

Transglutaminase 2 (human) (recombinant) (His) (stored in glycerol)

Ordering Information

[Order Online »](#)

| | |
|--------------|-----|
| NZY-T067-001 | 1mg |
|--------------|-----|

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

| | |
|-------------------|---|
| Handling | After opening, prepare aliquots and store at -20°C. Avoid freeze/thaw cycles. |
| Long Term Storage | -20°C |
| Shipping | Blue Ice |

Regulatory Status RUO - Research Use Only

Product Details

Activity >1500 U/mg [Activity is determined by measuring the rate of fluorescence enhancement after His₆-rhTG2-catalyzed monodansylcadaverine-incorporation into N,N-dimethylated casein according to Lorand et al., Anal. Biochem. 44 (221-231). 1U is defined as the increase in fluorescence intensity of 1 a.u./min (measured on a Cary eclipse fluorescence spectrophotometer, Varian; λ_{ex} = 332 nm, λ_{em} = 500 nm; band filter = 5 nm; detector strength = 600 V; temperature = 37°C, assay volume = 1 ml)].

Alternative Name Tissue-type Transglutaminase, TG2, TGase 2, Protein-glutamine- γ -glutamyltransferase

Application Notes 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. His₆-rhTG2 may also be used for

Formulation Liquid. In 10mM TRIS-HCl, pH 8.1, 150mM sodium chloride, 1mM EDTA , 1mM DTT and 25% Glycerol. His₆-rhTG2 is a Ca²⁺-dependent enzyme.
Activation: Add 10mM Ca²⁺ to activate His₆-rhTG2.

MW ~78kDa

Purity ≥90% (SDS-PAGE)

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