

**Background**

A mutant form of recombinant ubiquitin where lysine at position 6 has been replaced by arginine.

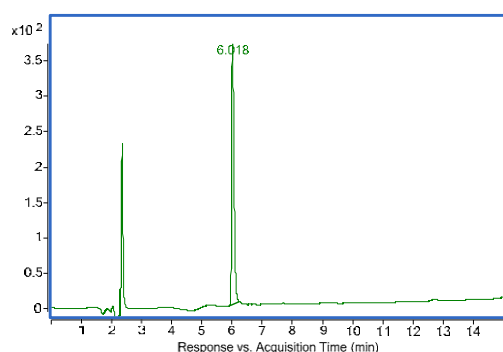
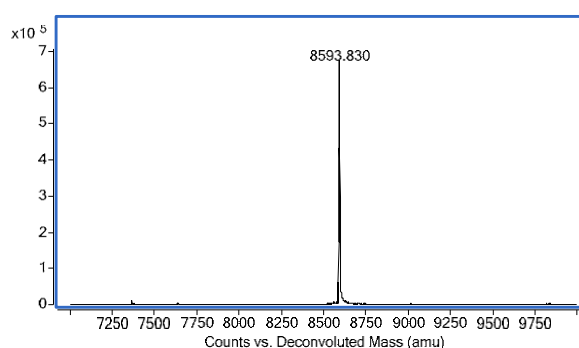
Ubiquitin is a small polypeptide that can be conjugated via its C-terminus to the amine group of a lysine residue on target proteins. This conjugation is referred to as monoubiquitination. Additional ubiquitin moieties can subsequently be conjugated to the initial ubiquitin using any one of the seven lysine residues on its surface. The formation of these ubiquitin chains is referred to as polyubiquitination.

**Application(s)**

- 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 Specifications**

|                     |  |
|---------------------|--|
| Tag                 | None   |
| Purity              | ≥ 95% by RP-HPLC AND SDS-PAGE                        |
| Molecular Weight    | 8,592.9 Da   |
| Quantity            | 1 mg at 4 mg/ml                                      |
| Species             | Human  |
| Expression System   | <i>E. Coli</i>                                       |
| Physical State      | Liquid   |
| Buffer              | 20 mM Tris, pH 7.5, 0.15 M NaCl, 10% glycerol        |
| Stability & Storage | Over 1 year -80°C. Avoid repeated freeze/thaw cycles |

**Product QC****RP-HPLC****Deconvoluted MS****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.

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