

PRODUCT INFORMATION



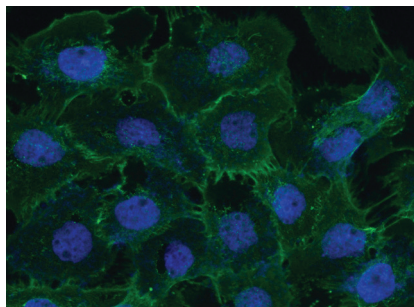
Tissue Factor/CD142 (human) Monoclonal Antibody

Item No. 43763

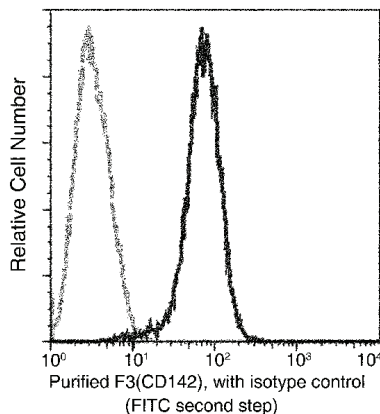
Overview and Properties

Contents:	This vial contains 100 µl of protein A purified monoclonal antibody.
Synonyms:	CD142, Cluster of Differentiation 142, Coagulation Factor III, TF, TFA, Tissue Thromboplastin
Immunogen:	Recombinant human TF
Cross Reactivity:	(+) Human
Species Reactivity:	(+) Human tissue factor; (-) HEK293 cell lysate
Form:	Liquid
Storage:	-80°C (as supplied)
Stability:	≥1 year
Storage Buffer:	0.2 µm filtered solution in PBS
Clone:	Monoclonal, Clone #05
Host:	Mouse
Isotype:	IgG2a
Applications:	ELISA, Flow Cytometry (FC), Immunocytochemistry (ICC), and Immunofluorescence (IF); the recommended starting dilution is 1:1,000-1:2,000 for ELISA, 1:25-1:100 for FC, and 1:20-1:100 for ICC and IF. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Images



Immunofluorescence staining of human tissue factor/CD142 in A431 cells. Cells were fixed with 4% PFA, permeabilized with 0.3% Triton X-100 in PBS, blocked with 10% serum, and incubated with Tissue Factor/CD142 (human) Monoclonal Antibody (1:60) at 4° overnight. Then cells were stained with the Alexa Fluor® 488-conjugated Goat Anti-mouse IgG secondary antibody (green) and counterstained with DAPI (blue). Positive staining was localized to cell membrane and cytoplasm.



Flow cytometric analysis of Tissue Factor/CD142 expression on A431 cells. Cells were stained with purified anti-Human Tissue Factor/CD142, then a FITC-conjugated second step antibody. The fluorescence histograms were derived from gated events with the forward and side light-scatter characteristics of intact cells.

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 10/28/2025

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

Tissue factor (TF), also known as CD142, is a type I transmembrane single-chain glycoprotein and the primary activator of the exogenous coagulation pathway.^{1,2} It is composed of an extracellular domain, which contains two fibronectin type III domains and interacts with factor VIIa (FVIIa), a transmembrane domain, and a cytoplasmic domain involved in signal transduction.³ TF is expressed by perivascular cells, including pericytes, adventitial fibroblasts, and vascular smooth muscle cells (VSMCs) and localizes to the cell surface.² Vascular injury exposes pre-formed TF-FVIIa complexes to its substrates, FX and FIX, initiating coagulation.^{2,3} TF levels and procoagulant activity are increased in atherosclerotic plaques isolated from patients with unstable coronary syndrome compared with those isolated from patients with stable coronary syndrome.² Overexpression of *F3*, the gene encoding TF, is positively correlated with tumor angiogenesis in advanced liver, breast, pancreatic, and lung cancers.¹ Cayman's Tissue Factor (human) Monoclonal Antibody can be used for ELISA, flow cytometry (FC), immunocytochemistry (ICC), and immunofluorescence (IF) applications.

References

1. Li, H., Yu, Y., Gao, L., *et al.* Tissue factor: A neglected role in cancer biology. *J. Thromb. Thrombolysis* **54(1)**, 97-108 (2022).
2. Grover, S.P. and Mackman, N. Tissue factor in atherosclerosis and atherothrombosis. *Atherosclerosis* **307**, 80-86 (2020).
3. Butenas, S. Tissue Factor Structure and Function. *Scientifica (Cairo)* 964862 (2012).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM