

CCR8 (human) (NT) polyclonal antibody

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

Citations: 5

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-210-762-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. 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 8, CD198
Application	ELISA, Flow Cytometry, IHC, WB
Formulation	Affinity purified antibody in 10mM KHPO ₄ , 140mM NaCl, 1mg/ml BSA and 0.1% sodium azide.
Host	Goat
Immunogen	Synthetic peptide corresponding to aa 12-36 (V12TDYYYPDIFSSPCDAELIQTNGKL36) of the N-terminal extracellular domain of human CCR8 (C-C chemokine receptor 8).
Purity Detail	Epitope-affinity purified IgG.
Recommendation Dilutions/Conditions	ELISA (1:50,000)Flow Cytometry (2-6µl per 10 ⁶ cells/50-100µl)Immunohistochemistry (1:125)Western blot (1:500)Optimal conditions must be determined individually for each application.
Species Reactivity	Human
UniProt ID	P51685
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee



Last modified: May 29, 2024

ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430

info@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989

info@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420

info@enzolifesciences.com

France
Phone: +33 472 440 655

info@enzolifesciences.com

Germany
Phone: +49 7621 5500 526

info@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488

Phone: +44 1392 825900
info@enzolifesciences.com