

## Human recombinant ubiquitin K63

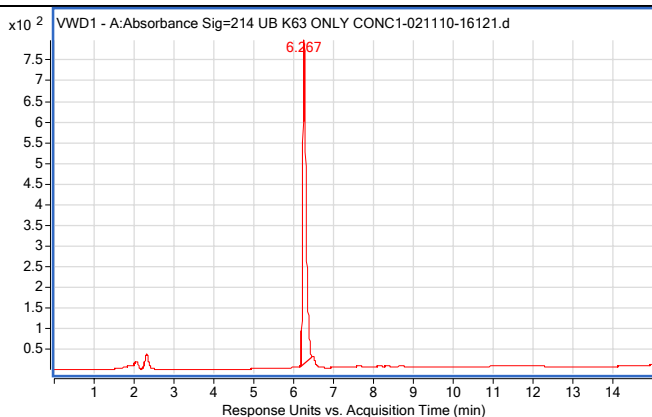
Cat. # SI208

**Background:** Human recombinant ubiquitin containing a single lysine at position 63 is produced in *E. coli* and purified to >95% homogeneity. This protein contains no extraneous tags.

**Application:** For use in conjunction with the E3LITE Customizable Ligase Assay and E2 Profiling Kit to investigate conjugation/ligation enzyme activity and/or drug discovery. Can also be used for conventional ubiquitin pathway research.

### Product Information

<b>Purity:</b>	≥ 95% by RP-HPLC
<b>Molecular Weight:</b>	8,732.9 Da
<b>Physical State:</b>	Liquid, 50mM Tris, pH 7.5, 0.15M NaCl, 10% glycerol
<b>Quantity:</b>	1 mg
<b>Solubility:</b>	>30 mg/ml
<b>Storage:</b>	-80° C. Avoid repeated freeze/thaw cycles



### RP-HPLC

### 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.
3. Marblestone, J.G., Kumar, K.G., Eddins, M.J., Leach C.A., Sterner, D.E., Mattern, M.R., Nicholson, B. (2010) Novel Approach for Characterizing Ubiquitin E3 Ligase Function. *J Biomol Screen* **15**:1220-1228.

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