

# Z-DON-Val-Pro-Leu-OMe

Site specific inhibitor of tissue transglutaminase.  $IC_{50} \sim 0.05 \mu M$ . Cell permeable at  $50 \mu M$ .

## Ordering Information

[Order Online »](#)

NZY-Z006-010	10mg
--------------	------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Handling	Keep cool and dry.
Long Term Storage	-20°C
Shipping	Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

Appearance	Pale yellow solid.
Formula	$C_{31}H_{44}N_6O_8$
MW	628.7
Purity	≥95% (HPLC)
Solubility	Soluble in 2% DMSO/aqueous buffer (>100µM) or in 2% v/v ethanol/aqueous buffers. Pre-dissolve e.g. 3.14 mg (5µmol) in 1 ml DMSO or 1 ml Ethanol – dilute by adding 49 ml buffer to obtain a 100 µM solution.
Technical Info / Product Notes	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)