PF-356231

MMP inhibitor

Specific non-chelating small-molecule inhibitor of MMP-12 (IC_{50} = 14 nM). At higher concentrations it inhibits other MMPs, but spares MMP-1, making it a valuable tool for investigation of MMP activity and function.

Citations: 2

View Online »

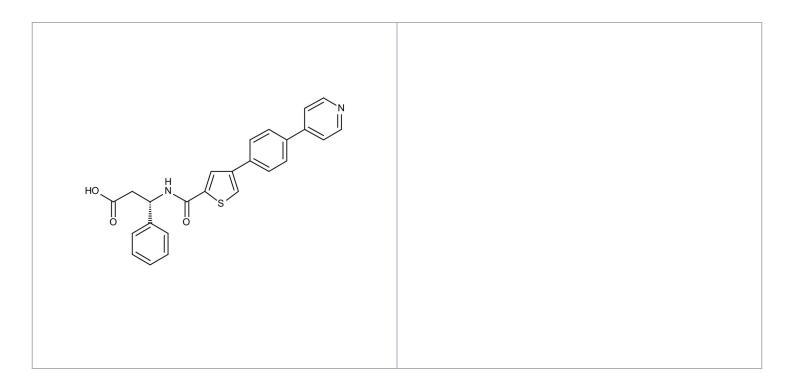
Ordering Information

Order Online »

BML-PI155-0005 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.

Long Term Storage -20°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Appearance Yellow solid.

CAS 766536-21-4

Couple Target MMP, MMP-12

Couple Type Inhibitor

 $\mathbf{Formula} \qquad \qquad \mathbf{C_{25}H_{20}N_{2}O_{3}S}$

MW 428.5

Purity ≥98% (TLC)

Soluble in DMSO (25mg/ml).

Last modified: May 29, 2024

