

## DUB-resistant, K48-linked Di-ubiquitin (Ub2)

Cat. # SI4812

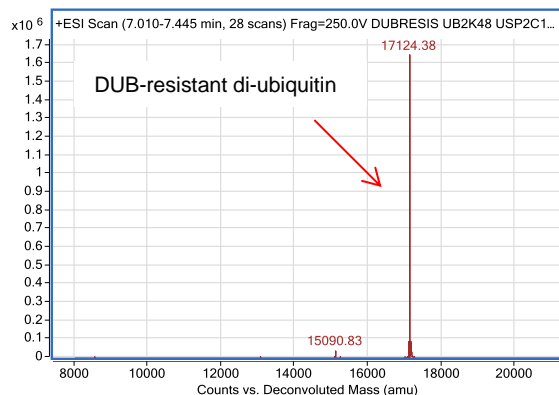
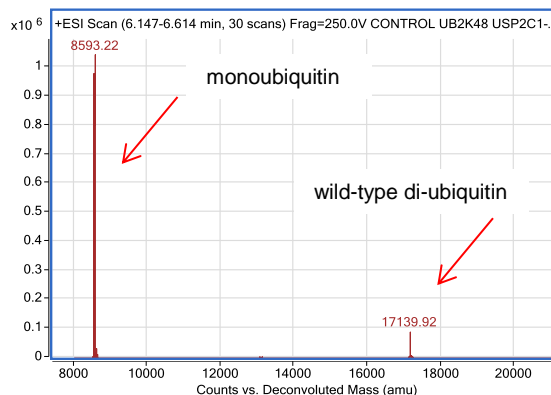
**Background:** DUB-resistant, K48-linked Di-ubiquitin (Ub2) contains an amino acid substitution near the C-terminus of the distal ubiquitin that renders it resistant to deubiquitylases (DUBs). DUB-resistant ubiquitin chains are ideal reagents for structural or binding studies with deubiquitylases or other ubiquitin binding proteins. They can also be used as inhibitors or to determine the linkage-specificity of DUBs. The product contains a K48R mutation in the distal ubiquitin and 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

### DUB-resistant, K48-linked Di-ubiquitin is not cleaved by USP2core

**DUB-resistant, K48-linked Di-ubiquitin (Ub2)****Wild-type di-ubiquitin (Ub2)**

**LC-MS analyses of DUB-resistant and wild-type di-ubiquitins treated with USP2core.** 25µg of DUB-resistant, K48-linked di-ubiquitin (Left panel; MW 17124.4) or 25µg of wild-type di-ubiquitin (Right panel; MW 17139.9) were treated with 0.5µM USP2core for 1 hour at room temperature and then subjected to LC-MS analysis. Neither loss of the di-ubiquitin peak or appearance of monoubiquitin were detected using the DUB-resistant di-ubiquitin whereas almost complete conversion of di-ubiquitin to monoubiquitin is observed with the wild-type di-ubiquitin.

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