

Ubiquitin-K48R K63R

Cat. # SI219

Background: This tag-free recombinant form of human ubiquitin is engineered to have a single substitution of arginine for lysine at position 48 and 63.

Application:

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

Product Information

Source	Human recombinant
Form	liquid
Molecular weight	8,592 Da
Purity	>=95% by RP-HPLC
Affinity Tag	None
Concentration	4 mg/mL
Buffer	20 mM Tris, pH 7.5, 0.15 M NaCl, 10% glycerol
Storage	Long term storage at -80°C. Avoid multiple freeze/thaw cycles.
Stability	1 year under recommended storage conditions

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