

WEB-2086

PAF receptor ligand

PAF receptor antagonist. Inhibits PAF-induced human platelet and neutrophil aggregation *in vitro* (IC_{50} =0.17 and 0.36 μ M respectively). *In vivo* it dose dependently blocks eosinophil activation and displays antiallergic and anti-inflammatory activity in various models. Induces differentiation and triggers apoptosis and in a variety of cell lines.

Citations: 9

[View Online »](#)

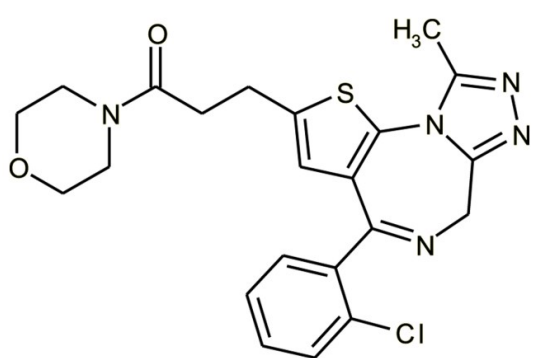
Ordering Information

[Order Online »](#)

BML-L137-0005	5mg
---------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Store dessicated.
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	Beige solid.
CAS	105219-56-5
Couple Target	PAF receptor
Couple Type	Ligand
Formula	$C_{22}H_{22}ClN_5O_2S$
Identity	Determined by 1H-NMR.
MW	456
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO or ethanol.

Last modified: May 29, 2024



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-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@enzolifesciences.com](mailto:info-fr@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)