

Ubiquitin-K63R

Cat. # SI218

Background: This tag-free recombinant form of human ubiquitin is engineered to have a single substitution of arginine for lysine at position 63. This mutant ubiquitin cannot form polyubiquitin chains through the lysine 63 linkage.

- Application:**
- 1) Synthesis of polyubiquitin chains that lack the lysine 63-linkage for structural or DUB linkage specificity studies
 - 2) Determination if polyubiquitin chain formation is through the lysine 63 linkage

Product Information

Purity:	≥ 95% by RP-HPLC and SDS-PAGE
Molecular Weight:	8,592 Da
Physical State:	Liquid, PBS
Quantity:	1mg
Solubility:	>30mg/ml
Storage:	-80°C. Avoid repeated freeze/thaw cycles

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.

All products are for research use only • not intended for human or animal diagnostic or therapeutic uses
Copyright © 2012 LifeSensors, Inc. All Rights Reserved