

CCR4 (mouse) polyclonal antibody

Mouse CCR4 is the homolog of human CCR4 receptor (85% identical), binding macrophage inflammatory protein-1 α (MIP-1 α) and RANTES. Mouse CCR4 transcripts are expressed in the thymus and in T-cell lines.

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

Citations: 7

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-210-621-C200	200 μ g
------------------	-------------

Manuals, SDS & CofA

[View Online »](#)

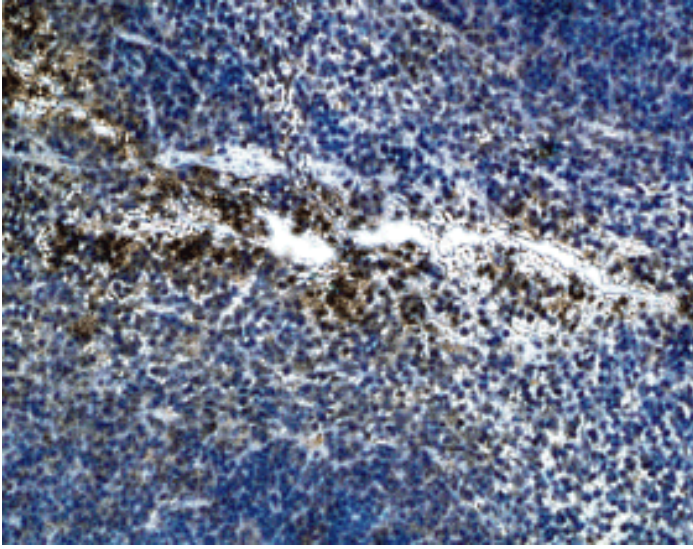


Figure: Immunohistochemistry on paraffin section of mouse spleen with CCR4 (mouse), pAb (Prod. No. ALX-210-621).

Handling & Storage

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

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	C-C chemokine receptor 4, CD194
Application	ELISA, Flow Cytometry, ICC, IHC (PS), WB
Crossreactivity	Does not cross-react with human CCR4.
Formulation	Liquid. In PBS containing 1mg/ml BSA and 0.1% sodium azide.
Host	Goat
Immunogen	Synthetic peptide corresponding to aa 8-26 (D8TTQDETVYNSYYFYESMP26C) of the N-terminal extracellular domain of mouse CCR4 (C-C chemokine receptor 4).
Purity Detail	Epitope-affinity purified F(ab`) ² fragment.
Quality Control	Immunocytochemistry on mouse spleen cells.
Recommendation Dilutions/Conditions	ELISA (1:100,000)Flow Cytometry (1-3µl per 2×10 ⁵ cells per 50µl)Immunocytochemistry (1:500)Immunohistochemistry (1:500)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	P51680

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

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