

Ubiquitin (human), (recombinant)

Ubiquitin is a 76 amino acid protein, which when covalently linked to proteins as a branched chain, serves as a tag marking proteins for a variety of cellular activities. Its primary sequence is remarkably conserved, with only three amino acid changes between the yeast and human sequences. Originally isolated from thymus and assumed to be a thymic hormone its primary role is now recognised as being in ATP-dependent protein degradation.

Citations: 15

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-UW8795-0005	5mg
-----------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	White powder.
CAS	79586-22-4
Formulation	Lyophilized.
MW	8565 Da
Purity	≥98% (SDS-PAGE)
Solubility	Soluble in water
Source	Produced in <i>E. coli</i> .
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](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-
be@enzolifesciences.com](mailto:info-be@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)