

Transglutaminase 1 (human) (recombinant) (His)

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. It may also be used for immunoprecipitation.

Ordering Information

[Order Online »](#)

NZY-T035-150	150 μ g
--------------	-------------

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 Add 10mM Ca²⁺ to activate TG1.> 8,000 U/mg [Activity is determined by measuring the rate of fluorescence enhancement after His6-rhTG1-catalyzed monodansylcadaverine-incorporation into N,N-dimethylated casein according to Lorand et al., Anal. Biochem. 44 (221-231). 1 U is defined as the increase in fluorescence intensity of 1 a.u./min (measured on a Cary eclipse fluorescence spectrophotometer, Varian; λ_{ex} =332nm, λ_{em} =500nm; band filter =5nm; detector strength =600V; temperature =37°C, assay volume =1ml)].

Alternative Name Keratinocyte Transglutaminase 1, TG1, TGase 1, Keratinocyte protein-glutamine- γ -glutamyltransferase

Appearance White solid.

Formulation Lyophilized from 50mM Tris-HCl, pH 8.0, containing 10mM Glutathione.

MW ~90kDa

Purity \geq 90% (SDS-PAGE)

Purity Detail Purified by Ni²⁺-affinity chromatography.

Reconstitution Reconstitute with water. Rotate vial gently until solid dissolves. After reconstitution the solution should be cooled on ice for short term storage.

Sequence Based on the TGM1-allele from I.M.A.G.E.-clone IRAKp961M1628 isolated from human skin squamous cell carcinoma. It is N-terminally fused to His₆-tag.

Source Produced in insect cells.



ENZO LIFE SCIENCES,
ENZO LIFE SCIENCES
(ELS) AG
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
ENZO LIFE SCIENCES
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
ENZO LIFE SCIENCES
Phone: +33 1 42 20 66 55
info-fr@enzolifesciences.com

Germany
ENZO LIFE SCIENCES
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