

# CD40 Recombinant monoclonal antibody (G28.5) (Rabbit IgGκ)

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-ABS447-0200 | 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                           | Rabbit   |
| Immunogen                      | Human erythrocyte-rosette-negative tonsillar lymphocytes.  |
| Isotype                        | IgGκ   |
| Purity Detail                  | Protein A affinity purified.   |
| Species Reactivity             | Human  |
| Technical Info / Product Notes | This chimeric rabbit antibody was made using the variable domain sequences of the original murine IgG1 format, for improved compatibility with existing reagents, assays and techniques. |
| 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)