

## MMP inhibitor

CTTHWGFTLC is a specific inhibitor of the gelatinases, MMP-2 and MMP-9, with an  $IC_{50}$  of 10-50  $\mu$ 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 »](#)

## Ordering Information

[Order Online »](#)

BML-PI136-0001	1mg
----------------	-----

## Manuals, SDS &amp; CofA

[View Online »](#)

## Handling & Storage

<b>Use/Stability</b>	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.
<b>Handling</b>	Protect from moisture.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

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



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-fr@enzolifesciences.com](mailto:info-fr@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)