

IgM (human), mAb (UCHB1)

Ordering Information

[Order Online »](#)

| | |
|--------------------|---------------------------|
| ANC-141-030 | Biotin - 100µg |
| ANC-141-020 | Purified - 100µg |
| ANC-141-820 | Preservative Free - 100µg |
| ANC-141-050 | R-PE - 120 tests |
| ANC-141-040 | FITC - 120 tests |

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

Clone UCHB1

Host Mouse

Immunogen Prolymphocytic leukemic B 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-141-020](#)

[ANC-141-030](#)

[ANC-141-040](#)

[ANC-141-050](#)

[ANC-141-820](#)



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