

# CCR6 polyclonal antibody

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

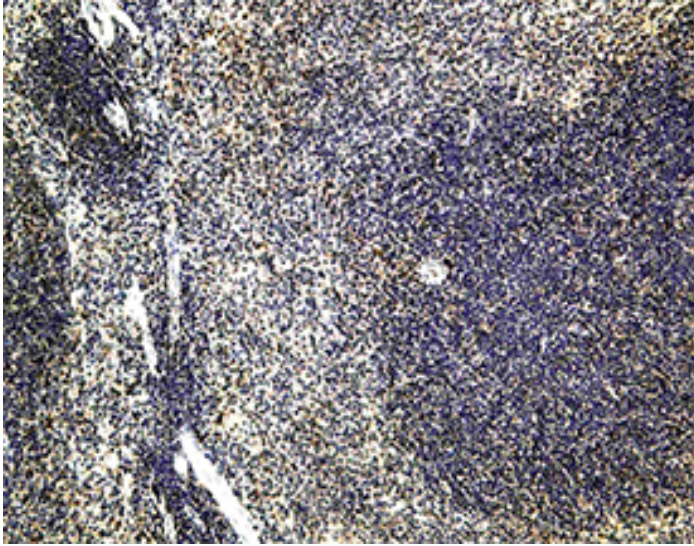
## Ordering Information

[Order Online »](#)

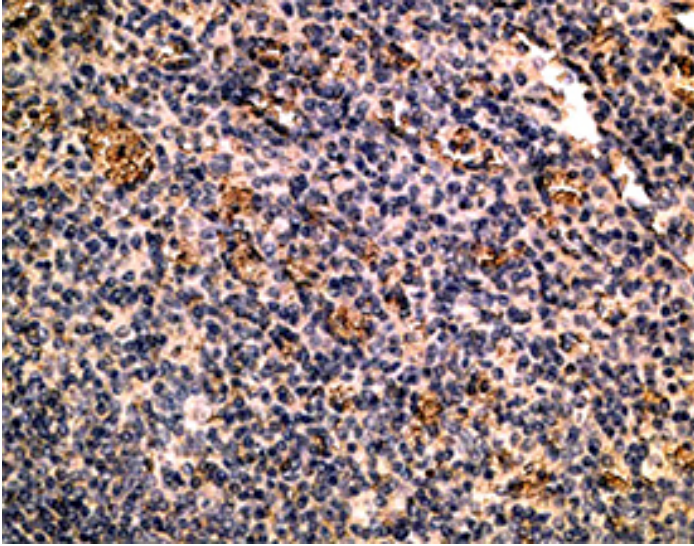
ALX-210-843-C200	200µg
------------------	-------

## Manuals, SDS & CofA

[View Online »](#)



Immunohistochemistry with CCR6, pAb (Prod. No. ALX-210-843) to human CCR6 on paraffin section of mouse spleen.



Immunohistochemistry with CCR6, pAb (Prod. No. ALX-210-843) to human CCR6 on paraffin section of human spleen.

## 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	C-C chemokine receptor 6, CD196
Application	ELISA, Flow Cytometry, ICC, IHC (PS)
Formulation	Liquid. In PBS containing 1mg/ml BSA and 0.1% sodium azide.
Host	Goat
Immunogen	Synthetic peptide corresponding to aa 18-46 (Y18FVSVNTSYYSVDSEMLLCSLQEVQRQFSR46) of human CCR6 (C-C chemokine receptor 6). Peptide sequence is <50% identical to other human chemokine receptors in this region.
Purity Detail	Epitope-affinity purified IgG.
Quality Control	Immunocytochemistry on human spleen or peripheral blood.
Recommendation Dilutions/Conditions	ELISA (1:75,000)Flow Cytometry (1-3µl per 106 cells/50-100µl)Immunocytochemistry (1:300)Immunohistochemistry (paraffin sections, 1:250)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human, Mouse
UniProt ID	P51681

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



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-be@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
fr@enzolifesciences.com](mailto:info-fr@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)