

Human recombinant ubiquitin K27

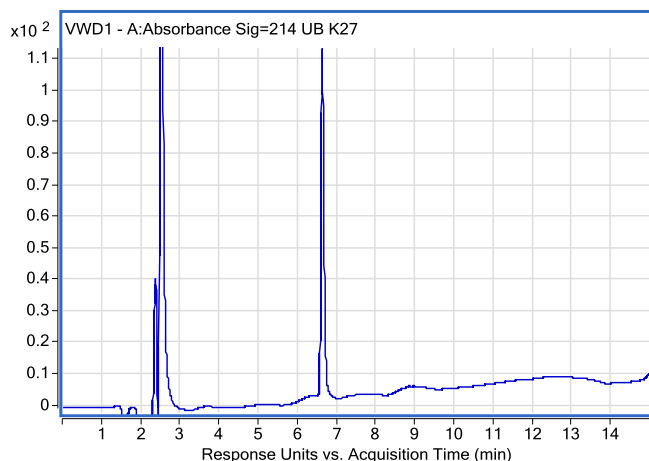
Cat. # SI204

Background: This tag-free recombinant form of human ubiquitin is engineered to have all available lysines mutated to arginines, except at position 27. This molecule, therefore, can only form polyubiquitin chains of K27 linkage type.

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

Purity:	≥ 95% by RP-HPLC
Molecular Weight:	8,732.9 Da
Physical State:	Liquid, PBS
Quantity:	1mg at 4 mg/mL
Solubility:	>30mg/ml
Storage:	-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* (In Press).

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