

CD35 Recombinant monoclonal antibody (8C12) (Rat IgG2ak)

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

Ordering Information

[Order Online »](#)

ENZ-ABS436-0200	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	CR1, C3BR, C4BR, KN, Complement receptor 1
Application	Flow Cytometry, IHC (FS), IP, WB
Clone	8C12
Formulation	Liquid. In PBS containing 0.02% Proclin 300.
Host	Rat
Immunogen	Purified murine CR1 protein.
Isotype	IgG2ak
Purity Detail	Protein A affinity purified.
Species Reactivity	Mouse
Specificity	Recognizes an epitope on CD35 (CR1, 190 kDa), but not on CD21 (CR2, 145-150 kDa), which are alternative splice variants of primary Cr2 gene transcripts.
UniProt ID	P17927

Worry-free Guarantee



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

This antibody is covered by our **Worry-Free Guarantee**.