



Ubiquitin Conjugating Kit

(HeLa lysate-based)

Catalog # BML-UW9915



Product Manual

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.**

Table of Contents



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.

Background	2
Kit Description	3
Suggested Application.....	3
Materials Supplied.....	4
Storage	5
Other Materials Needed	5
Ubiquitylation Assay	5
Analysis by Western Blotting	7
Example Procedure for Western Blotting.....	8
Example Results for Western Blotting	10
References.....	11
Contact Information	14

BACKGROUND

The covalent attachment of ubiquitin to proteins in the form of K⁴⁸-linked polyubiquitin chains (of at least four subunits in length) and their subsequent proteasomal degradation plays a fundamental role in the regulation of cellular function through biological events involving cell cycle, differentiation, immune responses, DNA repair, chromatin structure, and apoptosis¹⁻⁴. Attachment of polyubiquitin chains via linkage at alternative lysine residues has been implicated in a wide range of processes including DNA repair, translation, IκB kinase activation and endocytosis⁵. In addition, the attachment of mono-ubiquitin at single or multiple sites on target proteins, or of short ubiquitin chains, has been shown to regulate the location and activity of a diverse range of cellular proteins⁶.

Ubiquitinylation 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 specificity of ubiquitin ligation arises from the subsequent association of the E2-ubiquitin thioester with a substrate specific ubiquitin-protein isopeptide ligase (E3), which facilitates the formation of the isopeptide linkage between ubiquitin and its target protein. The HeLa S100 cytosolic lysate (Prod. No. BML-SW8750) provided contains the full HeLa cell complement of E1, E2s, 20S and 26S proteasome, deubiquitinating, and E3 enzymes. This lysate is ideal for demonstrating ubiquitin-proteasome mediated conjugation/ degradation of radiolabelled or immunodetectable substrates⁷⁻⁹. Addition of deubiquitinating enzyme inhibitors such as ubiquitin aldehyde (Prod. No. BML-UW8450), and proteasome inhibitor, such as epoxomicin (Prod. No. BML-PI127), help to facilitate the accumulation of ubiquitin-protein conjugates.

KIT DESCRIPTION

This kit provides the means of generating ubiquitin-conjugated lysate proteins. The HeLa S100 lysate facilitates controlled ubiquitin conjugation of endogenously-contained substrate proteins of interest through the ubiquitin cascade. Conjugate formation can be detected and monitored by Western blotting using the ubiquitin-conjugate specific antibody supplied and antibodies for specific target proteins. Modified proteins can then be subjected to further purification prior to their use in subsequent experiments if required. Proteasome/deubiquitylating enzyme inhibitor is provided in order to enhance yields of ubiquitin-conjugated proteins. Kit provides sufficient material for 20x50µl reactions.

SUGGESTED APPLICATION

1. Ubiquitylation of specific endogenous HeLa lysate proteins, followed by their immediate detection/analysis using antibodies to the protein(s) of interest, indicating a particular protein is a substrate for the ubiquitin-proteasome pathway.
2. Ubiquitylation of proteins of interest contained in exogenously added expression culture/cell extracts or tissue lysates/extracts, followed by their immediate detection/analysis or isolation/purification for use in subsequent experiments.
3. Modification of proteins using ubiquitin derivatives or mutants for improved detection/analysis or investigation of alternative (non-proteasomal) ubiquitin related pathways in subsequent experiments. For example, biotinylated-ubiquitin (Prod. No. BML-UW8705), for sensitive detection of low level ubiquitin-conjugates or methylated-ubiquitin (Prod. No. BML-UW8555) for mono-ubiquitin conjugate formation.

Note: The HeLa S100 fraction contains a wide range of Ub E3 ligases that may catalyze ubiquitin modification of the full range of available lysine residues of a specific protein of interest. Primary antibodies selected by the user for target protein detection should therefore not be specific for epitopes contain or are neighboring such modification sites.

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

MATERIALS SUPPLIED

1. **HeLa S100 fraction (S100)**, Catalog No. BML-SW8750-0100
 - * Use 10 μ L per 50 μ L reaction
 - * 200 μ L provided (2 x 100 μ L vials)
2. **10 \times Ubiquitin solution (Ub)**, Catalog No. BML-UW8795-0050
 - * Use 5 μ L per 50 μ L reaction
 - * 100 μ L provided (2 x 50 μ L vials)
3. **10 \times ATP-(Energy) Regeneration Solution (ATP-ERS)**, Catalog No. BML-EW9810-0100
 - * Use 5 μ L per 50 μ L reaction
 - * 100 μ L provided (1 x 100 μ L vial)
4. **10 \times Ubiquitylation Buffer**, Catalog No. BML-KW9885-0100
 - * Use 5 μ L per 50 μ L reaction
 - * 100 μ L provided (1 x 100 μ L vial)
5. **Ubiquitin aldehyde** (inhibitor), Catalog No. BML-UW8450-0050
 - * Use 1 μ L per 50 μ L reaction
 - * 50 μ L provided (1 x 50 μ L vial)
6. **Ubiquitin-conjugate specific antibody solution (FK2)**, Catalog No. BML-PW8810-0010
 - * Mono- and polyubiquitylated conjugates mouse mAb (FK2)
 - * 10 μ L of provided (1 x 10 μ L vial)
 - * Dilution of 1:100 to 1:1000 for Western blotting.

STORAGE

Store all components at -80°C for the highest stability. Components labeled with storage temperature other than -80°C can be stored at the temperature listed OR at -80°C. Avoid multiple freeze/thawing.

OTHER MATERIALS NEEDED

1. Microcentrifuge tubes
2. Dimethyl sulfoxide (DMSO)
3. Dithiothreitol (DTT) solution (50 mM in 20 mM Tris-HCl, pH7.5)
4. 2 × SDS-PAGE gel loading buffer For example: 0.25M Tris-HCl, pH 6.8, 4% SDS, 10% glycerol, 2% β-mercaptoethanol, 0.01% bromophenol blue.

UBIQUITINYLATION ASSAY

Two types of reaction are described, using the same basic assay set-up:

- 1) Ubiquitin modification of endogenous (HeLa S100 lysate) proteins
- 2) Ubiquitin modification of specific, exogenously-added target proteins

Ubiquitin modified proteins can then be analyzed immediately subjected to additional purification and/or used in subsequent experiments.

Assay protocol

Note: recommended total reaction volume = 50 μL.

Component	Lysate Protein-Ub	Lysate Protein-Ub (-ve control)
	Volume (μL)	Volume (μL)
dH2O	22	27
10x Ubiquitylation Buffer	5.0	5.0
HeLa S100 fraction	10	10
DTT (50mM)	2.0	2.0
Ubiquitin-aldehyde	1.0	1.0
10x Ubiquitin solution	5.0	5.0
10x ATP-ERS	5.0	-

Negative control reactions omitting ATP-ERS cofactor demonstrates that formation of ubiquitylated target protein/increased levels of endogenous ubiquitylated proteins is ATP dependent (required for E1 activation) and, hence, derived from the ubiquitin cascade.

Note: Suggested concentrations and assay incubation times are given as a starting point for such reactions and will require optimization by end-user.

Note: Incubation times will require optimization for ubiquitylation of specific target proteins. It is recommended to run a time course as an initial experiment and analyze with FK2 Ub-conjugate (BML-PW8810)/target protein specific antibodies as described in “Analysis by Western Blotting” to identify optimal incubation times.

Set-up assays/controls required as follows:

1. Add assay components to 0.5 mL 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 0.5-4 hours.
4. Store assay solutions at –20°C prior to additional processing/use in subsequent experiments.

For analysis of all/part of the crude assay solution by Western blotting:

5. Quench assays by addition of 50 µL 2x SDS-PAGE gel loading buffer followed by heating to 95°C for 10 minutes.
6. Proceed directly to “Analysis by Western blotting” or store at –20°C until ready

Note: DO NOT quench whole assay solution if additional purification or subsequent use of active/native ubiquitin-modified proteins is required.

ANALYSIS BY WESTERN BLOTTING

Summary of analysis steps

1. Separate proteins by SDS-PAGE.
2. Western transfer to PVDF membrane.
3. Block membrane with BSA/PBS-Tween solution.
4. Probe blot with either
 - i) FK2 ubiquitin-conjugate specific antibody supplied (BML-PW8810), or
 - ii) Appropriate target protein specific primary antibody, followed by a suitable secondary antibody.
5. Develop with Western Blotting detection reagents.

Note: Do not use milk in blocking/antibody binding solutions. 1-2% BSA in PBS or PBS-Tween is recommended.

Materials Needed

1. SDS-PAGE gels - User prepared or precast polyacrylamide gel (12% standard / 4-15% gradient).
2. Pre-stained SDS-PAGE molecular weight markers (e.g., SeeBlue® Plus2 Pre-stained Protein Standard)
3. PVDF membrane (e.g., Immobilon-P)
4. Target protein specific primary antibody.
5. Anti-mouse secondary antibody (HRP linked) for use with Ub-conjugate primary antibody FK2 (e.g., Goat anti-mouse IgG F(ab')₂, polyclonal antibody (HRP conjugate), Prod. No. ADI-SAB-100, or Anti-Mouse Polyvalent Immunoglobulins (G,A,M)-Peroxidase antibody produced in goat, Sigma, A0412).
6. Appropriate secondary antibody-HRP conjugate for use with chosen target protein specific primary antibody.
7. Western blotting detection reagents (e.g., ECL Reagent).
8. PBS solution 1X PBS.
9. PBS-Tween solution 1X PBS containing 0.2% Tween 20.
10. BSA/PBS-T blocking solution PBS-T containing 1-2% bovine serum albumin (BSA).

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

EXAMPLE PROCEDURE FOR WESTERN BLOTTING

Note: This protocol has been optimized using the materials indicated above. Using materials other than those listed may require additional optimization.

1. Apply ~10 μ L of each quenched reaction to the SDS-PAGE gel alongside selected molecular weight markers, electrophorese, and transfer protein to 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. Wash membrane for 3 x 10mins with PBS-T on a rocking platform at room temperature.

Detection

4. Dilute supplied FK2 ubiquitin-conjugate reactive primary antibody 1:100-1:1000 in BSA/PBS-T. (Alternatively dilute appropriate target protein specific primary antibody according to manufacturer's instructions (e.g., p53, monoclonal antibody (DO-1), at 1:1000)).
5. Incubate membrane with primary antibody solution for 1 hour at room temperature on a rocking platform, or overnight at 4°C.
6. Wash membrane for 3 x 10mins with PBS-T on a rocking platform.
7. Dilute appropriate secondary antibody according to the manufacturer's instructions (e.g., Anti-Mouse Polyvalent Immunoglobulins (G,A,M)-Peroxidase antibody produced in goat (Sigma, A0412) diluted 1:1000-1:5000 in BSA/PBS-T, following Ub-conjugate specific primary antibody)
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 6 x 10mins with PBS-T on a rocking platform.
10. Proceed to **step 17**.

Specific target protein detection (if required)

11. Dilute appropriate target protein specific primary antibody according to manufacturer's instructions.
12. Incubate membrane with target protein specific primary antibody solution overnight at 4°C on a rotor mixer.
13. Wash membrane for 3 x 10mins with PBS-T on a rocking platform.
14. Dilute appropriate secondary antibody in BSA/PBS-T according to the manufacturer's instructions.
15. Incubate membrane with secondary antibody solution for 1 hour at room temperature on a rocking platform, or as specified by the manufacturer.
16. Wash membrane for 6 x 10mins with PBS-T on a rocking platform.

Analysis

17. Prepare Western Blotting detection reagent according to the manufacturer's instructions.
18. Incubate membrane with detection reagent for appropriate time.
19. Detect emitted signal by luminography or CCD imaging instrument.

EXAMPLE RESULTS FOR WESTERN BLOTTING

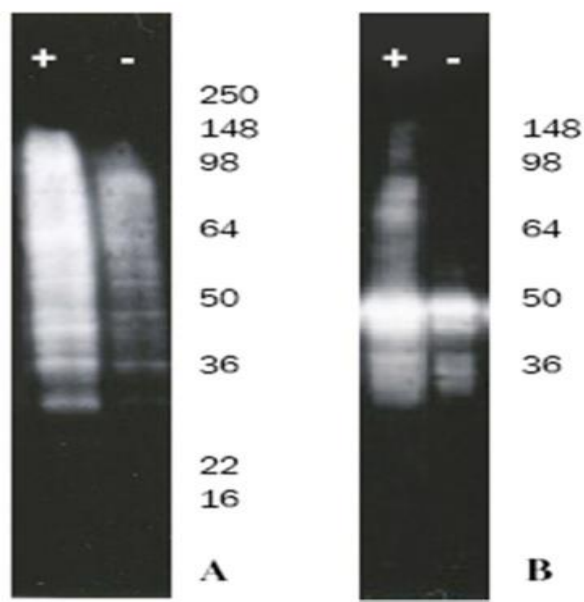


Figure: Western Blot of S100 ubiquitin conjugation assays of both endogenous lysate and exogenously added p53 proteins. Assays set-up and run as described in “Assay protocol” section. Ubiquitin-protein conjugate formation was detected by Western blotting of assays for A: General ubiquitinylation of endogenous HeLa S100 lysate proteins using the supplied FK2 ubiquitin-conjugate specific antibody (BML-PW8810) or B: specific modification of p53 present in HeLa S100 lysate using p53 specific monoclonal antibody DO-7 (BML-PW1095) as described in “Analysis by Western blotting”. Results demonstrate the utility of the HeLa conjugation kit for both the ubiquitin modification of endogenous HeLa S100 lysate proteins in general and of specific endogenous proteins of interest, such as p53. The elevated level (A) or formation (B) of ubiquitin modified proteins can be clearly seen in the +ve (ATP containing) assays. The lower level (A) or absence (B) of ubiquitin conjugated proteins in –ve control reactions (-ATP) demonstrates that their formation is ATP-dependent (required for E1 activation) and, hence, derived from the ubiquitin cascade.

REFERENCES

1. Hershko, A., and Ciechanover, A. "The ubiquitin system". *Annu. Rev. Biochem.* 67, 425–479, (1998)
2. Haas, A. L., and Siepmann, T. J. "Pathways of ubiquitin conjugation". *FASEB J.* 11, 1257–1268, (1997)
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, 1417–1423, (1999)
5. Pickart, C. M., and Fushman, D. "Polyubiquitin chains: polymeric protein signals". *Curr. Opin. Chem. Biol.* 8, 610-616, (2004)
6. Hicke, L. "Protein regulation by monoubiquitin". *Nat. Rev. Mol. Cell Biol.* 2, 195-201, (2001)
7. Hershko, A., Heller, H., Elias, S., and Ciechanover, A. "Components of ubiquitin-protein ligase system. Resolution, affinity purification, and role in protein breakdown". *J. Biol. Chem.*, 258, 8206-8214, (1983)
8. Hershko, A., Heller, H., Eytan, E., and Reiss, Y. "The protein substrate binding site of the ubiquitin-protein ligase system". (1986) *J. Biol. Chem.*, 261, 11992-11999, (1986)
9. Ganoth, D., Leshinsky, E., Eytan, E., and Hershko, A. "A multicomponent system that degrades proteins conjugated to ubiquitin. Resolution of factors and evidence for ATP-dependent complex formation". *J. Biol. Chem.*, 263, 12412-12419, (1988)
10. Cooper, H. J., Heath, J. K., Jaffray, E., Hay, R. T., Lam, T. T., and Marshall, A.G. "Identification of Sites of Ubiquitination in Proteins: A Fourier Transform Ion Cyclotron Resonance Mass Spectrometry Approach". *Anal. Chem.*, 76, 6982-6988 (2004)



Product Manual

NOTES



Product Manual

NOTES



Product Manual

**Global Headquarters
Enzo Life Sciences Inc.**

10 Executive Blvd
Farmingdale, NY 11735
(p) 1-800-942-0430
(f) 1-631-694-7501
(e) info-usa@enzolifesciences.com

Enzo Life Sciences (ELS) AG

Industriestrasse 17, Postfach
CH-4415 Lausen/ Switzerland
(p) +41/0 61 926 89 89
(f) +41/0 61 926 89 79
(e) info-ch@enzolifesciences.com

For local distributors and detailed product information visit us online:

www.enzolifesciences.com