# CCR2 (human) polyclonal antibody

This antibody is covered by our Worry-Free Guarantee.

**Ordering Information** 

Order Online »

ALX-210-821-C200

200µg

Manuals, SDS & CofA

**View Online »** 

# **Handling & Storage**

Handling For maximum product recovery after thawing, centrifuge the vial before opening the cap.

Avoid freeze/thaw cycles.

-20°C **Long Term Storage** 

**Shipping** Blue Ice

## Regulatory Status RUO - Research Use Only

### **Product Details**

**Alternative Name** C-C chemokine receptor 2, CD192

**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 17-

> 39 (G17EEVTTFFDYDYGAPCHKFDVKQ39) of the Nterminal extracellular domain of human CCR2 (C-C

chemokine receptor 2).

**Purity Detail** Epitope-affinity purified IgG.

**Recommendation Dilutions/Conditions** Immunocytochemistry (1:400)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** Human

**UniProt ID** P41597

### **Worry-free Guarantee**

This antibody is covered by our Worry-Free Guarantee



ENZO LIFE SCIENCES, Phone: 800.942.0430 info-

ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 usa@enzolifesciences.com eu@enzolifesciences.com

European Sales Office

Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France Phone: +33 472 440 655 fr@enzolifesciences.com

Phone: +49 7621 5500 526 de@enzolifesciences.com

UK & Ireland Phone (UK customers): 0845 601 1488 Phone: +44 1392 825900 uk@enzolifesciences.com