

PD-146176

Lipoxygenase inhibitor

A potent and selective 15-lipoxygenase inhibitor which lacks significant nonspecific antioxidant properties. It suppresses atherogenesis and limits atherosclerotic lesion development in the rabbit. It reduces oxidant stress-induced apoptosis in endothelial cells and inhibits proliferation in 15-LO overexpressing PC3 cells.

Citations: 6

[View Online »](#)

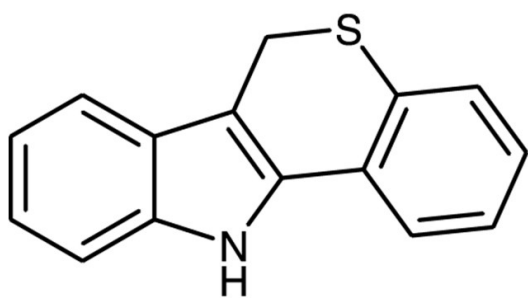
Ordering Information

[Order Online »](#)

BML-EI356-0010	10mg
BML-EI356-0050	50mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Handling Solutions should be stored at -20°C for up to 3 months.

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

CAS 4079-26-9

Couple Target Lipoxygenase

Couple Type Inhibitor

Formula $C_{15}H_{11}NS$

MW 237.3

Purity ≥98% (HPLC)

Solubility Soluble in DMSO (50mg/mL) and ethanol (10mg/mL).

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-eu@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)