

CD40 (human) monoclonal antibody (G28-5)

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

Citations: 1

[View Online »](#)

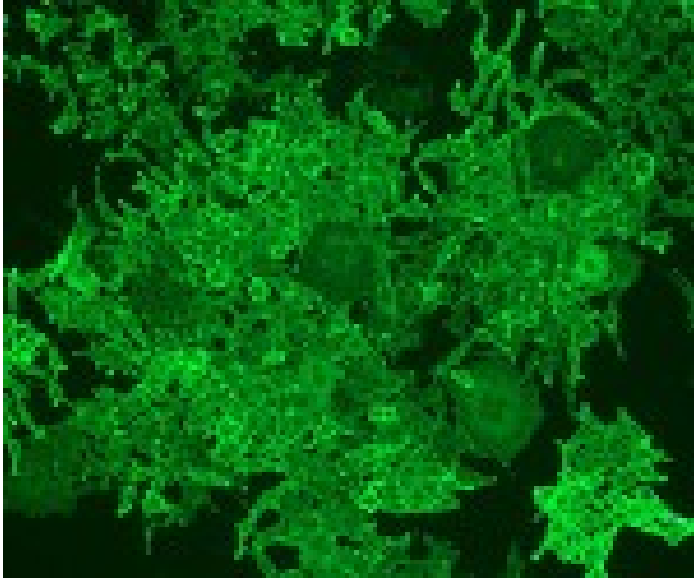
Ordering Information

[Order Online »](#)

ENZ-ABS130-0100	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Immunohistochemistry analysis of CD40-transfected L cells stained with CD40 (human) mAb, clone G28-5.

Handling & Storage

Short Term Storage -20°C

Long Term Storage -80°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Tumor necrosis factor receptor superfamily member 5, TNFRSF 5, B-cell surface antigen CD40, Bp50, CD40L receptor, CDw40

Application Flow Cytometry, FUNC, IHC (PS)

Application Notes Functional Application: cell activation.

Clone G28-5

Formulation Liquid. In TRIS-NaCl, pH 8.

Host Mouse

Immunogen Human erythrocyte-rosette-negative tonsillar lymphocyte.

Isotype IgG1k

Purity Detail Purified using QMA Hyper D ion exchange chromatography

Species Reactivity Human

UniProt ID P25942

Worry-free Guarantee This antibody is covered by our [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)