

Tissue Factor polyclonal antibody

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

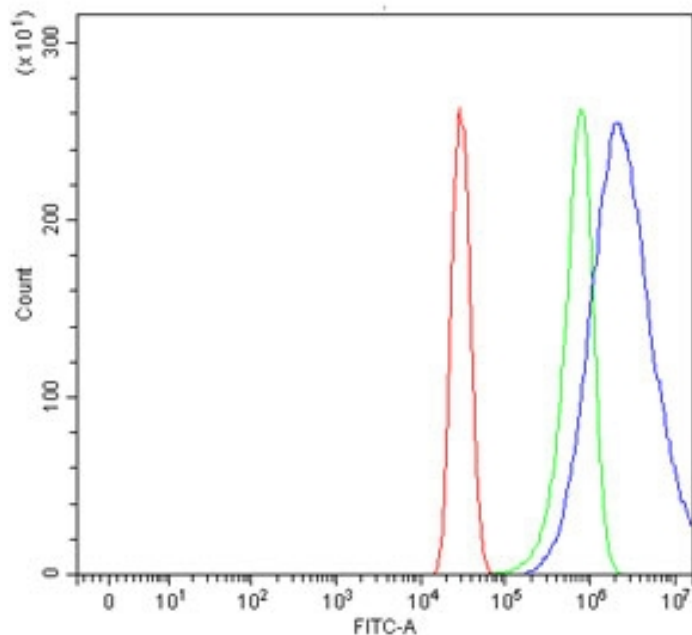
Ordering Information

[Order Online »](#)

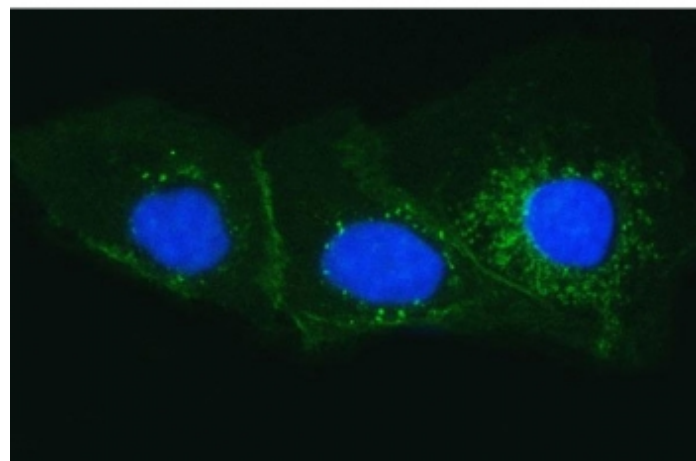
ENZ-ABS922-0100	100µg
-----------------	-------

Manuals, SDS & CofA

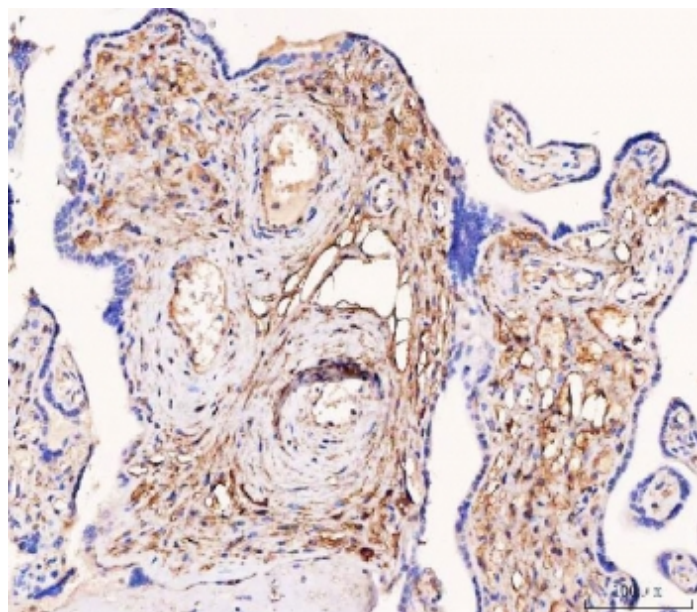
[View Online »](#)



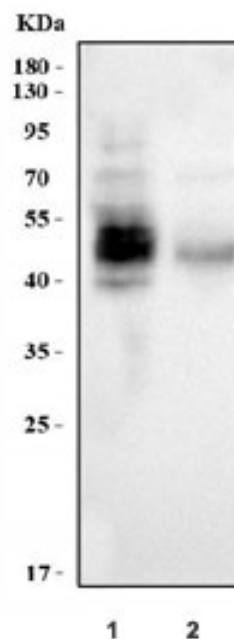
Flow cytometry testing of human A431 cells with Tissue Factor antibody at 1µg/million cells (blocked with goat sera); Red=cells alone, Green=isotype control, Blue= Tissue Factor antibody.



Immunofluorescent staining of FFPE human A431 cells with Tissue Factor antibody (green) and DAPI nuclear stain (blue). HIER: steam section in pH6 citrate buffer for 20 min.



IHC staining of FFPE human placental tissue with Tissue Factor antibody. HIER: boil tissue sections in pH8 EDTA for 20 min and allow to cool before testing.



Western blot testing of human 1) A431 and 2) HaCaT cell lysate with Tissue Factor antibody. Expected molecular weight: 33-50 kDa depending on glycosylation level.

Handling & Storage

Use/Stability	After reconstitution, the Tissue Factor antibody can be stored for up to one month at 4°C.
Handling	For long-term, aliquot and store at -20°C. Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	CD142, Coagulation factor III, Thromboplastin
Application	Flow Cytometry, IF, IHC (PS), WB
Formulation	Lyophilized from 1X PBS containing 2.5% BSA and 0.005% sodium azide.
Host	Rabbit
Immunogen	Human Tissue factor (aa 33-295).
Isotype	IgG
Purity Detail	Antigen affinity purified.
Recommendation Dilutions/Conditions	Flow cytometry (1-3µg/million cells)Immunofluorescence (5µg/ml)Immunohistochemistry (FFPE) (2-5µg/ml)Western blot (0.5-1µg/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human
UniProt ID	P13726
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)