

Uba6 (human), (recombinant)

UBA6 (ubiquitin-like modifier activating enzyme-6, also known as UbE1L2) is a novel ubiquitin activating enzyme that is also capable of activating the ubiquitin-like protein FAT10 (human leukocyte antigen F-associated transcript 10), which, similarly to ubiquitin, may serve as a signal for proteasomal degradation.

UBA6 exhibits 42% homology to the major ubiquitin activating enzyme UbE1 and contains all the key structural elements required for E1 enzyme activity. It has been shown to interact with a number of E2 and E3 enzymes, facilitating transfer of ubiquitin to substrate proteins such as p53 *in vitro*. Only one UBA6 specific E2 conjugating enzyme, known as Use1 or UbE2Z, which is capable of accepting activated FAT10, has been identified to date.

FAT10, is encoded by the major histocompatibility (MHC) class I locus, and its expression is induced by tumor necrosis factor α (TNF α) and interferon- γ (IFN γ). FAT10 is highly upregulated in hepatocellular carcinoma as well as in other gastrointestinal and gynecological cancers, though its precise biochemical and cellular functions have yet to be determined.

FAT10 forms covalent conjugates with cellular proteins through its C-terminal diglycine motif, though substrates have yet to be identified. Knockdown of Uba6 results in strong, but not complete diminution of FAT10 conjugation, an effect not observed with knockdown of UbE1. This finding indicates that Uba6 is both necessary and sufficient to activate FAT10 and thereby facilitate its conjugation.

Citations: 1

[View Online »](#)

Ordering Information

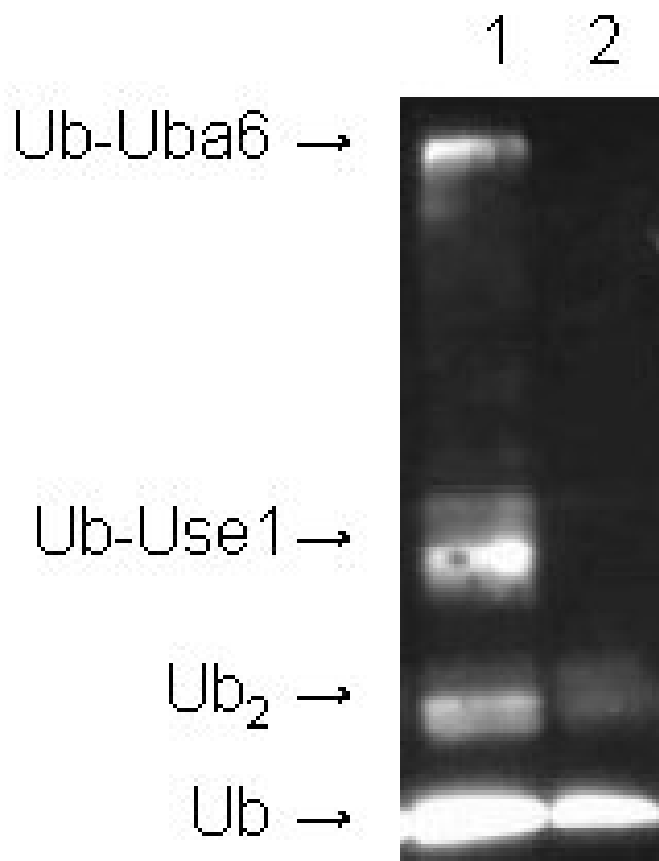
[Order Online »](#)

BML-UW0350-0050

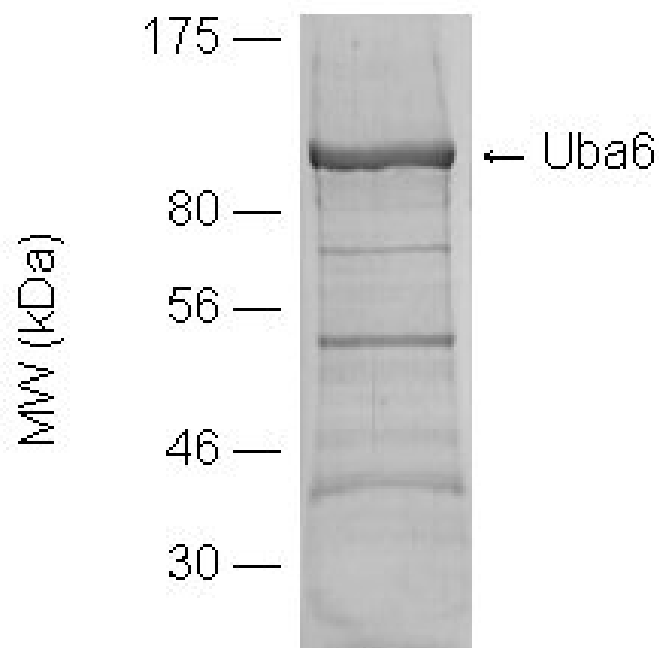
50 μ g

Manuals, SDS & CofA

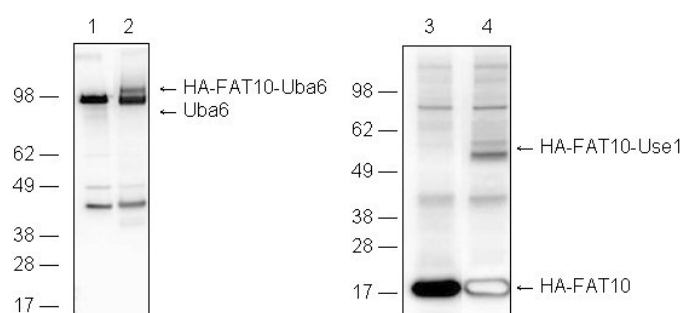
[View Online »](#)



Western blot of ubiquitin thioester assay showing Ub activation by Uba6 (BML-UW0350) and conjugation by Use1 in the presence (1) or absence (2) of Mg-ATP. Ubiquitin species detected using ubiquitin antibody BML-UG9510 at 1:1000 dilution.



SDS-PAGE image of Uba6 (3µg, BML-UW0350), coomassie stained.



Western blot of FAT10 thioester assay showing HA-FAT10 activation by Uba6 (BML-UW0350, 1-2) and conjugation by Use1 (3-4) in the absence (1,3) or presence (2,4) of Mg-ATP. Species of interest detected using either Uba6 antibody (BML-PW0525, 1-2) or anti-HA (3-4) at 1:1000 dilution.

Handling & Storage

Use/Stability	Stable for at least 6 months after receipt when stored at -80°C.
Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -80°C.
Short Term Storage	-80°C
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Ubiquitin-activating enzyme E1-like protein 2, Monocyte protein 4
Application Notes	For use in ubiquitin and FAT10 protein modification studies.
Formulation	Liquid. In 40mM Tris-HCl, pH 8.0, 110mM NaCl, 2.2mM KCl, 80 µg/mL FLAG peptide, 20% glycerol and 3mM DTT.
MW	~118kDa
Purity	≥65% (SDS-PAGE)
Source	Produced in baculovirus system
UniProt ID	A0AVT1

Last modified: May 29, 2024



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)