

PRODUCT INFORMATION



Tissue Factor/CD142 Extracellular Domain (human, recombinant)

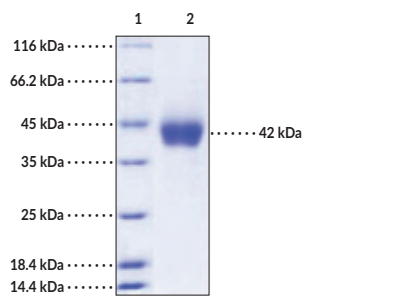
Item No. 43764

Overview and Properties

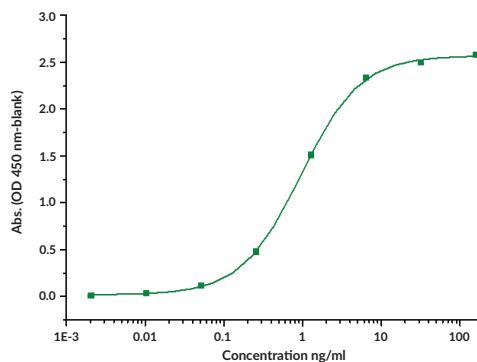
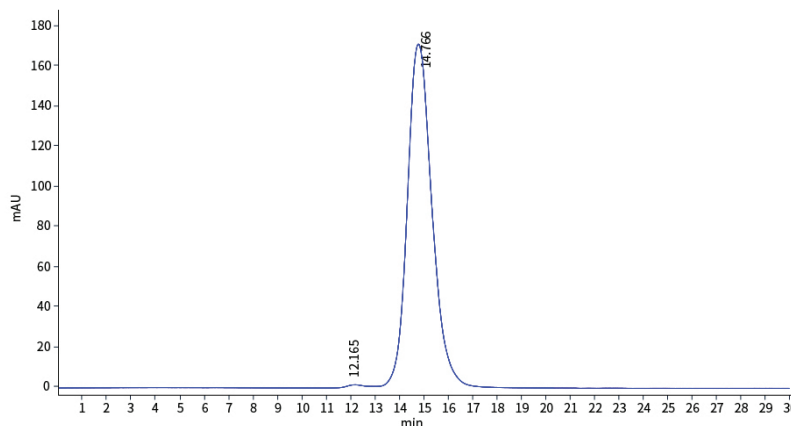
Synonyms:	CD142, Cluster of Differentiation 142, Coagulation Factor III, TF, TFA, Tissue Thromboplastin
Source:	Active recombinant human C-terminal His-tagged TF expressed in HEK293 cells
Amino Acids:	33-251
Uniprot No.:	P13726
Molecular Weight:	26.2 kDa
Storage:	-80°C (as supplied)
Stability:	≥1 year
Purity:	≥97% estimated by SDS-PAGE
Supplied in:	Lyophilized from sterile PBS, pH 7.4
Endotoxin Testing:	<1.0 EU/μg, determined by the LAL endotoxin assay
Bioactivity:	See figures for details

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Images



SDS-PAGE Analysis of Tissue Factor/CD142 Extracellular Domain. This protein has a predicted molar mass of 26.2 kDa. It has an apparent molecular weight of approximately 42 kDa by SDS-PAGE under reducing conditions.



Immobilized Tissue Factor/CD142 Extracellular Domain (human, recombinant) at 1 μg/ml (100 μl/well) binds to an anti-tissue factor/CD142 antibody EC₅₀ value of 0.9 ng/ml (routinely tested).

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, 12/23/2025

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/CD142 Extracellular Domain (human, recombinant) protein can be used for binding assays. This protein consists of 230 amino acids, has a calculated molecular weight of 26.2 kDa, and a predicted N-terminus of Ser33 after signal peptide cleavage. By SDS-PAGE, under reducing conditions, the apparent molecular mass of the protein is 42 kDa.

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