

Glutamate decarboxylase (65kDa isoform) (human) monoclonal antibody (83/2E6)

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

Citations: 4

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-803-054-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	Stable for at least 1 year when stored at +4°C.
Handling	Avoid freeze/thaw cycles.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	GAD65, GAD2
Application	ICC, IHC (FS), IP, Radioimmunoassay
Application Notes	GAD-binding of this antibody is competitively inhibited by human GAD-antibodies.
Clone	83/2E6
Crossreactivity	Does not cross-react with GAD67.
Formulation	Liquid. In PBS. pH 7.4, containing 0.9mg/ml sodium azide.
Host	Mouse
Immunogen	Recombinant human GAD65 (glutamate decarboxylase (65kDa isoform)).
Isotype	IgG1
Purity Detail	Ammonium sulphate precipitation.
Species Reactivity	Human, Rat

Specificity

Recognizes an epitope (aa 221-442) of human GAD65.
Binds to human and rat pancreas tissue but not mouse pancreas tissue. Not detectable on Bouin-fixed pancreatic tissues. No binding to viable rat islet cells.

UniProt ID

Q05329

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)