

## Ubiquitin-K6R

Cat. # SI212

---

**Background:** Recombinant ubiquitin having a single substitution of arginine for lysine at position 6.

Ubiquitin is a small polypeptide that can be conjugated via its C-terminus to amine groups of lysine residue on target proteins. This conjunction is referred to as monoubiquitylation. Additional ubiquitin moieties can be subsequently conjugated to this initial ubiquitin, utilizing any one (of seven) lysine residues on the surface of ubiquitin. The formation of these ubiquitin chains is referred to as polyubiquitylation.

**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

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