

Ubiquitin activating kit

For activation of ubiquitin for use in ubiquitinylation experiments. The kit provides the means of generating thioester linked, activated ubiquitin-E1 conjugates, utilizing the first step in the ubiquitin cascade, for investigation of ubiquitin activation, subsequent ubiquitin transfer to/interaction with E2 conjugating enzymes, and their use in the ubiquitinylation of E3 ligases and target substrate proteins. Ubiquitin antibody provided allows high sensitivity detection of ubiquitin and ubiquitin-conjugates.

The covalent attachment of ubiquitin to proteins (ubiquitinylation) 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 apoptosis. 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, in the presence of an Mg^{2+} cofactor, followed by its subsequent transfer to the active site cysteine of an 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

Citations: 5

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-UW0400A-0001

20 tests

Manuals, SDS & CofA

[View Online »](#)

- For use with wild type or mutant E2 enzymes in E1 initiated/mediated reactions
- Contains highly sensitive ubiquitin antibody for detection of ubiquitin and ubiquitin conjugates

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Notes

Uses:

1. Activation of Ubiquitin for conjugation to wild type or mutant E2 enzymes (user supplied) via thioester bond formation.
2. Use of Ub-E2 conjugates produced in the subsequent ubiquitin modification of specific target proteins in presence of a dedicated ubiquitin E3 ligase.
3. Investigation of ubiquitin activation by E1 activating enzyme.
4. Substitution of the wild type ubiquitin provided with ubiquitin mutants or derivatives (e.g. biotinylated-ubiquitin, Prod. No. BML-UW8705) allowing their activation and subsequent utility in the ubiquitin cascade.

Contents

20X Ubiquitin Activating Enzyme Solution (E1)
20X Ubiquitin Solution (Ub)
Ubiquitin Antibody Solution
20X Mg-ATP Solution
2X Non-reducing Gel Loading Buffer
10X Ubiquitylation Buffer

Quantity

20 tests

UniProt ID

P0CG47 (UBB), P0CG48 (UBC), P62979 (RPS27A),
P62987 (UBA52)

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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