

CD202b recombinant monoclonal antibody (116)

This antibody is covered by our [Worry-Free Guarantee](#).

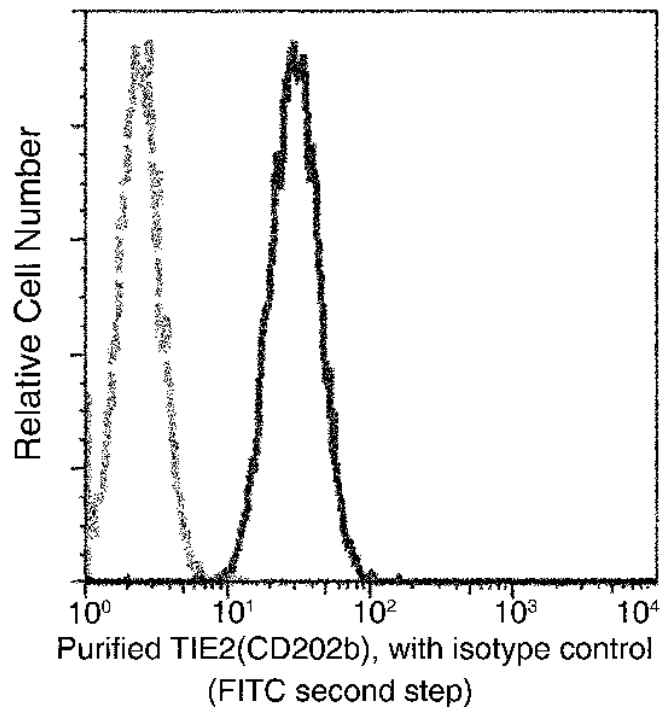
Ordering Information

[Order Online »](#)

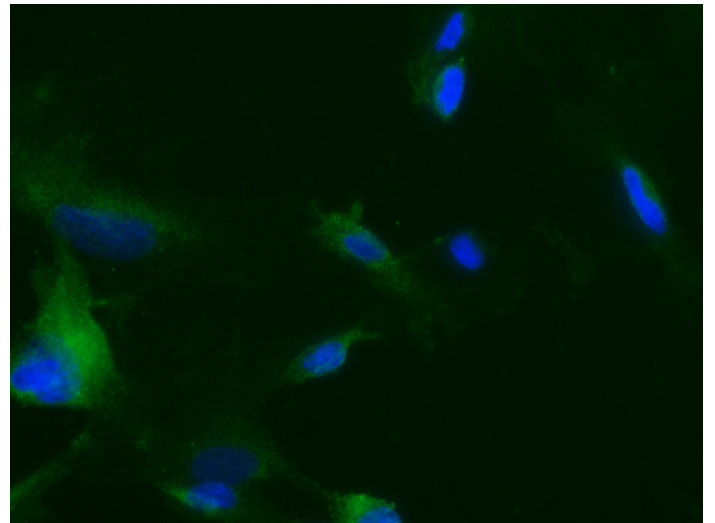
ENZ-ABS958-0100	100µl
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Flow cytometric analysis of Human TIE2(CD202b) expression on HUVEC cells. Cells were stained with purified anti-Human TIE2(CD202b), then a FITC-conjugated second step antibody. The histogram were derived from gated events with the forward and side light-scatter characteristics of intact cells.



Immunofluorescence staining of Human Tie2 in HUVEC cells. Cells were fixed with 4% PFA, permeabilized with 0.3% Triton X-100 in PBS, blocked with 10% serum, and incubated with rabbit anti-Human Tie2 monoclonal antibody (1:60) at 37 °C for 1 hour. Then cells were stained with the Alexa Fluor® 488-conjugated goat Anti-rabbit IgG secondary antibody (green) and counterstained with DAPI (blue). Positive staining was localized to cell membrane and cytoplasm.

Handling & Storage

Handling	Protected from light. Avoid freeze-thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Tie2, TEK, Angiopoietin-1 receptor
Application	ELISA, Flow Cytometry, ICC, IF
Clone	116
Formulation	Liquid. 0.2 µm filtered solution in PBS. Preservative free.
Host	Rabbit
Immunogen	Recombinant human Tie2 / CD202b / TEK protein.
Isotype	IgG
Purity Detail	Protein A purified.
Recommendation Dilutions/Conditions	ELISA (Capture) (1:250-1:2'000)Flow Cytometry (1:25-1:100)Immunocytochemistry (1:20-1:100)Immunofluorescence (1:20-1:100)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee .



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)