liquid, PBS with 10% glycerol
<b>Quantity:</b>	1mg
<b>Solubility:</b>	>30mg/mL
<b>Storage:</b>	-80° C. Avoid repeated freeze/thaw cycles



**RP-HPLC**



**Deconvoluted MS**

### References

1. Wilkinson, K. D. (2000). "Ubiquitination and deubiquitination: targeting of proteins for degradation by the proteasome." *Semin Cell Dev Biol* 11(3): 141-8.
2. Ciechanover, A. (2003) "The ubiquitin proteolytic system and pathogenesis of human diseases: a novel platform for mechanism-based drug targeting." *Biochem Soc Trans* 31(2): 474-81.

All products are for research use only • not intended for human or animal diagnostic or therapeutic uses  
Copyright © 2014 LifeSensors, Inc. All Rights Reserved