

PRODUCT INFORMATION



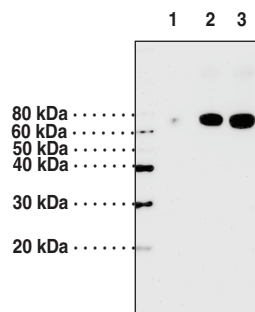
Fibrinogen (α chain) Polyclonal Antibody

Item No. 18033

Overview and Properties

Contents:	This vial contains 500 μ l of peptide affinity-purified antibody.
Synonyms:	FBG, FG, FGA
Immunogen:	Peptide from the C-terminal region of human fibrinogen (α chain)
Cross Reactivity:	(+) Fibrinogen (α chain); (-) Fibrinogen (β chain), fibrinogen (γ chain)
Species Reactivity:	(+) Human; other species not tested.
Uniprot No.:	P02671
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	\geq 1 year
Storage Buffer:	PBS, pH 7.2 with 50% glycerol and 0.02% sodium azide
Host:	Rabbit
Applications:	Western blot (WB); the recommended starting dilution is 1:200. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Image



Lane 1: Human Fibrinogen (Item No. 16088) (1 ng)
Lane 2: Human Fibrinogen (Item No. 16088) (5 ng)
Lane 3: Human Fibrinogen (Item No. 16088) (10 ng)

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, 04/01/2020

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

Fibrinogen is a hexameric glycoprotein that has roles in coagulation and hemostasis.^{1,2} It is comprised of two sets of A α , B β , and γ polypeptide chains encoded by *FGA*, *FGB*, and *FGG*, respectively, in humans.¹ Fibrinogen is synthesized in hepatocytes and secreted into the plasma. Following thrombin-mediated cleavage of N-terminal fibrinopeptides from the A α and B β chains, yielding the α and β chains, respectively, fibrinogen assembles into fibrin protofibrils and then mature fibers, which provide structure and viscoelasticity to blood clots.²⁻⁴ Mutations in *FGA*, *FGB*, or *FGG* have been found in patients with afibrinogenemia or hypofibrinogenemia.¹ Elevated plasma fibrinogen levels are associated with an increased risk of cardiovascular disease.⁵ Immune complexes containing citrullinated fibrinogen have been found in patients with anti-citrullinated protein antibody-positive rheumatoid arthritis.⁶ Cayman's Fibrinogen (α chain) Polyclonal Antibody can be used for Western blot applications. The antibody recognizes the α chain of fibrinogen at 70 kDa from human samples.

References

1. de Moerloose, P., Casini, A., and Neerman-Arbez, M. Congenital fibrinogen disorders: An update. *Semin. Thromb. Hemost.* **39(6)**, 585-595 (2013).
2. Pieters, M. and Wolberg, A.S. Fibrinogen and fibrin: An illustrated review. *Res. Pract. Thromb. Haemost.* **3(2)**, 161-172 (2019).
3. Mosesson, M.W. Fibrinogen and fibrin structure and functions. *J. Thromb. Haemost.* **3(8)**, 1894-1904 (2005).
4. Weisel, J.W. and Litvinov, R.I. Fibrin formation, structure and properties. *Subcell. Biochem.* **82**, 405-456 (2017).
5. Kamath, S. and Lip, G.Y.H. Fibrinogen: Biochemistry, epidemiology and determinants. *Q.J.M.* **96(10)**, 711-729 (2003).
6. Sokolove, J., Zhao, X., Chandra, P.E., *et al.* Immune complexes containing citrullinated fibrinogen costimulate macrophages via toll-like receptor 4 and Fc γ receptor. *Arthritis Rheum.* **63(1)**, 53-62 (2011).

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