

Aristolochic acid

Reversible noncompetitive PLA₂ inhibitor.

Phospholipase A₂ (PLA₂) inhibitor active against the enzymes found in many snake venoms as well as those of human platelets and synovial fluids. Inhibits ionophore-stimulated PLA₂ activity in human neutrophils. Has been shown to be nephropathic and carcinogenic.

Citations: 13

[View Online »](#)

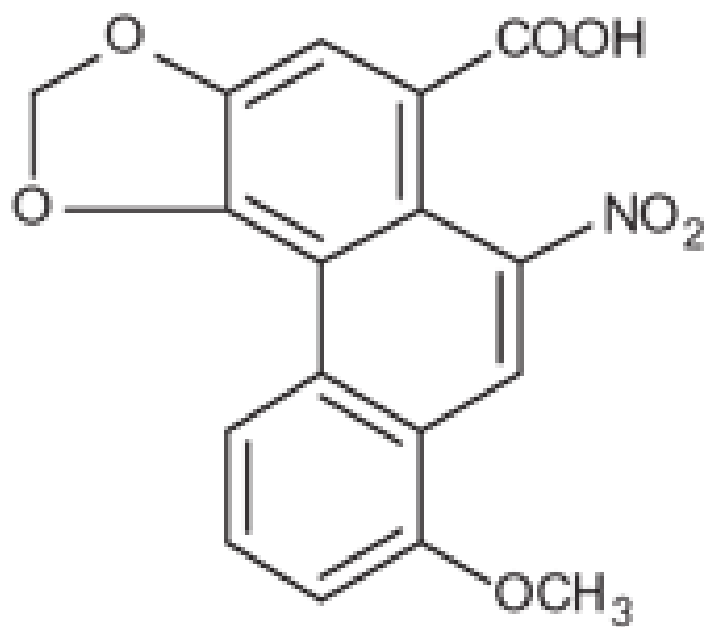
Ordering Information

[Order Online »](#)

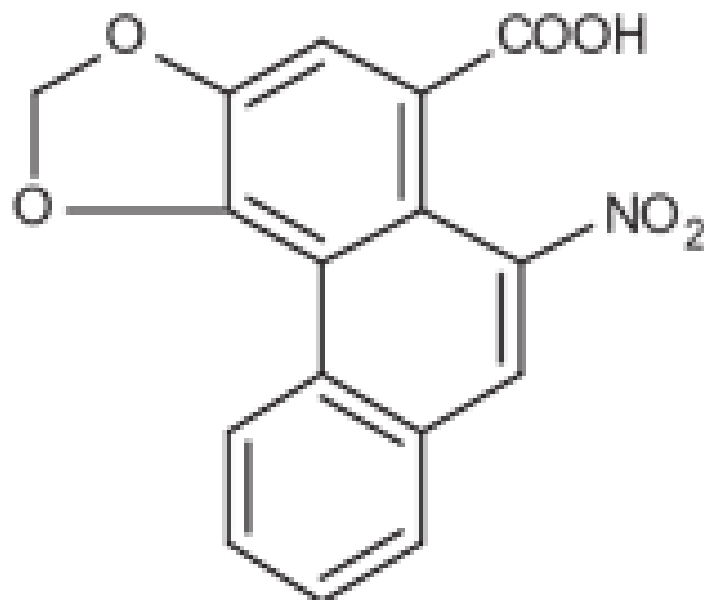
ALX-270-047-M025	25mg
------------------	------

Manuals, SDS & CofA

[View Online »](#)



Aristolochic acid I



Aristolochic acid II

Handling & Storage

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

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	8-Methoxy-6-nitrophenanthrol(3,4-d)-1,3-dioxide-5-carboxylic acid, 6-Nitrophenanthrol(3,4-d)-1,3-dioxide-5-carboxylic acid
Appearance	Yellow solid.
CAS	313-67-7 (Aristolochic acid I), 475-80-9 (Aristolochic acid II), 67123-64-2 (mixture)
Couple Target	PLA
Couple Type	Inhibitor
Formula	$C_{17}H_{11}NO_7$ (Aristolochic acid I) and $C_{16}H_9NO_6$ (Aristolochic acid II)
MI	14: 786
MW	341.3 (Aristolochic acid I) and 311.3 (Aristolochic acid II)
Purity	≥97% (~1:1 mixture of aristolochic acids I (A) and II (B))
RTECS	CF3325000
Solubility	Soluble in DMSO or 100% ethanol. Sparingly soluble in water.

Source Isolated from *Aristolochia*.



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-fr@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

Last modified: May 29, 2024