STT Negative Control for Gelatinase

MMP inhibitor

STTHWGFTLS is ten times less potent than, thus a negative control for CTTHWGFTLC (Prod. No. BML-PI136), which is a specific inhibitor of the gelatinases MMP-2 and MMP-9. Can be used in *in vitro* studies, zymography and *in situ* zymography, in tissue culture, and *in vivo*.

Citations: 3

View Online »

Ordering Information

Order Online »

BML-PI137-0005

5mg

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. After reconstitution, prepare aliquots and store at -20°C.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Appearance White to off-white solid.

Formulation Lyophilized.

MW 1136.2

Purity ≥95% (HPLC)

Sequence H-Ser-Thr-His-Trp-Gly-Phe-Thr-Leu-Ser-OH

Solubility Soluble in water (10mg/ml) or DMSO.

Source Synthetic.

Last modified: May 29, 2024



eu@enzolifesciences.com

uk@enzolifesciences.com