

GATA3 polyclonal antibody

This polyclonal antibody detects GATA3, a transcription factor involved in T-cell differentiation and epithelial development.

GATA3 is a transcription factor of the GATA family that regulates gene expression in various tissues, particularly in immune cells and breast epithelial cells. It is a critical regulator of T-helper 2 (Th2) differentiation and is commonly used as a biomarker in breast cancer and urothelial carcinoma.

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

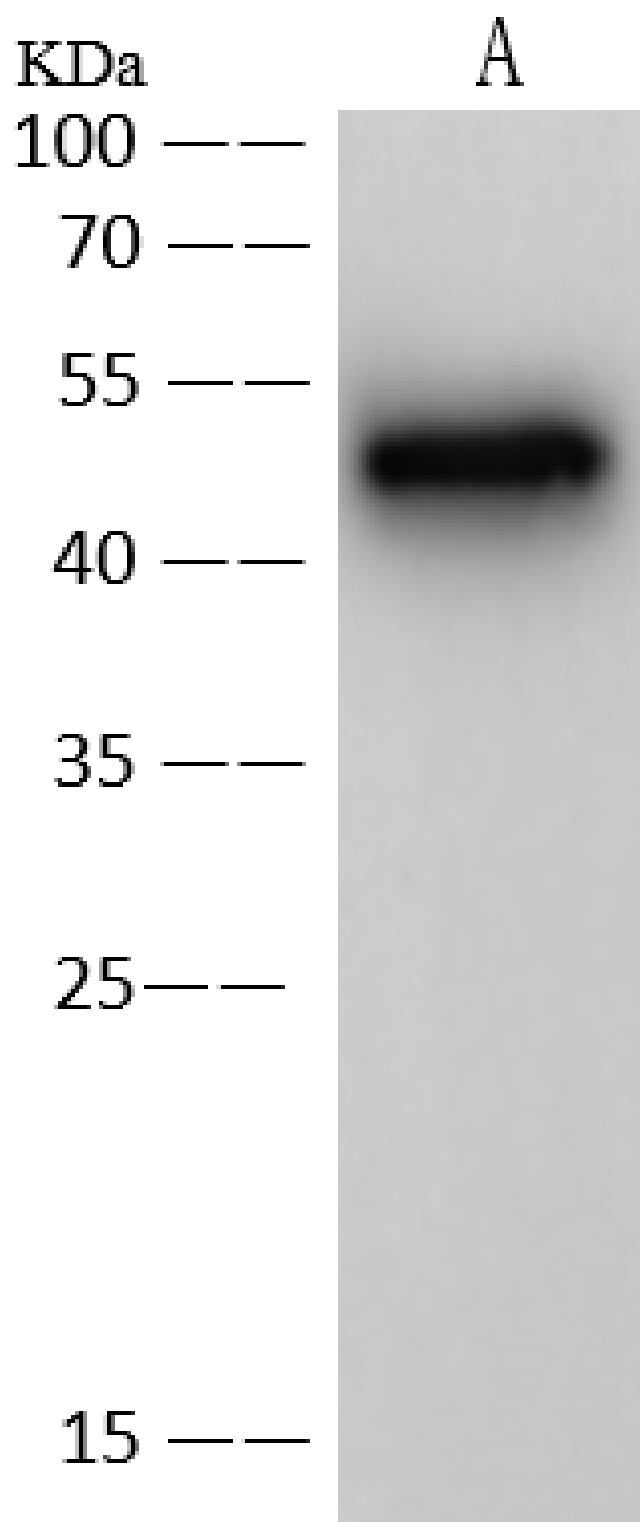
Ordering Information

[Order Online »](#)

ENZ-ABS1064-0100	100µl
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of mouse brain tissue using GATA3 polyclonal antibody (ENZ-ABS1063) at 1:500 dilution. Lane A: Mouse brain tissue lysate at 30 μ g per lane. Secondary antibody: Goat Anti-Rabbit IgG (H+L)/HRP at 1:10'000 dilution. Observed band size: 48 kDa

Handling & Storage

Handling Avoid freeze-thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Trans-acting T-cell-specific transcription factor GATA-3, GATA-binding factor 3

Application IP, WB

Crossreactivity Predicted to cross-react with mouse based on 100% sequence homology.

Formulation Liquid. 0.2 µm filtered solution in PBS. Carrier-free. Preservative-free.

Host Rabbit

Immunogen Synthetic peptide corresponding to the N-terminus of human GATA3.

Isotype IgG

Purity Detail Protein A & antigen affinity purified.

Recommendation Western Blot (1:500-1:2'000)

Dilutions/Conditions Immunoprecipitation (5-10 µL/mg of lysate)

Suggested dilutions/conditions may not be available for all applications.
Optimal conditions must be determined individually for each application.

Species Reactivity Human, Mouse

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)