F4/80 Recombinant monoclonal antibody (Recombinant CI: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 IgGĸ

Purity Detail Protein A affinity purified.

Species Reactivity Mouse

Technical Info / Product NotesThis 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

Last modified: May 29, 2024



info-

eu@enzolifesciences.com