

# Transglutaminase 7 (human) (recombinant) (His)

Human recombinant His<sub>6</sub>-TG7 is based on the TGM7-gene on plasmid pCRII-hTGz cl.14, corrected by the insertion of a C at position 1169. It is N-terminally fused to a hexahistidine-tag resulting in the encoded N-terminal amino acid sequence MHHHHHHDQVATLR.... His<sub>6</sub>-TG7 is produced in *E. coli* and purified by ion metal chelating chromatography to more than 90 % purity. The enzyme is Ca<sup>2+</sup>-dependent .

## Ordering Information

[Order Online »](#)

NZY-T011-250	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

Activity	≥200 U/mg [Activity is determined by measuring the rate of fluorescence enhancement after human recombinant His <sub>6</sub> -TG7-catalyzed monodansylcadaverine-incorporation into N,N-dimethylated casein according to L. Lorand et al., Anal. Biochem. 44, 221 (1971)].Activation: Add 10mM Ca <sup>2+</sup> to activate His <sub>6</sub> -TG7.
Alternative Name	TG7
Appearance	White lyophilized solid.
Application Notes	Human recombinant His <sub>6</sub> -TG7 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. The enzyme may also be used for
MW	~81kDa
Purity	≥80% (SDS-PAGE)
Reconstitution	Add the desired volume of water to the vial of lyophilized powder.
Source	Produced in <i>E. coli</i> .
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-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)

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