

## DUB-resistant, K11-linked Di-ubiquitin (Ub2)

Cat. # SI1112

- 
- Background:** DUB-resistant, K11-linked Di-ubiquitin (Ub2) contains an amino acid substitution near the C-terminus of the distal ubiquitin that renders it resistant to deubiquitylases (DUBs). DUB-resistant ubiquitin chains are ideal reagents for structural or binding studies with deubiquitylases or other ubiquitin binding proteins. They can also be used as inhibitors or to determine the linkage-specificity of DUBs. The product contains a wild-type C-terminus on the proximal ubiquitin.
- Application:** Structural or binding studies with deubiquitylases and other ubiquitin binding proteins. Inhibition or determination of linkage-specificity of DUBs
- 

### Product Information

<b>Purity:</b>	≥ 90% by RP-HPLC
<b>Molecular Weight:</b>	17,123.7 Da (calculated)
<b>Physical State:</b>	1mg/ml in 20 mM Tris pH 7.5, 0.15 M NaCl, 1 mM EDTA
<b>Quantity:</b>	100µg
<b>Storage:</b>	-80°C. Avoid repeated freeze/thaw cycles

---

All products are for research use only • not intended for human or animal diagnostic or therapeutic uses  
Copyright © 2013 LifeSensors, Inc. All Rights Reserved