

K48-linked tetra ubiquitin (Ub4)

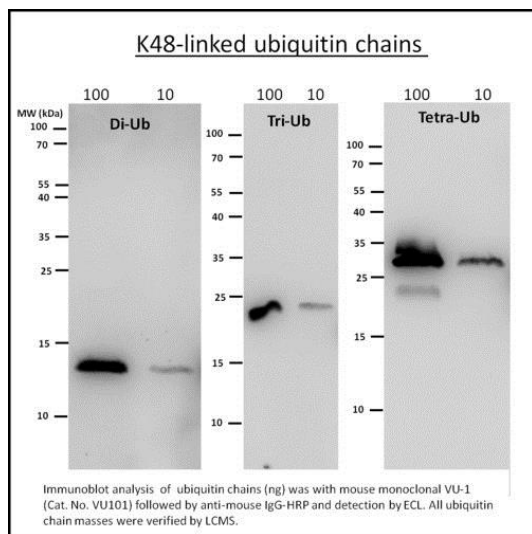
Cat. # SI4804

Background: Poly-ubiquitylation of target proteins through linkage at K48, is now the most thoroughly studied of the various chain linkages, and we once considered the hallmark of this post-translational modification. It is now clear that many, if not all, poly-Ub chain topologies likely play distinct and important roles in regulating cellular processes. Nevertheless, K48 linkage remains a critical pathway for the cells to maintain homeostasis through proteolytic degradation, and as such remains very intriguing for the study of DUBs that play a role in the degradation, as well as the proteasome itself.

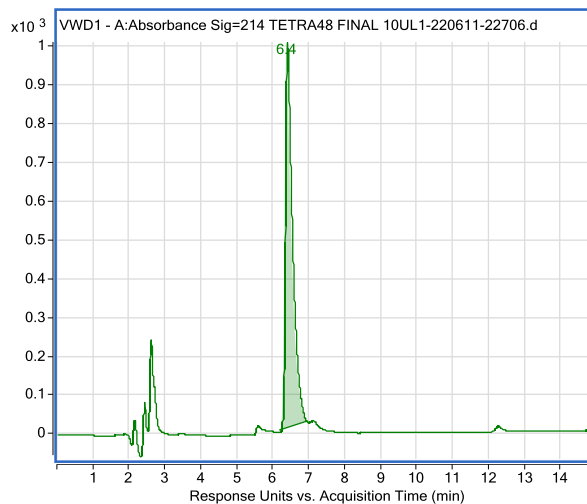
These tetra-ubiquitin chains generated from the enzymatic linkage (E2-25K) of wild-type ubiquitin through lysine 48. The most distal ubiquitin contains an arginine substitution for a lysine at position 48, limiting chain length.

Product Information

Purity:	≥ 95% by Western blot analysis
Molecular Weight:	34,233.4 Da
Physical State:	Liquid; 20mM Tris-HCl, pH 7.5, 0.15M NaCl, 1mM EDTA
Quantity:	25 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



Western blot analysis



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