CCR3 Recombinant monoclonal antibody (5E8-G9-B4)(Rabbit IgGk)

CCR3 is a chemokine receptor highly expressed of eosinophils, as well as Th2-Tcells, basophils, platelets, mast-cells, mononuclear phagocytes, airway epithelial cells, and hematopoietic progenitor cells (CD34+).

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

ENZ-ABS616-0200

200µg

Manuals, SDS & CofA

View Online »

Handling & Storage

Use/Stability Stable at 4°C for up to 3 months. For longer storage, aliquot and store at -20°C.

Handling Avoid freeze/thaw cycles.

Short Term Storage +4°C

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Flow Cytometry

Clone 5E8-G9-B4

Formulation Liquid. In PBS containing 0.02% Proclin 300.

Host Rabbit

Isotype IgGk

Purity Detail Protein A affinity purified.

Species Reactivity Human

Technical Info / Product NotesThis chimeric rabbit antibody was made using the variable

domain sequences of the original murine IgG2b format, for improved compatibility with existing reagents, assays and

techniques.

UniProt ID P51677

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

