

GAPDH monoclonal antibody (GA1R) (biotin conjugate)

Glyceraldehyde 3 phosphate dehydrogenase (GAPDH) is an enzyme involved in glycolysis. In addition to this role, it is involved in other cellular processes ranging from membrane fusion, neuronal apoptosis to cancer. GAPDH is expressed at high levels in most tissues and is therefore useful as protein loading control in Western blot analysis. Additionally, the antigen is conserved in most eukaryotes.

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

Citations: 2

[View Online »](#)

Ordering Information

[Order Online »](#)

ENZ-ABS277-0050	50µl
-----------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling Do not freeze.

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Glyceraldehyde-3-phosphate dehydrogenase

Application ELISA, Flow Cytometry, ICC, IF, IHC (PS), WB

Clone GA1R

Formulation Liquid. In PBS containing a proprietary stabilizer and 0.02% sodium azide.

Host Mouse

Immunogen Full-length recombinant GAPDH.

Isotype IgG1

Purity Detail Purified

Species Reactivity Chicken, E. coli, Hamster, Human, Insect, Mouse, Rabbit, Rat, Yeast

Specificity Recognizes GAPDH in multiple species, including human, mouse, rat, bacteria, chicken, hamster, insect, rabbit and yeast.

UniProt ID P04406

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-
fr@enzolifesciences.com](mailto:info-fr@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)

Last modified: May 29, 2024