CD40L Recombinant monoclonal antibody (hu5c8) (Human IgG1ĸ)

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

ENZ-ABS450-0200

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 CD40 ligand, Tumor necrosis factor ligand superfamily

member 5, TNFSF 5, CD154, tumour necrosis factor

related activation protein, TRAP

Application Flow Cytometry, FUNC, WB

Clone hu5c8 (Ruplizumab)

Formulation Liquid. In PBS containing 0.02% Proclin 300.

Host Human

Immunogen Human CD40L

IgG1ĸ

Purity Detail Protein A affinity purified.

Species Reactivity Human

Specificity The antibody binds to CD40L with an IC50 of 0.845 μg/ml.

UniProt ID P29965

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

.

Last modified: May 29, 2024



info-

eu@enzolifesciences.com