

TUBE 1 (Magnetic Beads)

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 polyubiquitin moieties over the single ubiquitin binding associated domain (UBA). In addition, TUBEs protect polyubiquitinated proteins from deubiquitination and proteasomal degradation, allowing for detection at relatively low abundance. These properties effectively allow TUBEs to “capture” protein in its polyubiquitinated state. TUBEs can be used to isolate, enrich, and identify ubiquitinated proteins from cell and tissue extracts. TUBE1 is based on UBAs from the protein ubiquilin. Magnetic beads coupled TUBE1 allow for efficient recovery of polyubiquitinated proteins in a single step without centrifuge. The magnetic TUBEs allows more complete removal of the unbound supernatant, which leads to lower background. Polyubiquitinated protein can easily be separated from TUBEs for proteomic studies or identification by Western blotting.

Ordering Information

[Order Online »](#)

LSI-UM-0401M-1000	1ml
-------------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Notes Protein isolation



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com