

Background

This tag-free recombinant form of human ubiquitin is engineered to have all available lysines mutated to arginines, except for the one at position 11. This molecule, therefore, can form only polyubiquitin chains of the K11 linkage type.

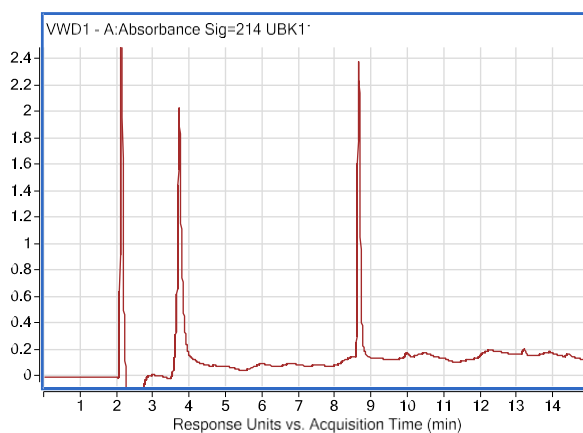
Application(s)

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
Molecular Weight	8,732.9 Da
Quantity	1 mg at 4 mg/mL
Species	Human
Expression System	<i>E. Coli</i>
Physical State	Liquid
Buffer	PBS
Solubility	>30 mg/ml
Stability & Storage	Over 1-year at -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) Wilkinson, K. D., Curr Opin Drug Discov Devel. 2000; 11(3): 141-8.
- 3) Ciechanover, A. Biochem Soc Trans. 2003; 31(2): 474-81.
- 4) Marblestone, JG et al., J Biomol Screen. 2010; 15(10):1220-8.

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