

CD28 (stimulating) (human), mAb (ANC28.1/5D10)

Ordering Information

[Order Online »](#)

ANC-177-820	Preservative Free - 100µg
-------------	---------------------------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application ELISA, Flow Cytometry, FUNC

Application Notes Functional Application (monomeric and cross linked ANC28.1 stimulates expression of IL-2 from CD28⁺ cells).

Clone ANC28.1/5D10

Host Mouse

Immunogen Human CD28 Ig fusion protein and HPB-MLT human tumor cells.

Isotype IgG1

Species Reactivity Human

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

[ANC-177-020](#)

[ANC-177-030](#)

[ANC-177-040](#)

[ANC-177-050](#)

[ANC-177-520](#)

[ANC-177-820](#)

[For functional application note please click here.](#)



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