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.

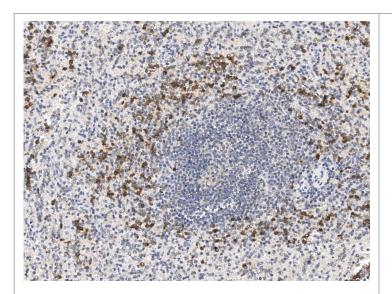
Ordering Information

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.

lgG lgG

Purity Detail Protein A and antigen affinity purified.

Recommendation Dilutions/Conditions Immunohistochemistry (paraffin sections) (1:200-1:1'000)

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

•



European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

Phone: +33 472 440 655 infofr@enzolifesciences.com Germany Phone: +49 7621 5500 526 infode@enzolifesciences.com UK & Ireland Phone (UK customers): 0845 601 1488 Phone: +44 1392 825900 infouk@enzolifesciences.com