

Microbial Transglutaminase Assay Kit (chromogenic) (hydroxamate)

Transglutaminases are a family of enzymes that catalyse the posttranslational modification of proteins by inserting an isopeptide bond within or between polypeptide chains (Folk, J. E. and Cole, P. W., 1966). These enzymes catalyse the acyl transfer reaction between the γ -carboxyamide group of peptide-bound glutamine residues and a variety of primary amines, particularly the ϵ -amino group of lysine (Lorand L. et al., 1962). The resulting cross link is of great significance, since it is highly stable and also resistant to mechanical and proteolytic degradation.

Ordering Information

[Order Online »](#)

NZY-Z009-001	1Kit
--------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Application Colorimetric detection

Application Notes Determination of microbial transglutaminase activity (MTG).
Kit is sufficient for 3 x 11 measurements.

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