

## Ubiquitin-K11R

Cat. # SI213

**Background:** This tag-free recombinant form of human ubiquitin is engineered to have a single substitution of arginine for lysine at position 11.

**Application:**

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

### Product Information

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