CCL22 (human) monoclonal antibody (349/6)

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

ALX-804-469-C250

250µg

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 000626

Last modified: May 29, 2024



info-

eu@enzolifesciences.com