CD40 Recombinant monoclonal antibody (G28.5) (Mouse IgG1κ)

48 kD Type I glycoprotein, a member of the TNFR superfamily primarily expressed on normal or malignant B cells or B cell lines, macrophages, follicular dendritic cells, endothelial cells, fibroblasts, and at low levels on plasma cells.

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

ENZ-ABS448-0200

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 TNFRSF 5, Tumor necrosis factor receptor superfamily

member 5, B-cell surface antigen CD40, Bp50, CD40L

receptor CDw40

Application Flow Cytometry, IHC, IP

Application Notes Functional Application: Activation

Clone G28.5

Formulation Liquid. In PBS containing 0.02% Proclin 300.

Host Mouse

Immunogen Human erythrocyte-rosette-negative tonsillar lymphocytes.

lgG1κ

Purity Detail Protein A affinity purified.

Species Reactivity Human

UniProt ID P25942

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

.

Last modified: May 29, 2024



info-

eu@enzolifesciences.com