UBIQAPTURE-Q® kit

Market-leading tool for isolation and detection of ubiquitinylated proteins.

For isolation and enrichment of ubiquitinylated proteins. The UBI-QAPTURE-Q[®] kit is an efficient tool for the selective isolation of ubiquitinylated proteins. The Kit facilitates the isolation of both mono- and poly-ubiquitinylated 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[®] 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 »

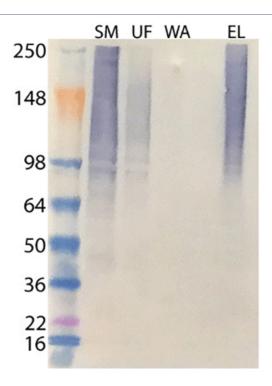
BML-UW8995A-0001

20 tests

Manuals, SDS & CofA

View Online »

- Efficient tool for the selective isolation of ubiquitinylated 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



Western Blot analysis of ubiquitin enrichment of lysate derived ubiquitinylated proteins. Ubiquitin enrichment experiment set-up and run as described in UBIQAPTURE-Q® kit (Prod. No. BML-UW8995A). Ubiquitin-protein conjugates present in starting material, unbound fraction and elution fraction were detected by western blotting using the provided ubiquitin-conjugate specific HRP-linked antibody (Prod. No. ENZ-ABS840HRP) at a dilution of 1:1000. Key: SM = Starting Material, UF = Unbound Fraction, WA = Wash Fraction, and EL = Elution Fraction.

Handling & Storage

Use/Stability Upon receipt, the control lysate (Prod. No. BML-UW0130) should be stored at -80°C. All

other components should be stored at 4°C.

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Notes

Uses:

- Isolation and detection of the full range of ubiquitinylated 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 ubiquitinylated protein conjugates of interest from particular cell/tissue lysates.
- Purification/pull down of ubiquitinylated proteins from cell free in vitro assays; for example to isolate Ub-p53 species from ubiquitinylation assays mediated by HeLa S100 lysate (Prod. No. BML-SW8750).
- Release of free proteins in their active/native form by cleavage of ubiquitin/ubiquitin chains from the UBI-QAPTURE-Q[®] matrix using deubiquitinylating enzymes such as USP2 (Prod. No. BML-UW9850).
- Release of ubiquitinylated proteins in their active/native form by elution from the UBI-QAPTURE-Q[®] 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

 $\label{eq:bound} \begin{array}{l} \textbf{UBI-QAPTURE-Q}^{\circledR} \ \textbf{matrix} \ (\text{Prod. No. BML-UW0125}) \ 0.5 \\ \text{ml settled resin provided; 50\% suspension.} \end{array}$

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

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

Provides sufficient material for 20 x 20µl reactions.

Quantity

Last modified: May 29, 2024



ENZO LIFE SCIENCES,

usa@enzolifesciences.com

Phone: 800.942.0430

info-