

VAS-2870

NADPH oxidase inhibitor

Novel inhibitor of NAD(P)H oxidase (Nox). In vascular smooth muscle cells VAS-2870 completely abolished PDGF-mediated NAD(P)H oxidase activation and ROS production (100% at 10 μ M). It inhibits PDGF-induced chemotaxis but not proliferation of VSMC. In mouse embryonic stem cells VAS-2870 blocked PDGF-BB-induced vasculogenesis. OxLDL-induced ROS formation in human endothelial cells was completely inhibited by VAS-2870 (100% at 5 μ M). VAS-2870 inhibits PMS-induced oxidative burst in HL-60 cells (IC₅₀=2 μ M).

Citations: 14

[View Online »](#)

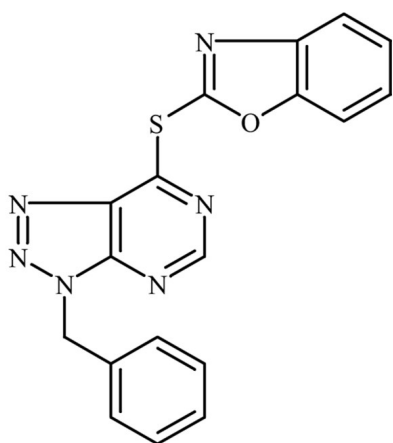
Ordering Information

[Order Online »](#)

BML-EI395-0010	10mg
BML-EI395-0050	50mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store solutions at -20°C for up to 3 months.
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	3-Benzyl-7-(2-benzoxazolyl)thio-1,2,3-triazolo(4,5-d)pyrimidine
Appearance	White to off-white solid.
CAS	722456-31-7
Couple Target	NADPH oxidase
Couple Type	Inhibitor
Formula	$C_{18}H_{12}N_6OS$
Identity	Determined by NMR.
MW	360.4
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (up to 10mg/ml) or DMF (up to 25mg/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)