

CCR7 polyclonal antibody

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

Ordering Information

[Order Online »](#)

ALX-210-844-C100	100µg
------------------	-------

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. 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 7, CD197
Application	Flow Cytometry, 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 2-22 (D2PGKPRKNVLVALLVIFQVC22) of mouse CCR7 (C-C chemokine receptor 7).
Purity Detail	Epitope-affinity purified IgG.
Quality Control	Immunohistochemistry on mouse spleen.
Recommendation	Flow Cytometry (1:10)Immunocytochemistry (1:500)Immunohistochemistry (paraffin
Dilutions/Conditions	sections, 1:250)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, Mouse
Technical Info / Product Notes	Sequence comparison: Mouse CCR7: DPGKPRKNVLVALLVIFQVC Human CCR7: DLGHPMKSVLVALLVIFQVC
UniProt ID	P47774



ENZO LIFE SCIENCES,
INC.
Phone: +1 942.042.0430
info-
usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
Phone: +41 61 926 8989
info-
eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 2 460 0000
info-
be@enzolifesciences.com

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

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

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

This antibody is covered by our **Worry-Free Guarantee**.