

Fluorescein-labeled K63 TUBE1 (FLR-K63 TUBE1)

Cat. # UM504F

Background: Fluorescein-labeled (FLR) K63 TUBE 1 binds K63-linked polyubiquitin with a 1000 to 10,000-fold preference for K63 over other linkage types and provides a sensitive and cost-effective tool for determining the abundance of K63-linked polyubiquitin in the cell, or identifying the linkage-type associated with a particular protein of interest. Based on an engineered peptide developed by [Sims et al.](#) [1], FLR-K63 TUBE 1 consists of multiple ubiquitin interacting motifs separated by rigid linkers, presumably increasing its affinity to extended K63 polyUb chains. FLR- K63 TUBE 1 is useful for detection and characterization of K63-linked polyubiquitylated proteins by 'far Western' blotting and *in situ* detection (histo- or cyto-chemical staining).

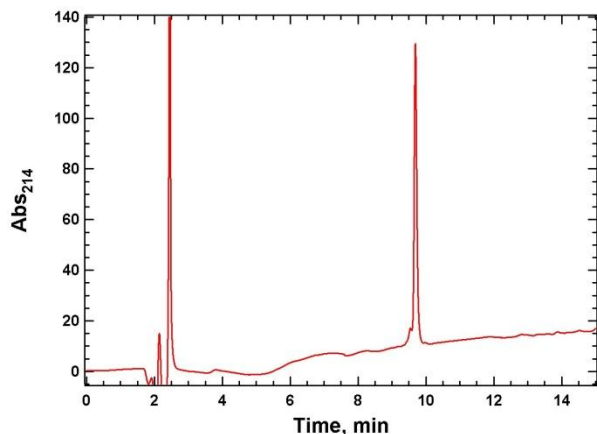
FLR-K63 TUBE 1 has one fluorescein molecule conjugated to a single, defined location within the protein.

Application:

- "Far Western" blot analysis
- Histo- or cyto-chemical staining. *N.b.* Sample fixation will require optimization.

Product Information

Purity:	≥ 95% by RP-HPLC
Molecular Weight:	10,222 Da
Physical State:	Liquid
Quantity:	50 µg
Buffer:	PBS
Storage:	-80° C. Avoid repeated freeze/thaw cycles



RP-HPLC

References

1. Sims, J.J., Scavone, F., Cooper, E.M., Kane, L.A., Youle, R.J., Boeke, J.D., and Cohen, R.E., Polyubiquitin-sensor proteins reveal localization and linkage-type dependence of cellular ubiquitin signaling. *Nat Methods*. **9**:303-309 (2012).

All products are for research use only • not intended for human or animal diagnostic or therapeutic uses

Copyright © 2009 LifeSensors, Inc. All Rights Reserved