

K63-linked di-ubiquitin (Ub2)

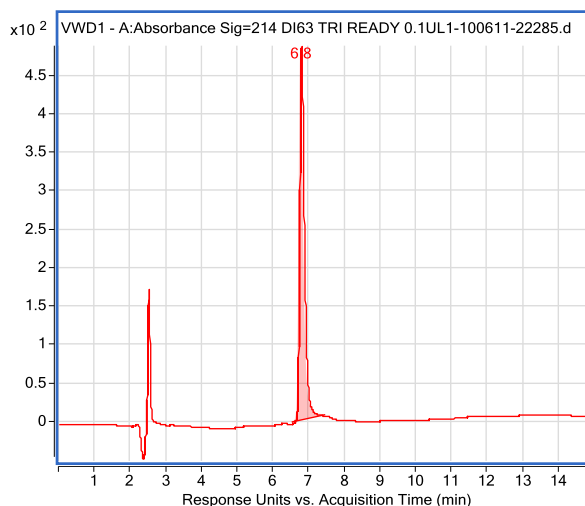
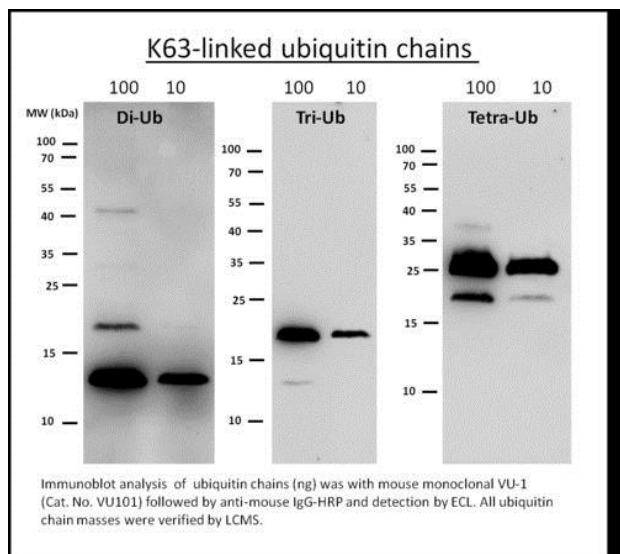
Cat. # SI6302

Background: Poly-ubiquitylation of target proteins through K63 has recently become the focus of intense study. Strikingly, the topology of this linkage and its apparent role in cellular processes are quite different from that of K48. Poly-Ub chains of this type appear to play a role in endocytic trafficking, DNA repair, neurodegeneration and more.

These di-ubiquitin chains are generated from the enzymatic linkage of wild-type ubiquitin through lysine 63. The most distal ubiquitin contains an arginine substitution for a lysine at position 63, limiting chain length.

Product Information

- Purity:** $\geq 95\%$ by Western blot analysis
- Molecular Weight:** 17,139.7 Da
- Physical State:** Liquid; 20mM Tris-HCl, pH 7.5, 0.15M NaCl, 1mM EDTA
- Quantity:** 50 μ g
- Solubility:** >1mg/mL
- Storage:** -80° C. Avoid repeated freeze/thaw cycles
- Concentration:** Lot dependent, please see Certificate of Analysis or vial label for actual concentration



RP-HPLC Profile

All products are for research use only • not intended for human or animal diagnostic or therapeutic uses
 Copyright © 2009 LifeSensors, Inc. All Rights Reserved