

Fibrinogen (human plasma)

Fibrinogen is a soluble glycoprotein found in the plasma, with a molecular weight of 340kDa. It comprises of three pairs of non-identical polypeptide chains (alpha, 63.5kDa, beta, 56kDa, and gamma, 47kDa chains) linked to each other by disulphide bonds. Low plasma fibrinogen concentrations are therefore associated with an increased risk of bleeding due to impaired primary and secondary hemostasis. Therefore Fibrinogen is an essential component of the blood coagulation system. Also it may play key roles in the process of atherosclerotic lesion formation, with subsequent effects on cardiovascular diseases. And increasing evidence from epidemiological studies suggests that elevated plasma fibrinogen levels are associated with an increased risk of ischaemic heart disease (IHD), stroke and other thromboembolism.

Citations: 2

[View Online »](#)

Ordering Information

[Order Online »](#)

ENZ-PRT157-0100	100mg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Short Term Storage -80°C

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Factor 1

Formulation Liquid. In 20mM sodium citrate-HCl, pH 7.4.

MW ~330 kDa

Purity ≥95%

Source Isolated from human plasma. Plasminogen depleted.

UniProt ID P02671 (α chain), P02675 (β chain), P02679 (γ chain)

Last modified: May 29, 2024



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)