

Ubiquitin conjugating kit (HeLa lysate-based)

For the generation of ubiquitin conjugated proteins. HeLa S100 lysate facilitates controlled ubiquitin conjugation of substrate proteins (exogenous or endogenous) of interest through the ubiquitin cascade. Conjugate formation can be detected and monitored by Western blotting using the highly sensitive ubiquitin-conjugate specific antibody supplied and/or antibodies for specific target proteins. Modified proteins can be subjected to further purification prior to their use in subsequent experiments if required. Suggested uses: Generation of ubiquitin conjugated proteins. Exogenous or endogenous HeLa lysate proteins (tagged/radio-labeled/immuno-detectable) can be ubiquitinated followed by immediate detection/analysis. Subsequent analysis could include proteasomal degradation, ubiquitin modification site mapping (by mass spectrometry), and the effect of ubiquitin modification on enzyme interactions, activity and function, Ubiquitination of proteins of interest from cell or tissue extracts, Modification of proteins using ubiquitin derivatives or ubiquitin mutants for improved detection, analysis or investigation of alternative (non-proteasomal) ubiquitin signaling pathways.

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-UW9915-0001	20 tests
-----------------	----------

Manuals, SDS & CofA

[View Online »](#)

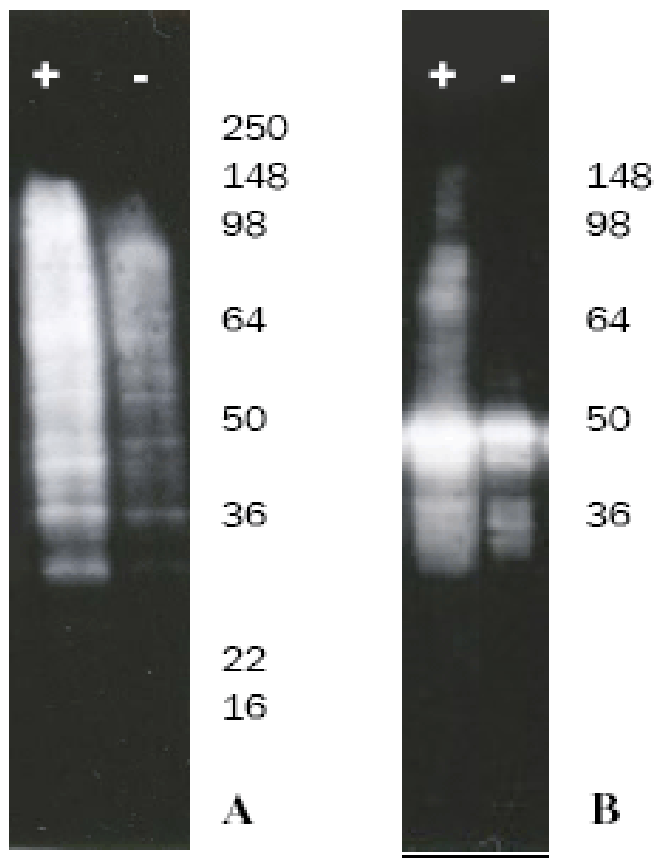


Figure: Western blot of S100 ubiquitin conjugation assays of both endogenous lysate and exogenously added p53 proteins. Assays set-up and run as described in “Assay protocol” section. Ubiquitin-protein conjugate formation was detected by Western blotting of assays for A: General ubiquitinylation of endogenous HeLa S100 lysate proteins using the supplied FK2 ubiquitin-conjugate specific antibody (PW8810) or B: specific modification of p53 present in HeLa S100 lysate using p53 specific monoclonal antibody DO-7 (Prod. No. 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 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.

Handling & Storage

Use/Stability	All kit components should be stored at -80°C to ensure stability and activity. Inhibitors are stable for up to 3 months at -20°C when dissolved in DMSO.
Handling	Avoid freeze/thaw cycles.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application Notes	Uses: <ol style="list-style-type: none">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.
Contents	<p>HeLa S100 fraction (Prod. No. BML-SW8750), 1 mg (200µl provided as 2×100µl vials), 5mg/ml</p> <p>10x Ubiquitin solution (UB) (Prod. No. BML-UW8795), 100µl (provided as 2X50µl vials), 1mg/ml</p> <p>10x ATP-(Energy) regeneration solutions (ATP-ERS) (Prod. No. BML-EW9810), 100µl</p> <p>10x Ubiquitylation buffer (Prod. No. BML-KW9885), 100µl</p> <p>Ubiquitin aldehyde (inhibitor) (Prod. No. BML-UW8450), 50 µl, 1mg/ml</p> <p>Ubiquitin-conjugate specific antibody solution (UBCJ2) (Prod. No. ENZ-ABS840), 10 µl, Mono- and polyubiquitylated conjugates, mAb</p>
Quantity	Provides sufficient material for 20 x 50µl 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-
fr@enzolifesciences.com](mailto:info-fr@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)