

CD32 (human), mAb (7.3)

Ordering Information

[Order Online »](#)

ANC-181-580

Fab - 100µg

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 FcγRII

Application Flow Cytometry

Application Notes Functional Application (blocks immune complex binding, as does the F(ab')₂ fragment). The F(ab')₂ fragment blocks non-antigen-specific binding of immunoglobulins to the FcγRII and FcγRI receptors *in vitro*. It is therefore a useful tool to reduce strong background staining of Fc-containing MAbs in FC and IHC. Also useful when recombinant Fc:fusion proteins are employed as detection tools for the detection of the cognate ligands.

Clone 7.3

Host Mouse

Immunogen Human K562 tumor cells and L cells transfected with human CD32 (FcγRII).

Isotype IgG1

Species Reactivity Human

Specificity Recognizes an epitope in domain 2 of all CD32 isoforms.

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

[ANC-181-020](#)

[ANC-181-030](#)

[ANC-181-050](#)

[ANC-181-520](#)

[ANC-181-580](#)

[ANC-181-820](#)



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