

## Fluoresceinated-K48-linked tetra ubiquitin (K48-Ub4-FLR)

Cat. # SI4804F

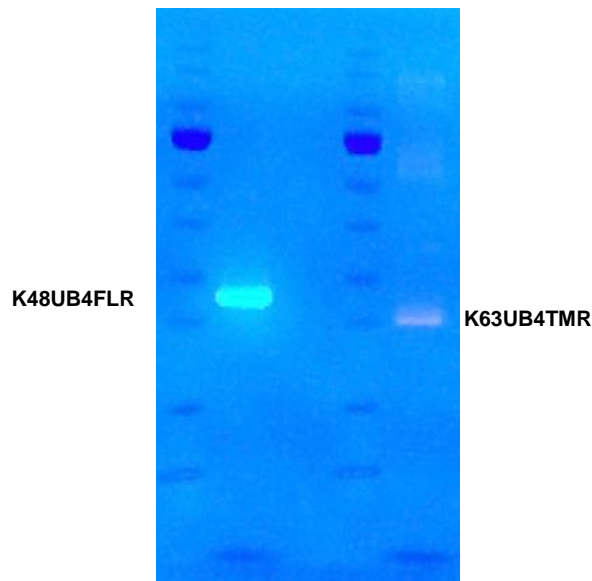
**Background:** Poly-ubiquitylation of target proteins through linkage at K48, is now the most thoroughly studied of the various chain linkages, and was once considered the hallmark of this post-translational modification. It is now clear that many, if not all, poly-Ub chain topologies likely play distinct 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 are generated by the enzymatic linkage of wild-type ubiquitin through lysine 48. The most distal ubiquitin contains a Lys48 to Arg substitution limiting chain length. The fluorescein moiety is attached to a single, non-lysine site in the proximal ubiquitin.

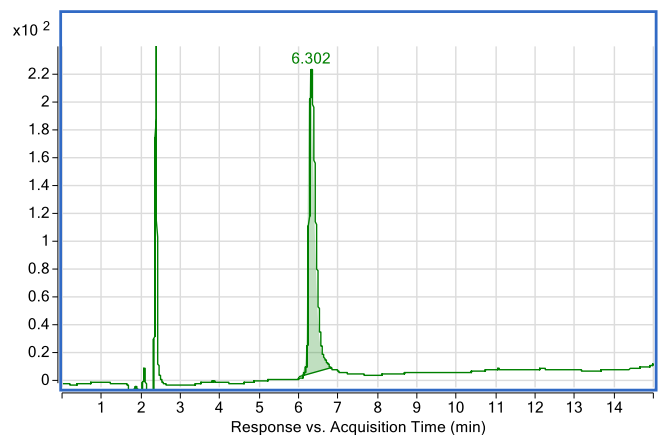
**Applications:** Optimization of pull-down conditions for anti-K48 TUBE and anti K63-TUBE reagents

### Product Information

<b>Purity:</b>	≥ 95% RP-HPLC
<b>Molecular Weight:</b>	34,678.3 Da
<b>Physical State:</b>	Liquid
<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 refer to 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 © 2017 LifeSensors, Inc. All Rights Reserved