

PRODUCT INFORMATION



Transglutaminase 2 (human, recombinant)

Item No. 23595

Overview and Properties

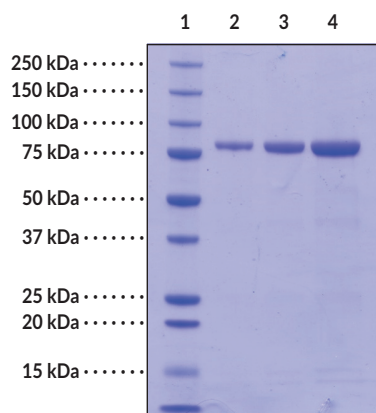
Synonyms: TG2, TGase 2, TGase C, TGase H, TGC, TGM2, Tissue Transglutaminase, Transglutaminase C, Transglutaminase H
Source: N-terminally His-tagged human TG2 protein (full length) purified from *E. coli*.
Amino acids: 2-687 (full length)
Uniprot No.: P21980
Molecular Weight: 79.4 kDa
Storage: -80°C (as supplied)
Stability: ≥1 year
Purity: *batch specific* (≥85% estimated by SDS-PAGE)
Supplied in: 50 mM HEPES, pH 7.2, with 150 mM sodium chloride, 1 mM DTT, 1 mM EDTA, and 10% glycerol

Protein

Concentration: *batch specific* mg/ml
Activity: *batch specific* U/ml
Specific Activity: *batch specific* U/mg
Unit Definition: One unit is defined as nmol of Z-glutamyl-hydroxamate-Gln produced per minute at pH 8.0 at 37°C.

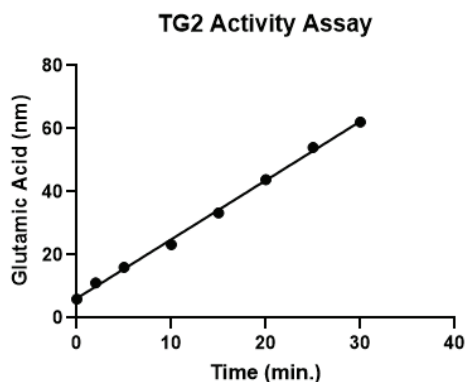
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Images



Lane 1: MW Markers
Lane 2: Transglutaminase 2 (1 µg)
Lane 3: Transglutaminase 2 (2 µg)
Lane 4: Transglutaminase 2 (4 µg)

Representative gel image shown; actual purity may vary between each batch.



TG2 activity was determined at 37°C using Cayman's Transglutaminase Activity Assay Kit (Item No 702250) with 5 µg of TG2 and 12 mM of Z-Gln-Gly substrate.

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 10/23/2025

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD

ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

Transglutaminase 2 (TG2) is the most abundant member of the transglutaminase enzyme family that is found in the intra- and extracellular spaces of various tissues.¹ It shares a common domain structure with other TGs that includes an N-terminal β -sandwich containing integrin and fibronectin binding sites, a catalytic core for acyl transfer reactions, and two C-terminal β -barrel domains with the second containing a phospholipase C binding sequence. Unlike other TGs, TG2 has a guanidine nucleotide binding site between its catalytic core and first β -barrel. TG2 catalyzes protein crosslinking in a calcium-dependent manner, creating an inter- or intramolecular bond between the ϵ -amino group of a lysine residue and the γ -carboxamide group of a glutamine residue that is highly resistant to proteolysis.² TG2 also exhibits calcium-independent enzyme activities, including GTPase, protein kinase, and disulfide isomerase activity *in vitro* and *in vivo*.³ Intracellular TG2 has important roles in protein stabilization, cytoskeletal regulation, and apoptosis.¹ It interacts with microtubule-associated protein tau-isoform Tau-F (Tau-4) and acetylcholinesterase, implicating TG2 in the pathology of neurodegenerative diseases. Extracellular TG2 has been linked to wound healing, receptor signaling, cell motility and adhesion, as well as stabilization of the extracellular matrix (ECM).³

References

1. Odi, B.O., and Coussons, P. Biological functionalities of transglutaminase 2 and the possibility of its compensation by other members of the transglutaminase family. *ScientificWorldJournal* 714561, (2014).
2. Porta, R., Esposito, C., Metafora, S., *et al.* Mass spectrometric identification of the amino donor and acceptor sites in a transglutaminase protein substrate secreted from rat seminal vesicles. *Biochemistry* 30(12), 3114-3120 (1991).
3. Belkin, A.M. Extracellular TG2: Emerging functions and regulation. *FEBS J.* 278(24), 4704-4716 (2011).