

Ubiquitin (human), (recombinant) (GST- tag)

GST-ubiquitin is activated by E1 and is a competent substrate for the E2 and E3 enzymes.

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-UW8620-0001	1mg
-----------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Notes

GST-ubiquitin is useful for studying ubiquitinylation *in vitro* and for the detection and purification of ubiquitinylated substrates.

Formulation

Liquid. In TBS, pH 7.5.

MW

38.5 kDa

Purity

≥98% (SDS-PAGE)

Source

Produced in *E. coli*.

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)