

Transglutaminase 2 (human) (recombinant) (His)

Ordering Information

[Order Online »](#)

NZY-T002-001	1mg
NZY-T002-250	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-γ-glutamyltransferase
Appearance	White solid.
Formulation	Lyophilized from 50mM NaH ₂ PO ₄ , pH 8, containing 150mM NaCl, maltodextrin and <0.1mM imidazole.
MW	~78kDa.
Purity	≥90% (SDS-PAGE).
Reconstitution	Reconstitute with 160μl water. For activation add 10mM Ca ²⁺ .
Source	Produced in <i>E. coli</i> .
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. 44 , 221 (1971)).
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