

CD40L (mouse) monoclonal antibody (MR1)

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

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-804-331-C100	100µg
ALX-804-331-C500	500µg

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name CD154, gp39, TNFSF 5

Application FUNC

Application Notes Functional Application: blocks T cell co-stimulation *in vivo*, promoting allograft acceptance and donor-specific tolerance to secondary skin grafts.

Clone MR1

Formulation Liquid. In PBS.

Host Hamster

Immunogen Mouse T-cells (Th1).

Isotype IgG1

Purity Detail Protein A-affinity purified.

Species Reactivity Mouse

UniProt ID P27548

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



ENZO LIFE SCIENCES,
Last modified: **May 29, 2024**
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-fr@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