

CD16/CD32

Recombinant monoclonal antibody (2.4G2) (Rat IgG2bκ)

This antibody is covered by our [Worry-Free Guarantee](#).

Ordering Information

[Order Online »](#)

ENZ-ABS411-0200	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

Alternative Name	Fcy III/II Receptor, Low affinity immunoglobulin γ Fc region receptor III, IgG Fc receptor III, Fc-γ RIII, FcRIII, Low affinity immunoglobulin γ Fc region receptor II, Fcy receptor IIB, Fc-γ RII, Fc-γ-RIIB, FcRIIIgG Fc receptor IIβ, Lymphocyte antigen 17, Ly-17
Application	Flow Cytometry, FUNC, IF, IHC, IP
Clone	2.4G2
Formulation	Liquid. In PBS containing 0.02% Proclin 300.
Host	Rat
Immunogen	Mouse BALB/c Macrophage J774 Cell Line.
Isotype	IgG2bk
Purity Detail	Protein A affinity purified.
Species Reactivity	Mouse
Specificity	Recognizes an epitope on the extracellular domains of the mouse FcγIII and FcγII, thereby inhibiting their ability to bind antibodies via their Fc region. Also binds the FcγI receptor (CD64) via its Fc domain.
UniProt ID	P08508 (CD16), P08101 (CD32)

Worry-free Guarantee

This antibody is covered by our [Worry-Free Guarantee](#)

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

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

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

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