

CD11b polyclonal antibody

The anti-CD11b antibody detects human CD11b, an integrin alpha subunit expressed on monocytes, macrophages, granulocytes, and microglia. This antibody is optimized for use in immunohistochemistry (IHC).

CD11b, also known as integrin alpha M, is a transmembrane protein that forms part of the integrin $\alpha M\beta 2$ complex (Mac-1 or CR3) when paired with the CD18 beta subunit. It plays a critical role in immune responses, mediating leukocyte adhesion, migration, and phagocytosis. CD11b is predominantly expressed on monocytes, macrophages, neutrophils, and microglia, making it a key marker for studying innate immunity and neuroinflammation. In the central nervous system, CD11b is a critical marker for activated microglia, the resident immune cells of the brain, and is widely used to study neuroinflammatory responses in neurodegenerative diseases such as Alzheimer's and Parkinson's disease. It is also involved in pathogen recognition, tissue remodeling, and inflammation, making it a target of interest in autoimmune disorders and infectious diseases.

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

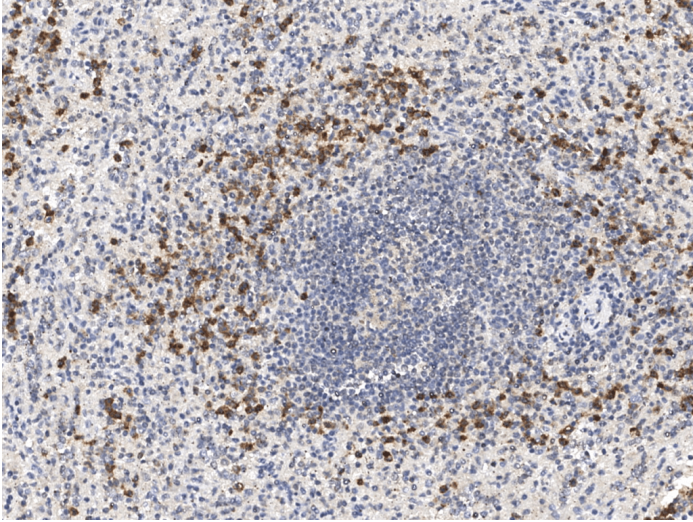
Ordering Information

[Order Online »](#)

ENZ-ABS1053-0100	100µl
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Immunohistochemical staining of human CD11b in human spleen with rabbit polyclonal antibody at 1:300 dilution, formalin-fixed paraffin embedded sections.

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	ITGAM, Integrin alpha M
Application	IHC (PS)
Formulation	Liquid. 0.2 µm filtered solution in PBS. Carrier-free. Preservative-free.
Host	Rabbit
Immunogen	Recombinant human CD11b / ITGAM / Integrin alpha M fragment.
Isotype	IgG
Purity Detail	Protein A and antigen affinity purified.
Recommendation	Immunohistochemistry (paraffin sections) (1:200-1:1'000)
Dilutions/Conditions	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

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com