

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

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](#).

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 Notes

This 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](#)



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
Last modified: **May 29, 2024**
usa@enzolifesciences.com

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info@enzolifesciences.com

France
Phone: +33 472 440 655
info@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info@enzolifesciences.com

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