

# Blocking peptide for Glutamic acid decarboxylase mAb (11) (Prod. No. BML- GC3208)

Immunization and blocking peptide for Glutamic Acid Decarboxylase (GAD) antibody (BML-GC3208). The tetradecapeptide corresponds to the N-terminal residues 4-17 of rat GAD65. The sequence is wholly conserved human GAD65.

## Ordering Information

[Order Online »](#)

|                 |      |
|-----------------|------|
| BML-GP3209-0050 | 50µg |
|-----------------|------|

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

**Use/Stability** As indicated on product label or CoA when stored as recommended.

**Long Term Storage** -20°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Appearance** Lyophilized solid.

**MW** 1416.43

**Purity** ≥99%

**Sequence** Pro-Gly-Ser-Gly-Phe-Trp-Ser-Phe-Gly-Ser-Glu-Asp-Gly-Ser

**Solubility** Soluble in Water.

**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)