

# CXCR4 (mouse) polyclonal antibody

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

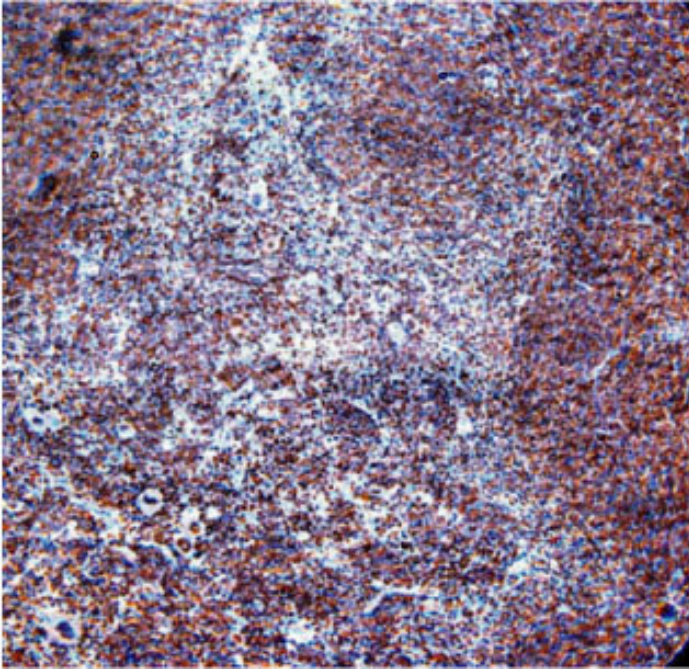
## Ordering Information

[Order Online »](#)

ALX-210-819-C200	200µg
------------------	-------

## Manuals, SDS & CofA

[View Online »](#)



**Figure:** Staining of CXCR4 in mouse spleen paraffin section using PAb to CXCR4 (mouse) (Prod. No. [ALX-210-819](#)).<sup>”””</sup>,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<sup>6</sup> 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 <a href="#">Worry-Free Guarantee</a> .



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)