

IBA1 recombinant monoclonal antibody (031)

The anti-IBA1 antibody is a highly specific and sensitive tool for detecting IBA1 (Ionized Calcium-Binding Adapter Molecule 1), a key protein expressed in microglia and macrophages. This antibody is optimized for use in immunohistochemistry (IHC), immunofluorescence (IF), and ELISA, making it an essential tool for studying microglial activation, neuroinflammation, and immune responses in the central nervous system.

IBA1 (Ionized Calcium-Binding Adapter Molecule 1) is a cytoplasmic protein highly expressed in microglia, the resident immune cells of the central nervous system (CNS), as well as in macrophages. It is a member of the EF-hand calcium-binding protein family and plays a crucial role in microglial activation, phagocytosis, and cell migration in response to injury or disease. IBA1 is widely used as a specific marker for microglia in studies of neuroinflammation and neurodegeneration. Its upregulation is associated with the activation of microglia during pathological conditions such as Alzheimer's disease, Parkinson's disease, multiple sclerosis, and stroke. IBA1 also participates in cytoskeletal reorganization, enabling microglia to adapt to the dynamic demands of the CNS environment.

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

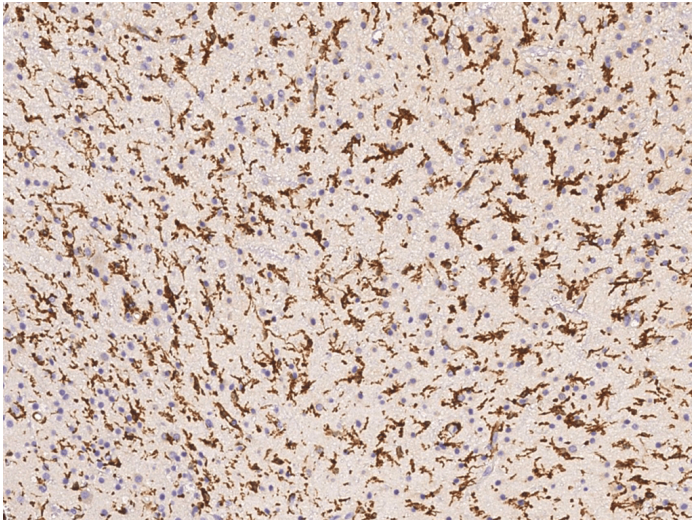
Ordering Information

[Order Online »](#)

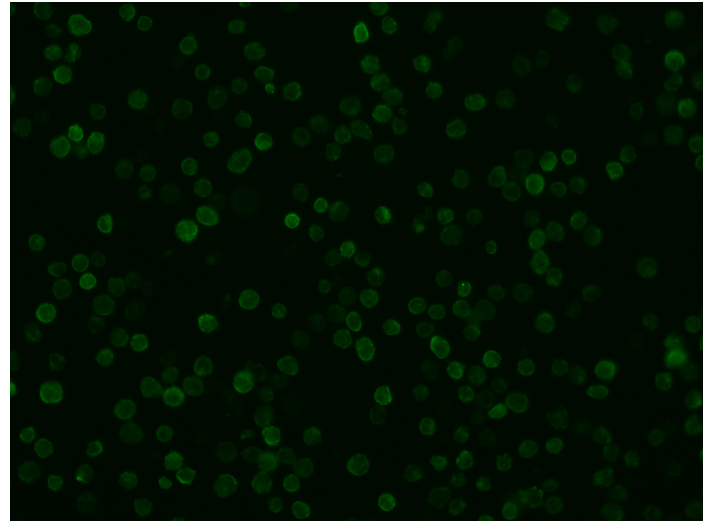
ENZ-ABS1052-0100	100µl
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Immunohistochemical staining of human Iba1 in human brain with rabbit monoclonal antibody (1:200, formalin-fixed paraffin embedded sections).



Immunofluorescence staining of Iba1 in K562 cells. Cells were fixed with 4% PFA, permeabilized with 0.1% Triton X-100 in PBS, blocked with 10% serum, and incubated with rabbit anti-human AIF1 monoclonal antibody (dilution ratio 1:60) at 4°C overnight. Then cells were stained with the Alexa Fluor®488-conjugated Goat Anti-rabbit IgG secondary antibody (green). Positive staining was localized to Cytoplasm and Cell membrane.

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 AIF1, Ionized Calcium-Binding Adapter Molecule 1, Allograft inflammatory factor 1

Application ELISA, ICC, IF, IHC (PS)

Clone 031

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

Host Rabbit

Immunogen Recombinant human IBA1 / AIF1 protein.

Isotype IgG

Purity Detail Protein A purified.

Recommendation ELISA (1:5'000-1:10'000) Immunohistochemistry (paraffin sections) (1:100-1:500)
Dilutions/Conditions Immunocytochemistry/Immunofluorescence (1:20-1:100)

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

Species Reactivity Human

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)