

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

Autophagy is a highly conserved bulk protein degradation pathway responsible for the turnover of long-lived proteins, disposal of damaged organelles, and clearance of aggregate-prone proteins. Autophagy is involved in various physiological or pathological processes, such as development, host defense response, cancer and neuronal degenerative diseases. It involves concerted action of more than 20 specific autophagy (ATG) proteins that mediate the formation of a double-membrane vesicle, the autophagosome, which engulfs the substrate and delivers it to the lysosome for degradation.

Formation and expansion of the pre-autophagosomal structure in yeast requires the attachment of the ubiquitin-like protein ATG8 *via* its C-terminal glycine to the amino group of phosphatidylethanolamine (PE), enabling its anchoring to the isolation membrane of the autophagosome. In mammals, ATG8 is represented by at least seven related proteins that fall into two subgroups, LC3- and GABARAP-like proteins. The free and PE linked versions of these proteins are often referred to as LC3-I and LC3-II respectively.

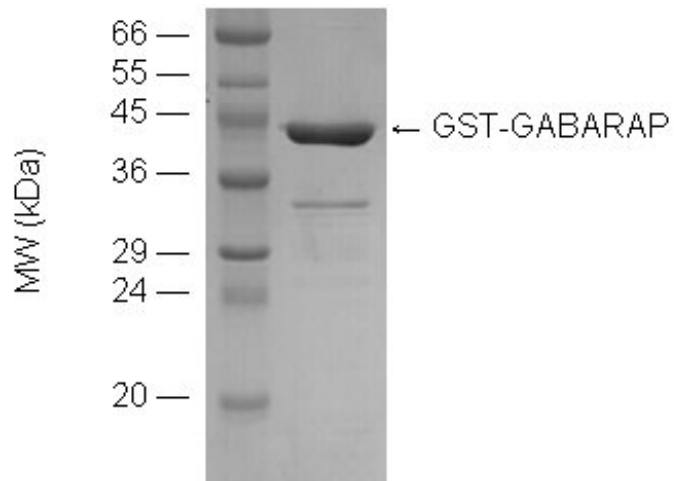
Ordering Information

[Order Online »](#)

BML-UW1175-0500	500µg
------------------------	-------

Manuals, SDS & CofA

[View Online »](#)



SDS-PAGE image (Coomassie stained) of GST-GABARAP (2.5 μ g, BML-UW1175).

Handling & Storage

Use/Stability	Stable for at least 12 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	Gamma-aminobutyric acid receptor-associated protein
Application Notes	For use in general and selective autophagy studies.
Formulation	Liquid. In PBS.
MW	~42.1kDa
Purity	≥90% (SDS-PAGE)
Source	Produced in <i>E. coli</i> .
UniProt ID	O95166



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)