

Otubain-1 (human), (recombinant) (His- tag)

Otubain-1 belongs to the ovarian tumor (OTU) domain class of cysteine protease deubiquitinating enzymes and has been implicated in mediating lymphocyte antigen responsiveness and may be generally involved in RNA processing and cell adhesion/morphology. Human otubain-1 has been shown to exhibit high linkage-specificity in vitro, cleaving Lys⁴⁸ (K⁴⁸)-linked polyubiquitin but not K⁶³-, K²⁹-, K⁶-, or K¹¹-linked polyubiquitin, or linear α -linked polyubiquitin. Cleavage is not limited to either end of a polyubiquitin chain, and both free and substrate-linked polyubiquitin are disassembled.

Ordering Information

[Order Online »](#)

BML-UW0680-0100	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)

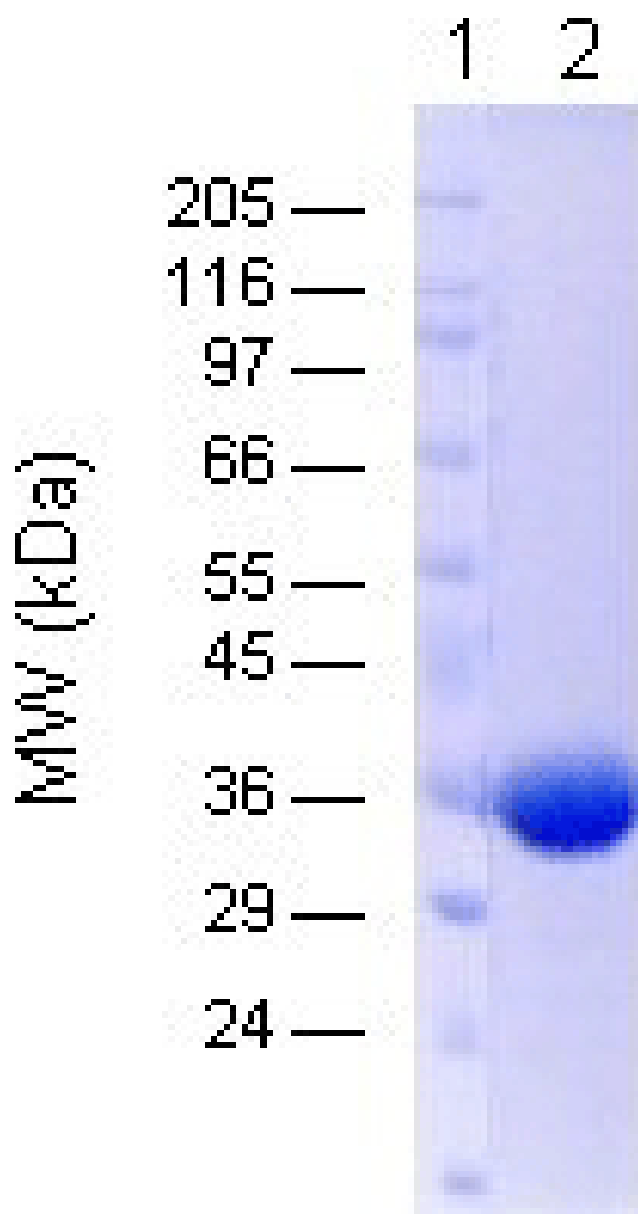


Figure 1: SDS-PAGE image of His₆-Otubain-1 (Coomassie). Lane 1: MW markers. Lane 2: Otubain-1 (human) (recombinant) (His) (Prod. No. BML-UW0680) (8μg).

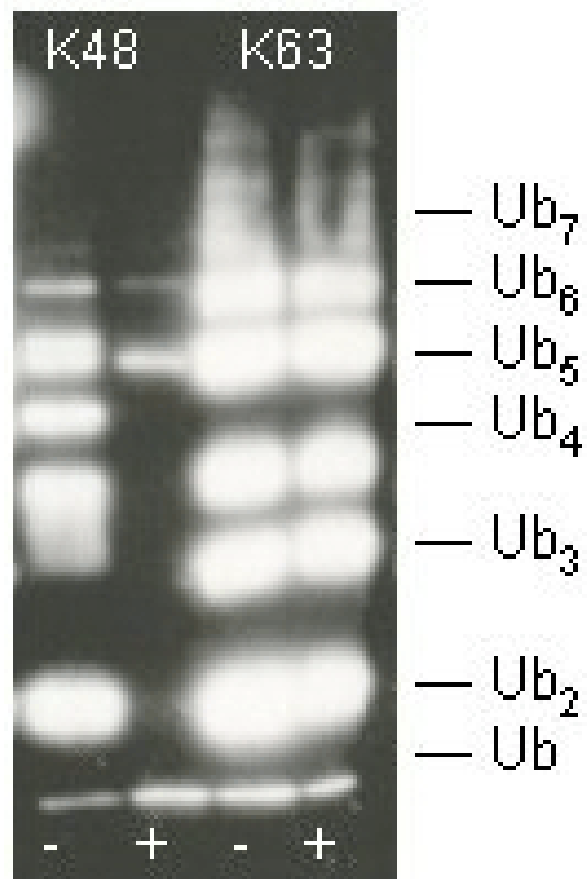


Figure 2: Deubiquitylation assay: Western Blot showing polyubiquitin chains (K⁴⁸-linked, Prod. No. BML-UW8860; K⁶³-linked, Prod. No. BML-UW9570) incubated with (+) or without (-) Otubain-1 (Prod. No. BML-UW0680)

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	OTUB1, Ubiquitin thioesterase OTUB1, Ubiquitin-specific-processing protease OTUB1, Deubiquitinating enzyme OTUB1
------------------	--

Application Notes

Uses:

1. Deubiquitinylation assays using polyubiquitin chains or ubiquitin-protein conjugates.
2. Studies of the K⁴⁸-linkage specific polyubiquitin chain binding and deubiquitinylation.
3. Investigation of the role of Otubain-1 in various biological processes.
4. Development of inhibitors for Otubain-1 deubiquitinylation activity.

Formulation	Liquid. In 50mM TRIS, pH 7.5, containing 50mM NaCl and 1mM DTT.
-------------	---

MW	~33kDa
----	--------

Purity	≥95% (SDS-PAGE)
--------	-----------------

Source	Produced in <i>E. coli</i> .
--------	------------------------------

UniProt ID	Q96FW1
------------	--------



Last modified: May 29, 2024
usa@enzolifesciences.com

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

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@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