

CD40-mulg Fusion Protein

Ordering Information

[Order Online »](#)

ANC-504-050	R-PE - 50 tests
ANC-504-060	APC - 50 tests

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name TNFRSF 5

Application Notes FUNC (Blocks binding of anti-human CD40 MAbs to Raji human tumor cells).

Source Produced in CHO cells. Soluble molecule consisting of the extracellular domain (173 aa) of human CD40 fused to the Fc portion of mouse IgG2a (232 aa).

Technical Info / Product Notes For the Original Manufacturer's Data Sheet please click on the corresponding product number:

[ANC-504-020](#)

[ANC-504-030](#)

[ANC-504-050](#)

[ANC-504-060](#)

[ANC-504-820](#)

UniProt ID P25942



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com