

Ubiquitin, (biotinylated)

Biotinylated ubiquitin (prepared by the action of biotinamidocaproic acid N-hydroxysuccinimide ester on ubiquitin) is useful for studying ubiquitinylation in vitro, where it represents an alternative to radio-iodinated ubiquitin. This product contains a number of multiply biotinylated ubiquitin species, which has proven itself to be a reagent of considerable utility particularly for carrying out E1/E2-mediated thiolester assays.

Citations: 3

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-UW8705-0100	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Appearance	White solid.
Application Notes	Biotinylated ubiquitin is activated by E1 and is a competent substrate for E2 and E3 enzymes. The biotinyl moiety also provides a potential route to affinity purification of labeled species.
Formulation	Lyophilized from Acetonitrile:water (1:1).
Purity Detail	Mass spectral analysis shows presence of [biotinyl]1-5-ubiquitin with [biotinyl]2-4-ubiquitin being the predominate species.
Solubility	In order to ensure full solubility it is recommended that the material be dissolved in 5µL DMSO prior to dilution with the buffer of choice to the desired working concentration.
Source	Produced in <i>E. coli</i> .
UniProt ID	P0CG47 (UBB), P0CG48 (UBC), P62979 (RPS27A), P62987 (UBA52)



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)