

PF-356231

MMP inhibitor

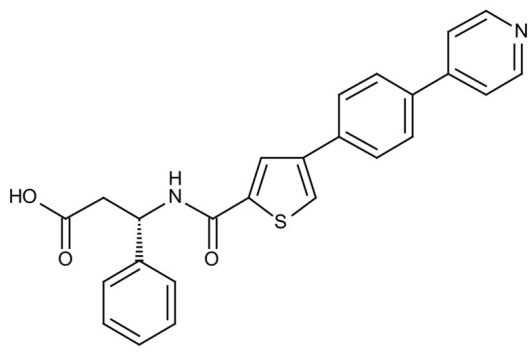
Specific non-chelating small-molecule inhibitor of MMP-12 ($IC_{50} = 14 \text{ nM}$). At higher concentrations it inhibits other MMPs, but spares MMP-1, making it a valuable tool for investigation of MMP activity and function.

Citations: 2

[View Online »](#)

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 |
| Formula | $C_{25}H_{20}N_2O_3S$ |
| MW | 428.5 |
| Purity | ≥98% (TLC) |
| Solubility | Soluble in DMSO (25mg/ml). |

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com