

CCL22 (human) monoclonal antibody (349/6)

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

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	Preservative may be added for storage at 4°C.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	MDC
Application	ELISA, WB
Clone	349/6
Formulation	Liquid. 0.2µm-filtered solution in PBS, pH 7.4. Contains no preservatives.
Host	Rat
Immunogen	Recombinant human CCL22 (macrophage-derived chemokine; MDC).
Isotype	IgG
Purity Detail	Protein G-affinity purified.
Recommendation Dilutions/Conditions	Western Blot (1-2µg/ml)ELISA (capture, 20µg/ml; use microtiter plates coated with this antibody in 0.05M carbonate/bicarbonate buffer, pH 9.6; substrate: 0.26mg/ml tetramethyl benzydine and 0.01% hydrogen peroxide stabilized in 0.05M citrate buffer, pH 3.8)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human
UniProt ID	O00626

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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