

CD134L (human), mAb (ANC10G1)

Ordering Information

[Order Online »](#)

ANC-400-030	Biotin - 100µg
ANC-400-020	Purified - 100µg
ANC-400-820	Preservative Free - 100µg
ANC-400-050	R-PE - 120 tests
ANC-400-040	FITC - 120 tests
ANC-400-060	APC - 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

Alternative Name OX40L, TNFSF 4, CD252

Application ELISA, Flow Cytometry, FUNC

Application Notes Functional Application (blocks binding of recombinant CD134 (OX40))

Clone ANC10G1

Host Mouse

Immunogen HUVEC (human umbilical vein endothelial cells) and recombinant soluble human CD134L (OX40L).

Isotype IgG1

Species Reactivity Human

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

[ANC-400-020](#)

[ANC-400-030](#)

[ANC-400-040](#)

[ANC-400-050](#)

[ANC-400-060](#)

[ANC-400-820](#)



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