

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.

UM501M was designed by coating polymeric high-capacity magnetic beads to allow superior enrichment of polyubiquitinated proteins while minimizing non-specific binding to proteins in tissue and cellular lysates.

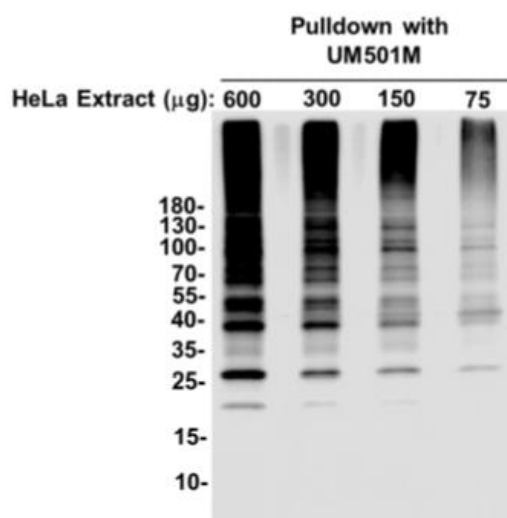
Application(s)

- Pull down of poly-ubiquitinated proteins from cell lines, tissues and organs
- Protection of poly-ubiquitinated proteins from both deubiquitination and degradation by the proteasome

Product Specifications

Affinity Tag	None
Purity	(prior to coupling) $\geq 95\%$ by SDS-PAGE
Quantity	1ml
Expression System	<i>E. coli</i>
Physical State	Liquid
Buffer	PBS, pH 7.2, 0.05% sodium azide
Stability	Over 1 year at +4 °C. Do not centrifuge, dry or freeze the beads.

Product QC



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.