

# Fluorescent Transglutaminase Assay Kit

Transglutaminases catalyze acyl transfer reactions from glutamine residues in proteins or peptides to primary amines. The transglutaminase-catalyzed covalent coupling of monodansylcadaverine into N,N-dimethylcasein produces a shift in intensity and wavelength of fluorescence of the dansyl group. The transglutaminase activity can be monitored online by measurement of the fluorescence (excitation wavelength 332 nm; emission wavelength 500 nm). The relative transglutaminase activity is shown by increase of fluorescence intensity over time.

## Ordering Information

[Order Online »](#)

NZY-T036-001	1Kit
--------------	------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

<b>Use/Stability</b>	The unopened reagents are stable until the expiration date printed on the box when stored at +4°C.
<b>Handling</b>	Protect from light.
<b>Long Term Storage</b>	+4°C
<b>Shipping</b>	Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Application Notes</b>	For the determination of human transglutaminase activity (hTG1, hTG2 and hTG6).
<b>Contents</b>	Substrate reagent: 2x 0.185mg monodansylcadaverine Buffer reagent: 2x 20ml lyophilized TRIS, pH 8.0, containing calcium chloride, sodium chloride, N,N-dimethylcasein and glutathione.
<b>Technical Info / Product Notes</b>	For the Original Manufacturer's data sheet please <a href="#">click here</a> .



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)