

## **Auto-ubiquitylation Kit**

Catalog #BML-UW0970-0001

For assessment of protein Ub E3 ligase activity

\* New antibody included.

## **USE FOR RESEARCH PURPOSES ONLY**

Unless otherwise specified expressly on the packaging, all products sold hereunder are intended for and may be used for research purposes only and may not be used for food, drug, cosmetic or household use or for the diagnosis or treatment of human beings. Purchase does not include any right or license to use, develop or otherwise exploit these products commercially. Any commercial use, development or exploitation of these products or development using these products without the express written authorization of Enzo Life Sciences, Inc. is strictly prohibited. Buyer assumes all risk and liability for the use and/or results obtained by the use of the products covered by this invoice whether used singularly or in combination with other products.

## **LIMITED WARRANTY; DISCLAIMER OF WARRANTIES**

These products are offered under a limited warranty. The products are guaranteed to meet all appropriate specifications described in the package insert at the time of shipment. Enzo Life Sciences' sole obligation is to replace the product to the extent of the purchasing price. All claims must be made to Enzo Life Sciences, Inc., within five (5) days of receipt of order. THIS WARRANTY IS EXPRESSLY IN LIEU OF ANY OTHER WARRANTIES OR LIABILITIES, EXPRESS OR IMPLIED, INCLUDING WARRANTIES OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, AND NONINFRINGEMENT OF THE PATENT OR OTHER INTELLECTUAL PROPERTY RIGHTS OF OTHERS, AND ALL SUCH WARRANTIES (AND ANY OTHER WARRANTIES IMPLIED BY LAW) ARE EXPRESSLY DISCLAIMED.

## **TRADEMARKS AND PATENTS**

Several Enzo Life Sciences products and product applications are covered by US and foreign patents and patents pending.

**FOR RESEARCH USE ONLY.  
NOT FOR USE IN DIAGNOSTIC PROCEDURES.**



Please read entire booklet before proceeding with the assay.



Carefully note the handling and storage conditions of each kit component.



Please contact Enzo Life Sciences Technical Support if necessary.

## Table of Contents

Background .....	2
Kit Description .....	3
Suggested Application.....	3
Kit Components.....	4
Storage .....	5
Other Materials Needed .....	5
Auto-Ubiquitylation Assay.....	5
Analysis by Western Blotting .....	7
Example Procedure for Western Blotting.....	8
Example Results for Western Blotting .....	9
References.....	11
Contact Information.....	10

## BACKGROUND

The covalent attachment of ubiquitin to proteins (ubiquitylation) plays a fundamental role in the regulation of cellular function through biological events involving cell cycle, differentiation, immune responses, DNA repair, chromatin structure, and apoptosis<sup>1,2,3,4</sup>.

Ubiquitylation is achieved through three enzymatic steps. In an ATP-dependent process, the ubiquitin activating enzyme (E1) catalyzes the formation of a reactive thioester bond with ubiquitin, followed by its subsequent transfer to the active site cysteine of a ubiquitin carrier protein (E2). The selectivity of the ubiquitin cascade for a particular substrate protein relies on the interaction between the E2 conjugating enzyme (of which a cell contains relatively few) and a ubiquitin-protein ligase (E3), of which over 600 have been identified to date<sup>2,5</sup>.

The E3s are a large, diverse group of proteins, characterized by one of several defining motifs. These include a HECT (homologous to E6-associated protein C-terminus), RING (really interesting new gene) or U-box (a modified RING motif without the full complement of Zn<sup>2+</sup>-binding ligands) domain. Whereas HECT E3s have a direct role in catalysis during ubiquitylation, RING and U-box E3s facilitate protein ubiquitylation. These latter two E3 types act as adaptor-like molecules. They bring an E2 and a substrate into sufficiently close proximity to promote the substrate's ubiquitylation. Although many RING-type E3s, such as MDM2 and c-Cbl, can apparently act alone, others are found as components of much larger multi-protein complexes, such as the anaphase-promoting complex.

Taken together, these multifaceted properties and interactions enable E3s to provide a powerful, and specific, mechanism for protein clearance within all cells of eukaryotic organisms utilizing the ubiquitin-proteasome system. The importance of E3s is highlighted by the number of normal cellular processes they regulate, and the number of diseases associated with their loss of function or inappropriate targeting.

E3 ligases also undergo auto-ubiquitylation, through modification of specific lysine residues within an individual ligase, providing a mechanism thought to be responsible for the regulation of the E3 enzyme itself<sup>5,6</sup>.

## KIT DESCRIPTION

This E3 ligase auto-ubiquitylation kit enables proteins to be tested for ubiquitin E3 ligase activity through assessment of their ability to undergo auto-ubiquitylation. Utilizing the first three steps in the ubiquitin cascade the kit facilitates ubiquitylation of known or putative E3 ligase enzymes followed by Western blot analysis using the highly sensitive reagents provided or using antibodies to the specific protein of interest (user supplied). A high integrity ubiquitin E3 ligase enzyme is also provided for use as a positive control.

The Kit provides sufficient material for approximately 10 auto-ubiquitylation assays.

## SUGGESTED APPLICATION

1. Qualitative assessment of an Ub E3 ligase enzyme's activity through its ability to auto-ubiquitylate.
2. Testing of proteins for auto-ubiquitylation activity allowing their identification as putative ubiquitin E3 ligases.
3. Ubiquitylation of substrate proteins (user provided) specific to a particular ubiquitin E3 ligase.

**Note:** Protocol provided covers applications 1-2. Assay set-up can be readily modified for alternative applications by inclusion, omission or substitution of specific components.

## KIT COMPONENTS

**1. Ubiquitin Activating Enzyme Solution, E1 (20X):**

*20X Ub Activating E1*

Ubiquitin activating enzyme E1 (human), (recombinant) (His-tag) (BML-UW9410-0025).

Use 2.5µL per 50µL reaction.

25µL provided, sufficient for 10 x 50µL reactions.

**2. UbcH5a (h)(rec.) (His-tag) (20X):**

*20X Ub Conjugating E2*

UbcH5a (human), (recombinant) (His-tag) (BML-UW9050-0025).

Use 2.5µL per 50µL reaction.

25µL provided, sufficient for 10 x 50µL reactions.

**3. Ubiquitin (10X):**

*10X Ubiquitin Solution*

Ubiquitin (BML-UW8795-0050).

Use 5µL per 50µL reaction.

50µL provided, sufficient for 10 x 50µL reactions.

**4. Hdm2 (h) (rec.) (catalytic RING domain) (GST-tag) (20X):**

*20X Control Ub Ligase E3*

Hdm2 (catalytic RING domain) (human), (recombinant) (GST-tag) (BML-UW0200-9090).

Use 2.5µL per 50µL reaction.

25µL provided, sufficient for 10 x 50µL reactions.

**5. Mg<sup>2+</sup>-ATP Solution (20X):**

*20X Mg<sup>2+</sup>-ATP Solution*

Mg<sup>2+</sup>/ATP (BML-EW9805-0025).

Use 2.5µL per 50µL reaction.

25µL provided, sufficient for 10 x 50µL reactions.

**Note:** Ensure Mg<sup>2+</sup>-ATP is fully dissolved by warming to room temperature and mixing by vortex prior to use.

**6. Ub E3 Ligase Buffer (10X):**

*10X Ub E3 Ligase Buffer (BML-KW0965-0050)*

Use 5µL per 50µL reaction

50µL provided, sufficient for 10 x 50µL reactions

**7. Ubiquitin mAb (P4G7-H11):**

*Ubiquitin Antibody Solution*

Ubiquitin, mouse monoclonal antibody (ADI-SPA-203-0025)

25µL provided. Dilution of at least 1:500-1:1000 recommended for Western blotting.

## STORAGE

All kit components should be stored at  $-80^{\circ}\text{C}$  to ensure stability and activity. Components with storage temperatures other than  $-80^{\circ}\text{C}$  can be stored at the temperature listed on the label OR at  $-80^{\circ}\text{C}$ . Avoid multiple freeze/thawing.

## OTHER MATERIALS NEEDED

1. Eppendorf tubes (0.5 mL)
2. 2x SDS-PAGE gel loading buffer (e.g. 0.25M Tris-Cl, pH 6.8, 4% SDS, 10% glycerol, 2%  $\beta$ -mercaptoethanol, 0.01% bromophenol blue).
3. DTT (Dithiothreitol) solution (50mM in dH<sub>2</sub>O)
4. Inorganic pyrophosphatase solution (IPP) (100U/mL in 20mM Tris-HCl, pH7.5) (e.g. pyrophosphatase, inorganic, Baker's Yeast, Sigma, I1643)
5. 10X PBS solution (VWR, J373) or 10X TBS solution
6. Tween 20 solution (Sigma, P-5927)

## AUTO-UBIQUINYLATION ASSAY

The protocol set out in this section describes the running of reactions to assess the auto-ubiquitylation activity of the control Ub E3 ligase enzyme provided and user supplied proteins for subsequent analysis by Western blotting.

Hdm2 RING domain (BML-UW0200) is provided as a control ubiquitin E3 ligase for use in auto-ubiquitylation assays.

### Assay protocol

**Note:** recommended total reaction volume = 50µL.

\* Adjust dH<sub>2</sub>O volume in accordance with available Ub E3 ligase protein concentration. A final assay concentration of 150-300nM is recommended as a starting point for Ub E3 ligase auto-ubiquitylation (e.g. use 2.5µL of 6µM Ub E3 ligase protein solution).

\*\* Inorganic pyrophosphatase solution (IPP) is recommended to be included in auto-ubiquitylation reaction but it is not absolutely necessary. Adjust dH<sub>2</sub>O volume if IPP is not included in reaction.

Component	Sample E3-Ub	Sample E3 (-ve control)	Hdm2-Ub (+ve control)	Hdm2 (-ve control)
	Volume/ $\mu$ L			
dH <sub>2</sub> O	21.5*	24.0*	19.0	21.5
10x Ub E3 ligase buffer	5.0	5.0	5.0	5.0
20x Ub E1	2.5	2.5	2.5	2.5
20x E2	2.5	2.5	2.5	2.5
10x Ubiquitin	5.0	5.0	5.0	5.0
20x E3 control (6 $\mu$ M)	-	-	2.5	2.5
*Sample E3 protein	X	X	-	-
50mM DTT	1.0	1.0	1.0	1.0
20x Mg <sup>2+</sup> -ATP	2.5	-	2.5	-
IPP ** (100U/mL)	10	10	10	10

Negative control reactions omitting Mg<sup>2+</sup>-ATP (BML-EW9805-0025) cofactors demonstrate formation of auto-ubiquitylated proteins is ATP-dependent (required for E1 activation, also known as one of the critical characters of ubiquitin activating enzyme E1) and hence derived from the ubiquitin cascade.

### Set-up assays/controls required as follows:

1. Add assay components to 0.5 mL Eppendorf tube(s) in order shown in table above. Keep all enzymes on ice throughout.
2. Mix tube contents gently.
3. Incubate at 37°C for 60 minutes.
4. Quench assays by addition of 50 $\mu$ L 2x SDS-PAGE gel loading buffer followed by heating to 95°C for 5 minutes or 70°C for 10 minutes.  
**Note:** This step removes all Ub thioester linked species (Ub-E1/Ub-E2) so only isopeptide linked Ub-E3 species are detected using ubiquitin antibody/Western blotting.
5. Proceed directly to “Analysis by Western blotting” or store at -20°C until ready.



## ANALYSIS BY WESTERN BLOTTING

### Summary of analysis steps

1. Separate proteins by SDS-PAGE.
2. Western transfer to nitrocellulose or PVDF membrane.

**Note:** Western blotting conditions appropriate for the transfer of large proteins may be required to ensure good transfer of Ubiquitylated-E3 protein to PVDF membrane. For example, use BSN transfer buffer 48mM Tris, pH9.2, 39mM glycine with 10% MeOH and 0.0375% SDS.

3. Block membrane with BSA/PBS-T solution.
4. Probe blot with either:
  - a) ubiquitin antibody supplied or
  - b) appropriate target protein specific primary antibody in conjunction with suitable secondary antibodies.
5. Develop with Western blotting detection reagents.

**Note:** Do NOT use milk in blocking/antibody binding solutions. Please use 5% BSA in PBS-Tween or TBS-Tween instead.

### Materials Needed

1. SDS-PAGE gels - User prepared (10% standard / 4-15% linear gradient)
2. Pre-stained SDS-PAGE molecular weight markers (e.g. See Blue™ Plus 2, Thermo Fisher Scientific, LC5925)
3. PVDF membrane or nitrocellulose membrane (e.g. EMD Millipore, Immobilon-P or GE Healthcare, 10600002)
4. Anti-mouse IgG secondary antibody (HRP linked) (e.g. Goat anti-mouse IgG F(ab')<sub>2</sub>, pAb (HRP conjugate), Prod. No. ADI-SAB-100).
5. (If required) Target protein specific primary antibody (user supplied) and appropriate secondary antibody-HRP conjugate.
6. Western blotting detection reagents (e.g. ECL Reagent).
7. PBS solution 1X PBS.
8. PBS-T solution 1X PBS containing 0.2% Tween 20.
9. BSA/PBS-T blocking solution PBS-T containing 5% bovine serum albumin (BSA).

**Note:** TBS-T can be used as an alternative to PBS-T if required.

**EXAMPLE PROCEDURE FOR WESTERN BLOTTING**

1. Apply 10 - 20 $\mu$ L of each quenched reaction to the SDS-PAGE gel alongside selected molecular weight markers, electrophorese, and transfer protein to nitrocellulose or PVDF membrane according to standard procedures.
2. Remove membrane from the transfer unit and block with BSA/PBS-T blocking buffer for 1 hour at room temperature on a rotor mixer.

**Note:** Drying PVDF membrane prior to blocking, as per Manufacturers' instructions, may considerably enhance results.

3. Check for completion of transfer as indicated by bands of pre-stained protein standard evident on membrane.

**Ubiquitin-conjugate detection**

4. Dilute supplied ubiquitin antibody 1:500 or 1:1000 in BSA/PBS-T.
5. Incubate membrane with ubiquitin antibody solution at room temperature for 1-2 hours or overnight at 4°C on a rotor mixer.
6. Wash membrane for 3 x 10mins with PBS-T on a rocking platform.
7. Dilute appropriate anti-mouse IgG secondary antibody (HRP-linked) according to the manufacturer's instructions (e.g. Prod. No. ADI-SAB-100, 1:3000 in BSA/PBS-T).
8. Incubate membrane with secondary antibody solution for 1 hour at room temperature on a rocking platform, or as specified by the manufacturer.
9. Wash membrane for 4 x 10mins with PBS-T on a rocking platform.
10. (optional) Wash membrane for 1x10mins with 1X PBS (without Tween) on a rocking platform.
11. Proceed to **step 20**.

**Specific target protein detection (if required)**

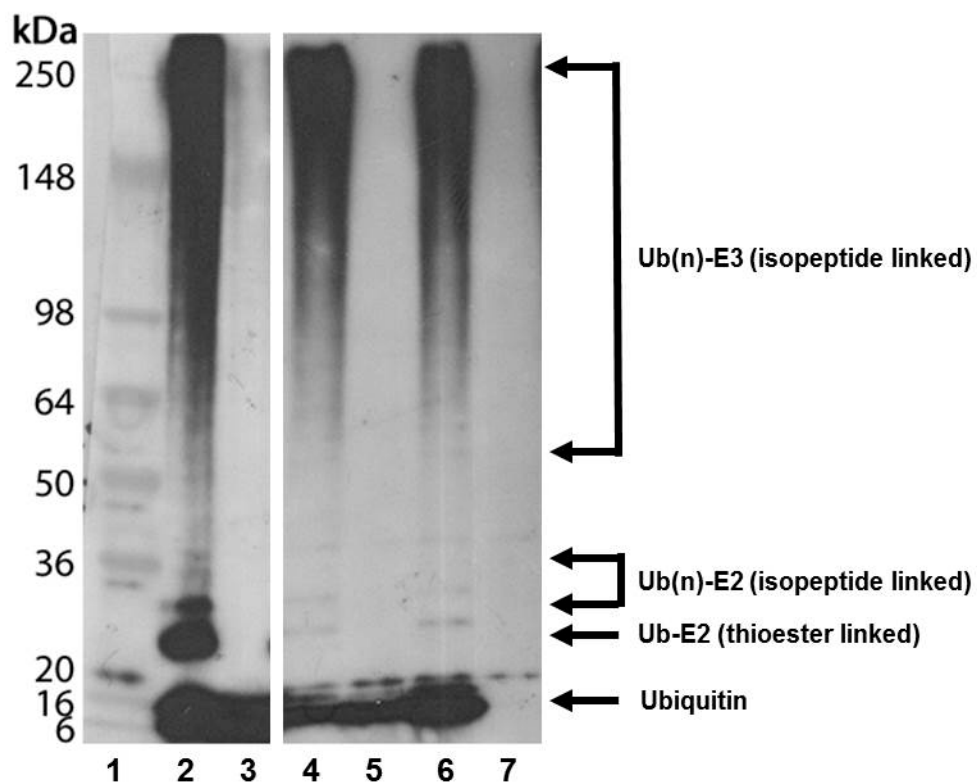
12. Dilute appropriate target protein specific primary antibody according to manufacturer's instructions.
13. Incubate membrane with target protein specific primary antibody solution overnight at 4°C on a rotor mixer.
14. Wash membrane for 3 x 10mins with PBS-T on a rocking platform.

15. Dilute appropriate secondary antibody according to the manufacturer's instructions in BSA/PBS-T.
16. Incubate membrane with secondary antibody solution for 1 hour at room temperature on a rocking platform, or as specified by the manufacturer.
17. Wash membrane for 4 x 10mins with PBS-T on a rocking platform.
18. (optional) Wash membrane for 1x10mins with 1X PBS (without Tween) on a rocking platform.
19. Proceed to **step 20**.

## Analysis

20. Prepare western blotting detection reagent according to the manufacturer's instructions.
21. Incubate membrane with detection reagent for appropriate time.
22. Detect emitted signal by luminography or CCD imaging instrument.

## EXAMPLE RESULTS FOR WESTERN BLOTTING



Lane 1	Pre-stained Protein Standard
Lane 2	Hdm2-Ub, (+) control, (non-reduced)
Lane 3	Hdm2, (-) control (no Mg-ATP), (non-reduced)
Lane 4	Hdm2-Ub, (+) control, (reduced)
Lane 5	Hdm2, (-) control (no Mg-ATP), (reduced)
Lane 6	Hdm2-Ub, (+) control, (reduced)
Lane 7	Hdm2, (-) control (no Ubiquitin), (reduced)

**Figure:** Western blot analysis of control Ub E3 ligase Hdm2 RING domain auto-ubiquitylation assays. Auto-ubiquitylation assays set-up and run as described in “Assay protocol”. Ubiquitylated E3 ligase species were detected by Western blotting as described in “Analysis by western blotting”, using the provided ubiquitin antibody (ADI-SPA-203) at a dilution of 1:1000 dilution.

Results demonstrate auto-ubiquitylation of the control Hdm2 RING domain ligase under the given assay conditions. Please note E2-Ub (thioester linked) is visible in non-reduced condition (left, lane 2), in which  $\beta$ -ME and DTT were omitted from SDS-PAGE gel loading buffer.

**REFERENCES**

1. Haas, A.L. and Siepmann, T. J. Pathways of ubiquitin conjugation. *FASEB J.* 11, 1257-1268 (1997)
2. Hershko, A. and Ciechanover, A. The ubiquitin system. *Annu.Rev.Biochem.* 67, 425-479 (1998)
3. Pickart, C.M. Mechanisms underlying ubiquitination. *Annu.Rev.Biochem.* 70, 503-533 (2001)
4. Strous, G.J. and Govers, R. The ubiquitin-proteasome system and endocytosis. *J.Cell Sci.* 112 ( Pt 10), 1417-1423 (1999)
5. Amemiya, Y., Azmi, P., and Seth, A. Autoubiquitination of BCA2 RING E3 ligase regulates its own stability and affects cell migration. *Mol.Cancer Res.* 6, 1385-1396 (2008)
6. Wu-Baer, F., Ludwig, T., and Baer, R. The UBXN1 protein associates with autoubiquitinated forms of the BRCA1 tumor suppressor and inhibits its enzymatic function. *Mol.Cell Biol.* (2010)

## NOTES

---

## NOTES

---



# Product Manual

## **GLOBAL HEADQUARTERS**

Enzo Life Sciences Inc.  
10 Executive Boulevard  
Farmingdale, NY 11735  
Toll-Free: 1.800.942.0430  
Phone: 631.694.7070  
Fax: 631.694.7501  
info-usa@enzolifesciences.com

## **EUROPE**

Enzo Life Sciences (ELS) AG  
Industriestrasse 17  
CH-4415 Lausen  
Switzerland  
Phone: +41/0 61 926 89 89  
Fax: +41/0 61 926 89 79  
info-ch@enzolifesciences.com

For local distributors and detailed product information visit us online:

[www.enzolifesciences.com](http://www.enzolifesciences.com)