

# Glutamic acid decarboxylase monoclonal antibody (11)

This antibody is covered by our [Worry-Free Guarantee](#).

## Ordering Information

[Order Online »](#)

BML-GC3208-0100	100µl
-----------------	-------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage      -20°C

Shipping      Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name      GAD

Application      IHC, WB

Formulation      Liquid. In PBS containing 10mM sodium azide.

Host      Mouse

Immunogen      The hybridoma secreting the antibody to GAD65 was generated by fusion of splenocytes from a non-obese diabetic (NOD) mouse which had received a single subdiabetogenic injection of streptozotocin. Effectively, Prod. No. BML-GC3208 is a mouse autoantibody to GAD.

Isotype      IgG1

Purity Detail      Partially purified.

Species Reactivity      Human, Monkey, Mouse, Rat

Specificity      Recognizes an N-terminal epitope. The antibody consistently labels neuronal GAD in the CNS and GAD-containing cells in mammalian (not mouse) pancreas.

UniProt ID      P48320

Worry-free Guarantee      This antibody is covered by our [Worry-Free Guarantee](#)

.

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)