

CD40 (mouse) monoclonal antibody (FGK45) (preservative free)

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: 21

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-805-046PF-C100	100µg
ALX-805-046PF-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 sterile PBS.
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-
fr@enzolifesciences.com](mailto:info-fr@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)