

Background

A mutant form of recombinant ubiquitin where lysine at position 27 has been replaced by arginine.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)

- Investigation of E1-E2-E3 mediated ubiquitin ligation.
- Investigation of DUB linkage specificity.

Product Specifications

Tag	None
Purity	≥ 95% by RP-HPLC
Molecular Weight	8,592 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

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.