

Transglutaminase 2 (dog) (recombinant) (His)

Ordering Information

[Order Online »](#)

NZY-T072-001	1mg
NZY-T072-250	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 >1500 U/mg [Activity is determined by measuring the rate of fluorescence enhancement after His₆-cITG2-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-cITG2 may also be used for

Formulation Lyophilized from 10mM TRIS-HCl, pH 8.1, 150mM sodium chloride, 1mM EDTA and 5mM DTT. Sample contains maltodextrin. His₆-rhTG2 is a Ca²⁺-dependent enzyme. Activation: Add 10mM Ca²⁺ to activate His₆-cITG2.

MW ~78kDa

Purity \geq 90% (SDS-PAGE)

Reconstitution Add water to the vial of lyophilized powder. Rotate vial gently until solid dissolves. After reconstitution the solution should be stored frozen in working aliquots.

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