CCL1 (human) monoclonal antibody (8P16-N4-L21)

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

ALX-804-489-C100

100µg

Manuals, SDS & CofA

View Online »

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name I-309

Application IHC (FS)

Clone 8P16-N4-L21

Formulation Liquid. In PBS.

Host Mouse

Immunogen Recombinant human CCL1 (I-309).

lgM lgM

Purity Detail DEAE purified.

Recommendation Dilutions/Conditions Immunohistochemistry (frozen sections, 5µg/ml)Suggested

dilutions/conditions may not be available for all

applications. Optimal conditions must be determined

individually for each application.

Species Reactivity Human

UniProt ID P22362

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

