

# SB-3CT

## MMP inhibitor

Competitive, mechanism-based inhibitor of matrix metalloproteinases 2 and 9 (MMP-2 and MMP-9). It possesses  $K_i$  values in the nanomolar range against the gelatinases (14nM for MMP-2; 600nM for MMP-9), but in the micromolar range against other metalloproteinases (206 $\mu$ M for MMP-1, 15 $\mu$ M for MMP-3, 96 $\mu$ M for MMP-7, and 4 $\mu$ M for ADAM17/TACE). This inhibitor has also been used *in vitro*, *in vivo*, and in tissue culture. *In vivo*, it is metabolized to an even more potent gelatinase inhibitor.

Citations: 10

[View Online »](#)

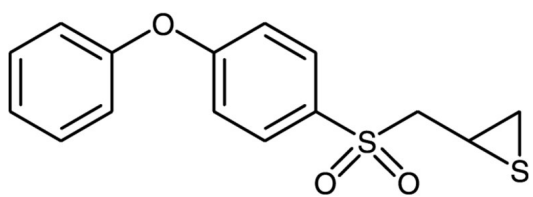
## Ordering Information

[Order Online »](#)

BML-EI325-0001	1mg
BML-EI325-0005	5mg

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

**Use/Stability** As indicated on product label or CoA when stored as recommended.

**Long Term Storage** -20°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Appearance** White solid.

**CAS** 292605-14-2

**Couple Target** MMP, MMP-2, MMP-9

**Couple Type** Inhibitor

**Formula**  $C_{15}H_{14}O_3S_2$

**MW** 306.4

**Purity** ≥98% (TLC)

**Solubility** Soluble in DMSO (at least 30mM) or 100% ethanol.



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)