

GY Y 4137

H₂S donor

Water-soluble, slow-releasing hydrogen sulfide (H₂S) donor. Shows vasodilator and antihypertensive activity in rats in both acute (L-NAME-induced) or chronic (spontaneously hypertensive) hypertension models. Inhibits LPS-induced release of various proinflammatory mediators (such as IL-1 β , IL-6, TNF- α , nitric oxide (NO) and PGE2) in RAW264.7 cells. Increases the synthesis of the anti-inflammatory chemokine IL-10 through NF- κ B, ATF-2 and HSP27 dependent pathways in RAW264.7 cells.

Citations: 8

[View Online »](#)

Ordering Information

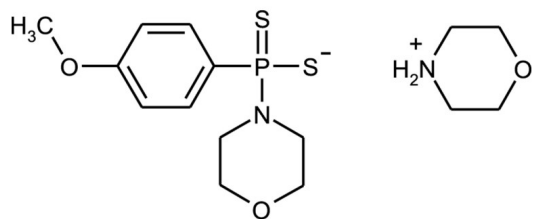
[Order Online »](#)

ALX-430-163-M010

10mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light and moisture.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	4-Methoxyphenyl(morpholino)phosphinodithioate morpholinium salt
Appearance	White solid.
CAS	106740-09-4
Formula	$C_{11}H_{15}NO_2PS_2 \cdot C_4H_{10}NO$
Identity	Identity determined by 1H-NMR.
MW	288.3 . 88.1
Purity	≥95%
Solubility	Soluble in water (30mg/ml) or DMSO (15mg/ml).
Source	Synthetic.

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-
be@enzolifesciences.com](mailto:info-be@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)