

Background

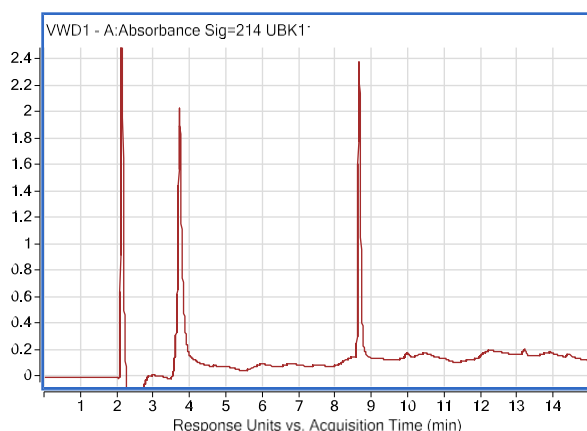
This tag-free recombinant form of human ubiquitin is engineered with all available lysines mutated to arginines. This molecule does not participate in polyubiquitin chain formation and can, therefore, be used as a control in studies of ubiquitination.

Application(s)

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

Product Specifications

Tag	None
Purity	≥ 95% by RP-HPLC & SDS-PAGE
Molecular Weight	8,760.9 Da
Quantity	1 mg
Species	Human
Expression System	<i>E. Coli</i>
Physical State	Liquid
Buffer	PBS
Solubility	>30 mg/ml
Storage	-80° C. Avoid repeated freeze/thaw cycles

Product QC**RP-HPLC****References**

1. Buneeva O, Medvedev A. Int J Mol Sci. 2022;23(7):3705.
2. Ohtake F., et al., Proc Natl Acad Sci U S A. 2018;115(7): E1401-E1408.
3. Marblestone, JG et al., J Biomol Screen. 2010; 15(10):1220-8.
4. Ciechanover, A. Biochem Soc Trans. 2003; 31(2): 474-81.
5. Wilkinson, K. D., Curr Opin Drug Discov Devel. 2000; 11(3): 141-8.

All products are for research use only • Not intended for human or animal diagnostic or therapeutic uses
Copyright © 2025 LifeSensors, Inc. All Rights Reserved