

## Background

Based on protein domains known to possess an affinity for ubiquitin, Tandem Ubiquitin Binding Entities (TUBEs) have been developed for the isolation and identification of ubiquitinated proteins. TUBEs display up to a 1000-fold increase in affinity for poly-ubiquitin moieties over the single ubiquitin-binding associated domain (UBA). In addition, TUBEs display a protective effect on polyubiquitinated proteins, allowing for detection at relatively low abundance. These properties effectively "capture" proteins in their polyubiquitin state. [UM502M](#) was designed by coating high-capacity magnetic beads to allow superior enrichment of poly-ubiquitinated proteins along with minimizing non-specific binding to proteins in tissue and cellular lysates.

The affinity of TUBE 2 for K63 linked tetra-ubiquitin is approximately equal to K48 linked tetra-ubiquitin (5-10nM).

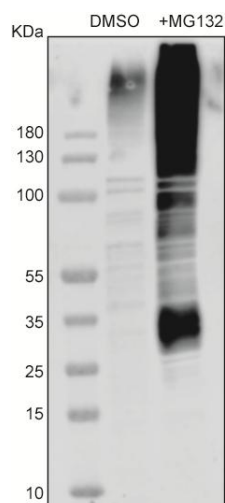
## Application(s)

- Pull down of poly-ubiquitinated proteins from cell lines, tissues, and organs
- Protection of polyubiquitinated proteins from both deubiquitination and proteasomal degradation
- [Bead-based ELISA](#) for studying ubiquitinated proteome
- Label-free [mass spectrometry for ubiquitinated proteomic analysis](#)

## Product Specifications

Affinity Tag	None
Purity	(prior to coupling) > 95% estimated by SDS-PAGE
Quantity	1 ml
Expression System	<i>E. coli</i>
Physical State	Liquid
Storage	+4°C. Avoid storage at lower temperatures

## Product QC



**Enrichment of total ubiquitinated proteome using UM502M.** 100  $\mu$ L of UM502M beads were added to 300  $\mu$ g of cell lysates derived from HeLa cells treated with either DMSO or 1  $\mu$ M MG-132 for 4 hours. The data represent overnight enrichment of both DMSO- and MG-132-treated lysates using UM502M at 4 °C on an end-to-end rotator. A characteristic increase in high-molecular-weight ubiquitin smears was observed in the MG-132-treated samples, indicating robust pull-down of the polyubiquitinated proteome. The enriched beads were resuspended in 30  $\mu$ L of 1X Laemmli sample buffer and loaded onto a 10% SDS-PAGE gel. The western blot membrane was probed using [anti-ubiquitin \(VU101\)](#).

## References

1. Garadi Suresh H et al., Mol Cell, 2024;84(12):2337-2352.
2. Kadimisetty K., et al., Methods Mol Biol, 2021;2365:185-202.
3. Hjerpe, R, et al., EMBO Rep., 2009; 10,1250-1258.

All products are for research use only • Not intended for human or animal diagnostic or therapeutic uses  
Copyright © 2025 LifeSensors, Inc. All Rights Reserved