

DUB-resistant, K63-linked Di-ubiquitin (Ub2)

Cat. # SI6312

-
- Background:** DUB-resistant, K63-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

Purity:	≥ 90% by RP-HPLC
Molecular Weight:	17,123.7 Da (calculated)
Physical State:	1mg/ml in 20 mM Tris pH 7.5, 0.15 M NaCl, 1 mM EDTA
Quantity:	100µg
Storage:	-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