

Glu-Glu monoclonal antibody (Glu-Glu)

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

Ordering Information

[Order Online »](#)

| | |
|-----------------|-------|
| ENZ-ABS161-0500 | 0.5ml |
|-----------------|-------|

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application IF, IP, WB

Application Notes This antibody is effective in immunofluorescence, immunoblotting, and immunoprecipitation of tagged proteins.

Clone Glu-Glu

Formulation Liquid. Phosphate-buffered solution + 0.03% Thimerosal.

Host Mouse

Immunogen Peptide sequence CEEEEYMPME derived from the polyoma virus medium T antigen.

Isotype IgG1κ

Purity Detail Purified by affinity chromatography.

Recommendation Western Blot (1:1,000)Immunofluorescence (1:1,000)Immunoprecipitation
Dilutions/Conditions (1:150)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

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



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)