

Factor XIII (A subunit) (human) (recombinant) (His)

Ordering Information

[Order Online »](#)

NZY-T027-200	200µg
--------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	After reconstitution, prepare aliquots and store at -20°C. Avoid freeze/thaw cycles.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Protein-glutamine γ -glutamyltransferase, Fibrin Stabilizing Factor
Appearance	White solid.
Formulation	Lyophilized from 20mM Tris-HCl, pH 7.5, containing 150mM NaCl, 1mM EDTA, 1mM DTT, maltodextrin.
MW	83kDa monomer, 166kDa homodimer.
Purity	$\geq 95\%$ (SDS-PAGE)
Reconstitution	Reconstitute with water.
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

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com