

K48-linked di-ubiquitin (Ub2)

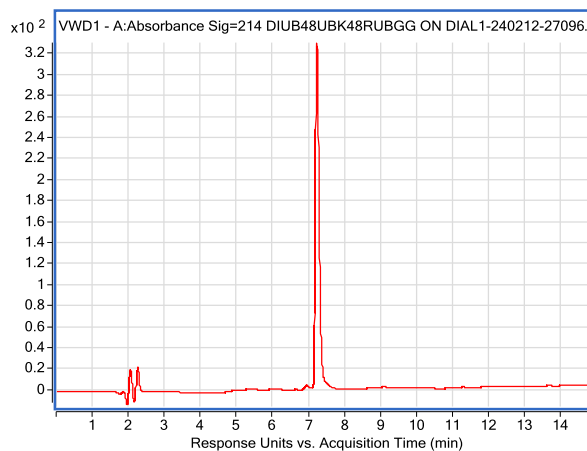
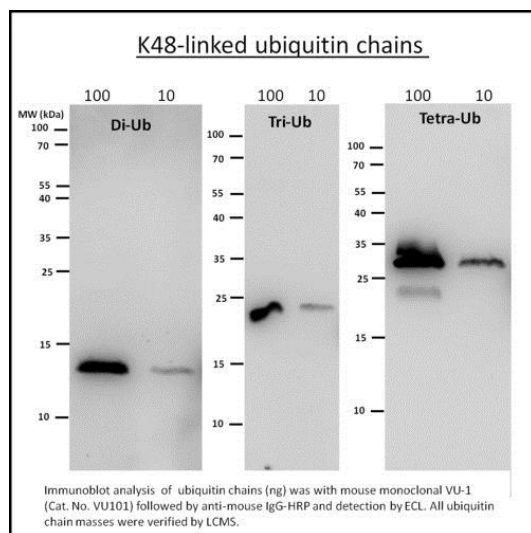
Cat. # SI4802

Background: Poly-ubiquitylation of target proteins through lysine 48 is the most thoroughly studied of the various chain linkages and once considered the hallmark of ubiquitylation. It is now clear that many, if not all, poly-Ub chain topologies play distinct and important roles in regulating cellular processes. Nevertheless, polyubiquitylation through lysine 48 is considered the primary signal for proteasome-mediated degradation.

K48-linked di-ubiquitin chains are generated from the enzymatic linkage (E2-25K) of wild-type ubiquitin through lysine 48. The most distal ubiquitin contains an arginine substitution for lysine at position 48, limiting chain length.

Product Information

Purity:	≥ 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:	100 µg
Solubility:	>1mg/mL
Storage:	-80° C. Avoid repeated freeze/thaw cycles
Concentration:	Lot dependent, please refer to Certificate of Analysis or vial label for actual concentration



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