

CTT

MMP inhibitor

CTTHWGFTLC is a specific inhibitor of the gelatinases, MMP-2 and MMP-9, with an IC_{50} of 10-50 μ M. Can be used in *in vitro* studies, zymography and *in situ* zymography, in tissue culture, and *in vivo*. Also available, negative control peptide STTHWGFTLS.

Citations: 1

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at -20°C for up to 1 year. Store solutions at -20°C for up to 3 months.
Handling	Protect from moisture.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Gelatinase inhibitor
Appearance	White to off-white solid.
Formulation	Lyophilized.
MW	1167.4
Purity	≥95% (HPLC)
Sequence	H-Cys-Thr-Thr-His-Trp-Gly-Phe-Thr-Leu-Cys-OH, cyclic peptide
Solubility	Soluble in water at 1mg/ml.
Source	Synthetic.

Last modified: May 29, 2024



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