## CCR6 polyclonal antibody

This antibody is covered by our Worry-Free Guarantee.

**Ordering Information** 

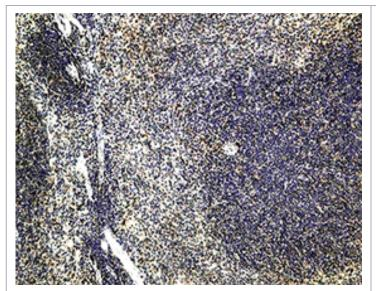
Order Online »

ALX-210-843-C200

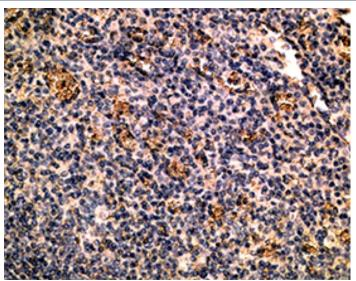
200µg

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

(Y18FVSVNTSYYSVDSEMLLCSLQEVRQFSR46) 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

Last modified: May 29, 2024



info-

eu@enzolifesciences.com