

CD40 (mouse) monoclonal antibody (FGK45)

Widely used stimulatory MAb to CD40. Shown to indirectly activate natural killer (NK) cells, producing significant antitumor and antimetastatic effects. Effective in boosting immune responses against infectious agents and can potentially be used to treat chronic autoimmune inflammatory processes.

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

Citations: 29

[View Online »](#)

Ordering Information

[Order Online »](#)

| | |
|------------------|-------|
| ALX-805-046-C100 | 100µg |
| ALX-805-046-C500 | 500µg |

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

| | |
|--------------------|---------------------------|
| 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 |
| Application | ELISA, Flow Cytometry, FUNC, WB |
| Application Notes | Functional Application: activates B and NK cells <i>in vivo</i> and <i>in vitro</i> . |
| Clone | FGK45 |
| Formulation | Liquid. In PBS containing 0.02% sodium azide. |
| Host | Rat |
| Immunogen | Recombinant mouse CD40 fusion protein. |
| Isotype | IgG2a |
| Purity Detail | Protein G-affinity purified. |
| Source | Purified from hybridoma tissue culture supernatant. |
| Species Reactivity | Mouse |
| UniProt ID | P27512 |
| 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)