

# CD134L (human), mAb (ANC10G1)

## Ordering Information

[Order Online »](#)

ANC-400-050	R-PE - 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](mailto:info-usa@enzolifesciences.com)

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

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

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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