

# CD40 (mouse) monoclonal antibody (FGK45) (biotin conjugate)

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: 11

[View Online »](#)

## Ordering Information

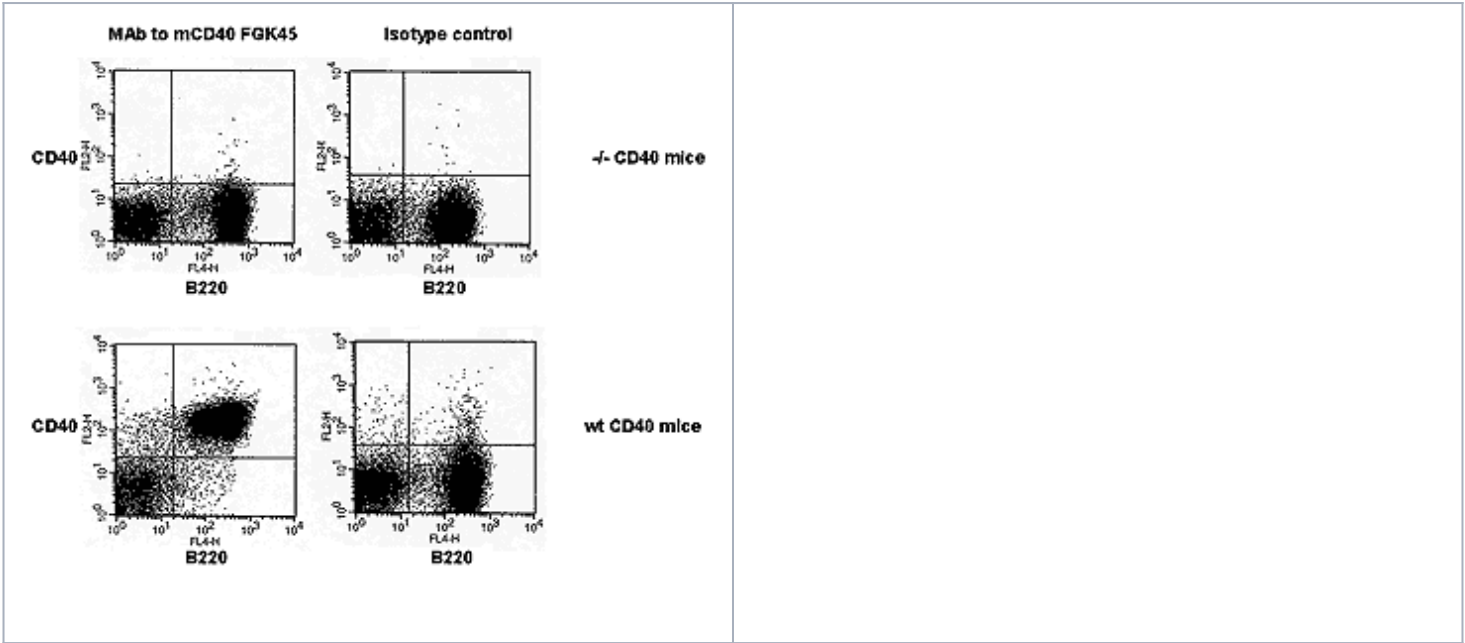
[Order Online »](#)

ALX-805-046B-C050

Biotin - 50µg

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Handling	Do not freeze. Protect from light.
Short Term Storage	+4°C
Long Term Storage	+4°C
Shipping	Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

Alternative Name	TNFRSF 5
Application	ELISA, Flow Cytometry
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 <a href="#">Worry-Free Guarantee</a> .



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)