

Blocking peptide for Glutamic acid decarboxylase pAb (Prod. No. BML- GC3008)

Immunization and blocking peptide for Glutamic Acid Decarboxylase (GAD) antibodies. The pentadecapeptide corresponds to the C-terminal residues [Cys+]572-585 of rat GAD65. The sequence is wholly conserved at the C-terminus of rat GAD67, mouse, pig and human GAD65/67, cat GAD67, and in human GAD-2.

Ordering Information

[Order Online »](#)

BML-GP3009-0050	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	1790.9
Sequence	Cys-Asp-Phe-Leu-Ile-Glu-Glu-Ile-Glu-Arg-Leu-Gly-Gln-Asp-Leu
Solubility	Soluble in Water.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com