

F4/80 Recombinant monoclonal antibody (Recombinant Cl:A3-1) (Rabbit IgGκ)

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

Ordering Information

[Order Online »](#)

ENZ-ABS638-0200	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	EGF-like module, Cell surface glycoprotein F4/80, EGF-like module receptor 1, EMR1 receptor, Em Gpf4, TM7, Ly-72
Application	Flow Cytometry, IHC
Clone	Recombinant Cl:A3-1
Formulation	Liquid. In PBS containing 0.02% Proclin 300.
Host	Rabbit
Immunogen	Thioglycollate stimulated peritoneal macrophages of mouse origin.
Isotype	IgGk
Purity Detail	Protein A affinity purified.
Species Reactivity	Mouse
Technical Info / Product Notes	This chimeric rabbit antibody was made using the variable domain sequences of the original Rat IgG2b format, for improved compatibility with existing reagents, assays and techniques.

UniProt ID Q61549



Worry-free Guarantee

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

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +32 3 466 0420
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
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