

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](#).

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 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
[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)