

## TAMRA-K63-linked tetra-ubiquitin (K63-Ub4-TMR)

Cat. # SI6304T

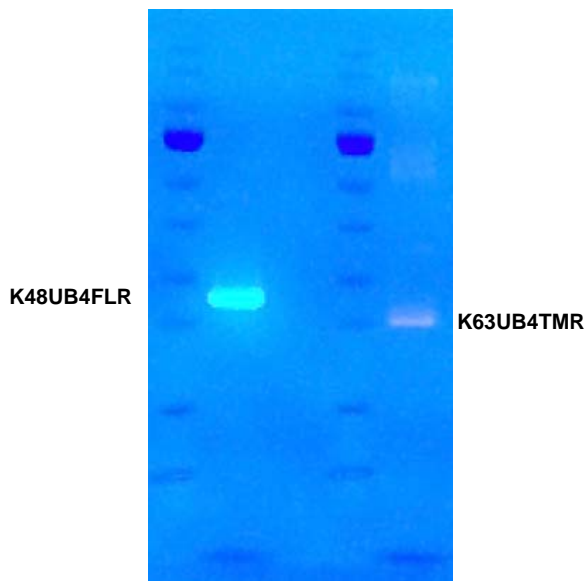
**Background:** Poly-ubiquitylation of target proteins through K63 has recently become the focus of intense study. Strikingly, the topology of this linkage and its apparent role in cellular processes are quite different from that of K48. Poly-Ub chains of this type appear to play a role in endocytic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

These tetra-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. The TAMRA moiety is attached to a single, non-lysine site in the proximal ubiquitin.

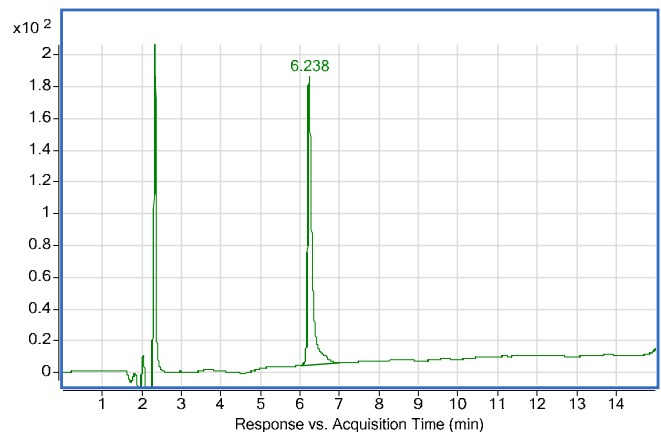
**Applications:** Optimization of pull-down conditions for anti-K48 TUBE and anti K63-TUBE reagents

### Product Information

<b>Purity:</b>	≥ 95% by RP-HPLC
<b>Molecular Weight:</b>	34,802.0 Da
<b>Physical State:</b>	Liquid; 20mM Tris-HCl, pH 7.5, 0.15M NaCl, 10% glycerol
<b>Quantity:</b>	25 µg
<b>Solubility:</b>	>1mg/mL
<b>Storage:</b>	-80° C. Avoid repeated freeze/thaw cycles
<b>Concentration:</b>	Lot dependent, please see Certificate of Analysis or vial label for actual concentration



**SDS-Gel, UV transillumination**



**RP-HPLC Profile**

All products are for research use only • not intended for human or animal diagnostic or therapeutic uses  
 Copyright © 2015 LifeSensors, Inc. All Rights Reserved

All products are for research use only • not intended for human or animal diagnostic or therapeutic uses  
Copyright © 2015 LifeSensors, Inc. All Rights Reserved