

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.
Isotype	IgG1κ
Purity Detail	Protein A affinity purified.
Species Reactivity	Human
UniProt ID	P25942

Worry-free Guarantee



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)

This antibody is covered by our **Worry-Free Guarantee**.