## **PRODUCT** INFORMATION



**Coagulation Factor X (human, recombinant)** 

Item No. 40252

### **Overview and Properties**

Synonyms:	Factor 10, Factor X, FX, Stuart Factor, Stuart-Prower Factor, Thrombokinase
Source:	Recombinant human C-terminal His-tagged coagulation factor X expressed in insect cells
Amino Acids:	1-488 (full length)
Uniprot No.:	P00742
Molecular Weight:	52.8 kDa
Storage:	-80°C (as supplied)
Stability:	≥1 year
Purity:	≥97% estimated by SDS-PAGE
Supplied in:	Lyophilized from sterile 20 mM Tris, 300 mM sodium chloride, 10% glycerol, pH 7.5,
	5% trehalose, 5% mannitol, and 0.01% Tween-80
Endotovin Testing	< 10 Ell/us determined by the LAL endetexin assay

**Endotoxin Testing:** < 1.0 EU/ $\mu$ g, determined by the LAL endotoxin assay Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

#### Images



Lane 1: MW Markers Lane 2: Coagulation Factor X

SDS-PAGE Analysis of Coagulation Factor X. This protein has a calculated molecular weight of 52.8 kDa. It has an apparent molecular weight of approximately 48 and 22 kDa by SDS-PAGE under reducing conditions due to glycosylation.

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, 09/17/2024

#### 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

Coagulation Factor X is a vitamin K-dependent plasma serine protease involved in the intrinsic, extrinsic, and common pathways of the blood coagulation cascade.<sup>1,2</sup> It is composed of an N-terminal light chain that contains a signal sequence, a propeptide, carboxylated glutamic acid residues (Gla domain), and two EGF domains and a C-terminal heavy chain that contains the activation peptide and a chymotrypsin-like catalytic domain.<sup>1</sup> Coagulation Factor X is produced as an inactive precursor, which is activated to Factor Xa following signal peptide cleavage, glycosylation, and further cleavage by the Factor VIIIa-IXa or tissue factor-FVIIa complexes of the intrinsic or extrinsic pathways, respectively. Factor Xa, in complex with Factor Va, converts prothrombin to thrombin. Factor Xa also cleaves severe acute respiratory syndrome coronavirus 2 (SARS-CoV-2) spike glycoprotein, also known as surface glycoprotein, and prevents SARS-CoV-2 viral entry in vitro, and Factor X levels are increased in the lungs of patients with COVID-19.<sup>3</sup> Knockout of F10, the gene encoding Factor X, in mice induces embryonic lethality and fatal internal bleeding of surviving neonates.<sup>4</sup> Mutations in F10 result in Factor X deficiency, a rare bleeding disorder characterized by CNS and gastrointestinal bleeding, as well as hemarthroses and hematomas.<sup>1</sup> Cayman's Coagulation Factor X (human, recombinant) protein consists of 468 amino acids, has a calculated molecular weight of 52.8 kDa, and a predicted N-terminus of Asn32 after signal peptide cleavage. By SDS-PAGE, under reducing conditions, the apparent molecular masses are 48 and 22 kDa, corresponding to the cleaved heavy and light chains, respectively, due to glycosylation.

#### References

- 1. Camire, R.M. Blood coagulation factor X: Molecular biology, inherited disease, and engineered therapeutics. J. Thromb. Thrombolvsis 52(2), 383-390 (2021).
- Schenone, M., Furie, B.C., and Furie, B. The blood coagulation cascade. Curr. Opin. Hematol. 11(4), 2. 272-277 (2004).
- 3. Dong, W., Wang, J., Tian, L., et al. Factor Xa cleaves SARS-CoV-2 spike protein to block viral entry and infection. Nat. Commun. 14(1), 1936 (2023).
- Dewerchin, M., Liang, Z., Moons, L., et al. Blood coagulation factor X deficiency causes partial embryonic lethality and fatal neonatal bleeding in mice. Thromb. Haemost. 83(2), 185-190 (2000).

#### 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