

Anti-M1 TUBE, FLAG

Cat. # UM606

Background: LifeSensors has developed TUBE (Tandem Ubiquitin Binding Entity) technology for the detection, characterization and isolation of polyubiquitylated proteins from cells and tissue extracts. TUBEs capitalize on the linkage of multiple ubiquitin interacting motifs (UIMs) to generate reagents with high affinity for polyubiquitin. TUBEs have 100 to 1000-fold higher affinity for polyubiquitin chains compared to monomer ubiquitin binding domains (UBDs). TUBEs both stabilize and bind to ubiquitylated proteins, serving as an indispensable tool for ubiquitologists. Our anti-M1 (linear) TUBE has high selectivity for M1-linked polyubiquitin over the more common K48- and K63-linked polyubiquitin chains, making it a powerful tool for studying the biological consequences of this ubiquitin linkage type.

Application:

- Isolation and enrichment of M1-polyubiquitinated proteins from cell and tissue extracts
- Isolation of ubiquitylated proteins for proteomic studies

Product Information

Affinity tag:	FLAG
Purity:	≥ 90% by SDS-PAGE
Molecular Weight:	23.1 kDa
Physical State:	Liquid
Quantity:	200 µg
Storage:	-80° C. Avoid repeated freeze/thaw cycles

References

1. Hjerpe, R., F. Aillet, F. Lopitz-Otsoa, V. Lang, P. England, and M.S. Rodrigues, *Efficient protection and isolation of ubiquitylated proteins using tandem ubiquitin-binding entities*. EMBO Rep, 2009. **10**: p. 1250-1258.
2. Aillet, F., F. Lopitz-Otsoa, R. Hjerpe, M. Torres-Ramos, V. Lang, and M.S. Rodriguez, *Isolation of ubiquitylated proteins using tandem ubiquitin-binding entities*. Meth Mol Biol, 2012. **832**: p. 173-183.

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