

Factor XIIIa (A Subunit) (human) (recombinant) (active) (His)

Homodimer (α_2) composed of two chains held together by non covalent bonds. After activation of the zymogen by Thrombin and Ca^{2+} to its active form (α_2^* , Factor XIIIa), Factor XIIIa catalyzes the formation of covalent bridges (ϵ -(γ -glutamyl)lysine bonds) between fibrin units to increase the elasticity of the clot network. The resulting cross-linked fibrin is insoluble and resistant to lysis.

Ordering Information

[Order Online »](#)

NZY-T070-001	1mg
NZY-T070-200	200 μ g

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Fibrin Stabilizing Factor

Application Notes Catalyzes acyl transfer reactions from glutamine residues in proteins or peptides to primary amines, e.g. the formation of ϵ -(γ -glutamyl) lysine bonds between proteins by transferring the acyl group of a peptide-bound glutamine residue to the primary amino group of a peptide-bound lysine residue.

Formulation Lyophilized from 50mM TRIS-HCl, pH 7.4, 300mM sodium chloride, 1mM CHAPS and 3mM Ca^{2+} . Sample contains sucrose.

MW ~78kDa

Purity $\geq 95\%$ (SDS-PAGE)

Reconstitution Add the desired volume of water (e.g. 400 μL) to the vial of lyophilized powder. Rotate vial gently until solid dissolves. After reconstitution the solution should be cooled on ice for short term storage.

Source Produced in insect cells.

Technical Info / Product Notes For the Original Manufacturer's data sheet please [click here](#).



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)