

# CCR8 (human) (NT) polyclonal antibody

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

Citations: 5

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-210-7620	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 <sub>4</sub> , 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 <sup>6</sup> 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 <a href="#">Worry-Free Guarantee</a>



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430

[info@enzolifesciences.com](mailto:info@enzolifesciences.com)  
Last modified: May 29, 2024

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

[info@enzolifesciences.com](mailto:info@enzolifesciences.com)  
[eu@enzolifesciences.com](mailto:eu@enzolifesciences.com)

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

[info@enzolifesciences.com](mailto:info@enzolifesciences.com)  
[be@enzolifesciences.com](mailto:be@enzolifesciences.com)

France  
Phone: +33 472 440 655

[info@enzolifesciences.com](mailto:info@enzolifesciences.com)  
[fr@enzolifesciences.com](mailto:fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526

[info@enzolifesciences.com](mailto:info@enzolifesciences.com)  
[de@enzolifesciences.com](mailto:de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900

[info@enzolifesciences.com](mailto:info@enzolifesciences.com)  
[uk@enzolifesciences.com](mailto:uk@enzolifesciences.com)