

Transglutaminase 5 (human), (recombinant) (His)

Ordering Information

[Order Online »](#)

NZY-T086-100	100µg
--------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Protein-glutamine γ -glutamyltransferase 5, Transglutaminase X, TGX, Tgase X, Tgase-5

Appearance Clear liquid.

Formulation Diluted in 6M urea, pH9, containing 80mM TRIS-HCl and 0.5 M sodium chloride.

MW ~78kDa

Purity $\geq 90\%$ (SDS-PAGE, visually)

Source Produced in *E. coli*, inclusion body preparation.

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