

CXCR4 (mouse) polyclonal antibody

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

Manuals, SDS & CofA

[View Online »](#)

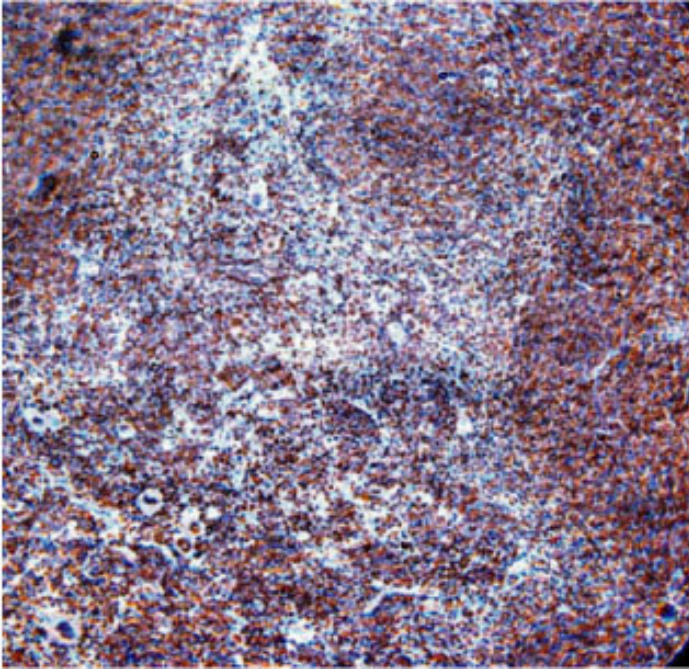


Figure: Staining of CXCR4 in mouse spleen paraffin section using PAb to CXCR4 (mouse) (Prod. No. [ALX-210-819](#)).^{”””},CXCR4 (mouse) polyclonal antibody Immunohistochemistry,CXCR4 (mouse) polyclonal antibody Immunohistochemistry,yes ALX-210-838,210-838-n2.gif,<http://static.enzolifesciences.com/fileadmin/files/image/210-838-n2.gif>,**Figure: (B)** Detection of caspase-9 processing during apoptosis. The antiserum detects procaspase-9 (46kDa) and the intermediate cleavage products of 37kDa and 35 kDa. No cross-reactivity with other caspases is observed.**Method:** Jurkat cells were treated with 2µM staurosporine (Prod. No. [ALX-380-014](#)). After 4 hours cell lysates were prepared and separated by SDS-Page (2×10^6 cells/lane) under reducing conditions. Proteins were immunoblotted with rabbit anti-caspase-9 (1:1'000). Following incubation with peroxidase-conjugated secondary antibodies caspase-9 processing was detected by enhanced chemoluminescent staining. The closed arrow indicates procaspase-9 (46kDa)

Handling & Storage

Handling	For maximum product recovery after thawing, centrifuge the vial before opening the cap. Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	CXC chemokine receptor 4, CD184
Application	ICC, IHC (PS), WB
Formulation	Liquid. In PBS containing 1mg/ml BSA and 0.1% sodium azide.
Host	Goat
Immunogen	Synthetic peptide corresponding to aa 14-40 (Y14SEEVGSGDYDSNKEPCFRDENVHFNR40) of the N-terminal extracellular domain of mouse CXCR4 (CXC chemokine receptor 4).
Purity Detail	Epitope-affinity purified IgG.
Quality Control	Immunocytochemistry on mouse spleen cells.
Recommendation Dilutions/Conditions	Immunocytochemistry (1:300)Immunohistochemistry (paraffin sections, 1:200)Western Blot (1:1,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Mouse
UniProt ID	O08565
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee .

Last modified: May 29, 2024



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