OX40L Recombinant monoclonal antibody (R4930) (Human IgG1ĸ)

Human OX40L, the ligand of OX40, stimulatory receptor that is expressed on activated T cells, natural killer (NK) cells and natural killer T (NKT) cells.

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

ENZ-ABS556-0200

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 TNFSF 4, Tumor necrosis factor superfamily member 4,

OX40 ligand, TAX transcriptionally-activated glycoprotein

1, TXGP1, gp34, CD252

Application Flow Cytometry, FUNC

Clone R4930 (Oxelumab)

Formulation Liquid. In PBS containing 0.02% Proclin 300.

Host Human

Immunogen Human OX40L

lgG1κ

Purity Detail Protein A affinity purified.

Species Reactivity Human

UniProt ID P23510

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 infousa@enzolifesciences.com

ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
infoeu@enzolifesciences.com

European Sales Office

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

Phone: +33 472 440 655 infofr@enzolifesciences.com

Phone: +49 7621 5500 526 infode@enzolifesciences.com UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com