

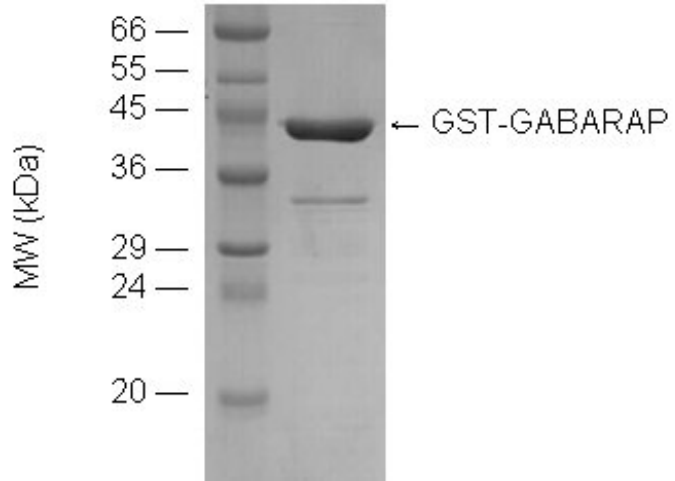
# GABARAP (human), (recombinant) (GST- tag)

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.

**Manuals, SDS & CofA**

[View Online »](#)



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

## Handling & Storage

<b>Use/Stability</b>	Stable for at least 12 months after receipt when stored at -80°C.
<b>Handling</b>	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -80°C.
<b>Short Term Storage</b>	-80°C
<b>Long Term Storage</b>	-80°C
<b>Shipping</b>	Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

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

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)