# 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 »

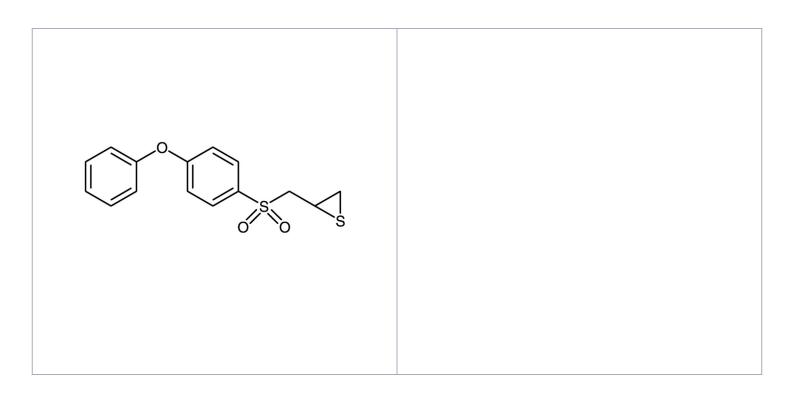
## **Ordering Information**

**Order Online** »

BML-El325-0001	1mg
BML-El325-0005	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.

Last modified: May 29, 2024