

Magnetic Beads-M1 TUBE

Cat. # UM406M

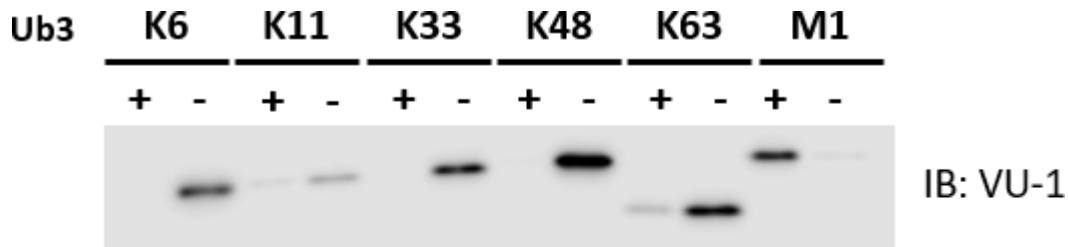
Background: Based on protein domains known to possess an affinity for ubiquitin, Tandem Ubiquitin Binding Entities (TUBEs) have been developed for the isolation and identification of ubiquitylated proteins. TUBEs display up to a 1000-fold increase in affinity for poly-ubiquitin moieties over the single ubiquitin binding associated domain (UBA). In addition, TUBEs display a protective effect on polyubiquitylated proteins, allowing for detection at relatively low abundance. These properties effectively "capture" protein in its polyubiquitin state.

Magnetic M1 TUBE is based on UBA from the protein ubiquilin. The affinity of M1 TUBE for M1 linked tetra-ubiquitin is approximately 10-fold higher than for K48 linked chains.

- Application:**
- Pull down of poly-ubiquitylated proteins from cell lines, tissues, and organs
 - Protection of poly-ubiquitylated proteins from both deubiquitylation and degradation by the proteasome

Product Information

Affinity tag:	N/A
Purity:	(prior to coupling) > 95% by SDS-PAGE
Molecular Weight:	not applicable
Physical State:	Liquid
Quantity:	1mL magnetic beads
Solubility:	not applicable
Storage:	+4 °C. Avoid storage at lower temperatures.



** + refers to the bound fraction

** - refers to the unbound fraction

References

Hjerpe, R, Aillet, F, Lopitz-Otsoa, F, Lang, V, England, P, and Rodriguez, MS., [Efficient protection and isolation of ubiquitylated proteins using tandem ubiquitin-binding entities](#). *EMBO Rep.* **10**,1250-1258 (2009).

All products are for research use only • not intended for human or animal diagnostic or therapeutic uses
 Copyright © 2009 LifeSensors, Inc. All Rights Reserved