CCR8 (mouse) polyclonal antibody

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

ALX-210-845-C200

200µg

Manuals, SDS & CofA

View Online »

Handling & Storage

Handling 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 8, CD198

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 7-33

(P7NVTMTDYYPDFFTAPCDAEFLLRGSM33) of mouse

CCR8 (C-C chemokine receptor 8).

Purity Detail Epitope-affinity purified IgG.

Quality Control Immunocytochemistry on mouse spleen cells.

Recommendation Dilutions/Conditions Immunocytochemistry (1:500)Immunohistochemistry

(paraffin sections (1:400))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 P56484

Worry-free Guarantee

This antibody is covered by our Worry-Free Guarantee



ENZO LIFE SCIENCES, INC. Phone: 800.942.0430 infousa@enzolifesciences.com European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

Phone: +33 472 440 655
infofr@enzolifesciences.com

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

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