

# UBIQAPTURE-Q<sup>®</sup> kit

**Market-leading tool for isolation and detection of ubiquitinated proteins.**

For isolation and enrichment of ubiquitinated proteins. The UBI-QAPTURE-Q<sup>®</sup> kit is an efficient tool for the selective isolation of ubiquitinated proteins. The Kit facilitates the isolation of both mono- and poly-ubiquitinated proteins (independent of lysine residue chain linkage) from cell extracts, tissue lysates and *in vitro* assay solutions through use of a high-binding affinity matrix. Captured proteins can be analyzed by Western blotting using the highly sensitive ubiquitin-conjugate specific antibody provided, using antibodies to specific proteins of interest, or eluted from the matrix for subsequent biochemical characterization. The UBI-QAPTURE-Q<sup>®</sup> matrix supplied with the kit has superior binding characteristics compared to other commercially available matrices and is compatible with a wide range of lysate buffers and cell/tissue samples from a variety of species.

Citations: 46

[View Online »](#)

## Ordering Information

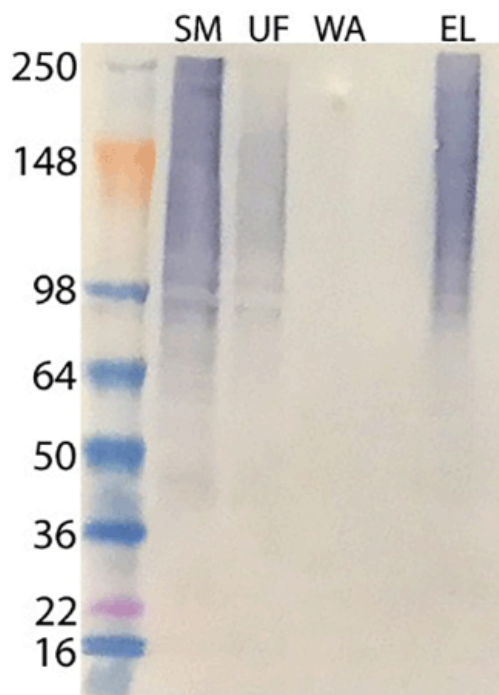
[Order Online »](#)

BML-UW8995A-0001	20 tests
------------------	----------

## Manuals, SDS & CofA

[View Online »](#)

- Efficient tool for the selective isolation of ubiquitinated proteins
- High binding affinity matrix, able to bind both monoubiquitinated and polyubiquitinated proteins
- Highly compatible with a wide range of lysate buffer, different sample types (cells, tissues) and a broad range of species



## Handling & Storage

<b>Use/Stability</b>	Upon receipt, the control lysate (Prod. No. BML-UW0130) should be stored at -80°C. All other components should be stored at 4°C.
<b>Long Term Storage</b>	-80°C
<b>Shipping</b>	Dry Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

### Application Notes

#### Uses:

1. Isolation and detection of the full range of ubiquitinated protein conjugates (mono-/multi-/poly-ubiquitin modified, lysine linkage independent) from specific cell/tissue lysates of interest, especially low abundance (regulatory) proteins involved in the ubiquitin-proteasome pathway.
2. Capture and analysis of specific ubiquitinated protein conjugates of interest from particular cell/tissue lysates.
3. Purification/pull down of ubiquitinated proteins from cell free in vitro assays; for example to isolate Ub-p53 species from ubiquitination assays mediated by HeLa S100 lysate (Prod. No. BML-SW8750).
4. Release of free proteins in their active/native form by cleavage of ubiquitin/ubiquitin chains from the UBI-QAPTURE-Q<sup>®</sup> matrix using deubiquitinating enzymes such as USP2 (Prod. No. BML-UW9850).
5. Release of ubiquitinated proteins in their active/native form by elution from the UBI-QAPTURE-Q<sup>®</sup> matrix using, for example, high salt buffer.

**Note:** Protocol provided covers applications 1-3. Assay set-up can be readily modified for alternative applications by inclusion, omission or substitution of specific components. Kit is not suitable for use with purified proteins.

### Contents

**UBI-QAPTURE-Q<sup>®</sup> matrix** (Prod. No. BML-UW0125) 0.5 ml settled resin provided; 50% suspension.

**Control ubiquitinated-protein lysate (Ub-lysate)** (Prod. No. BML-UW0130) 500 µg (100µl), 5mg/ml in 50mM HEPES, pH 7.6, 1mM DTT.

**Mono- and polyubiquitinated conjugates monoclonal antibody (UBCJ2) (HRP conjugate)** (Ubiquitin-conjugates specific HRP-linked antibody solution) (Prod. No. ENZ-ABS840HRP) 25µl.

### Quantity

Provides sufficient material for 20 x 20µl reactions.



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

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

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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