

# Transglutaminase 2 (human) (recombinant) (His)

## Ordering Information

[Order Online »](#)

NZY-T022-001	1mg
NZY-T022-250	250µg

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

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

**Regulatory Status** RUO - Research Use Only

## Product Details

Alternative Name	Tissue-type Transglutaminase, TG2, TGase 2, Protein-glutamine- $\gamma$ -glutamyltransferase
Appearance	White solid.
Formulation	Lyophilized from 10mM Tris-HCl, pH 8.1, containing 150mM NaCl, 1mM EDTA, 5mM DTT, and maltodextrin.
MW	~78kDa.
Purity	≥90% (SDS-PAGE).
Reconstitution	Reconstitute with water. For activation add 10mM Ca <sup>2+</sup> .
Source	Produced in insect cells.
Specific Activity	>750U/mg protein. One unit is defined as the increase of fluorescence intensity of 1 a.u./min. Activity is determined by measuring the rate of fluorescence enhancement after TG2-catalyzed monodansylcadaverine-incorporation into N, N-dimethylated casein (L. Lorand, et al.; Anal. Biochem. <b>44</b> , 221 (1971)).
Technical Info / Product Notes	For the Original Manufacturer's data sheet please <a href="#">click here</a> .



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)