

Artemisinin

iNOS inhibitor

Powerful antimalarial agent. Inhibits angiogenesis by down-regulating HIF-1 α and VEGF expression in mouse embryonic stem cells. Crosses the blood-brain barrier. Inhibitor of human iNOS (NOS II).

Citations: 8

[View Online »](#)

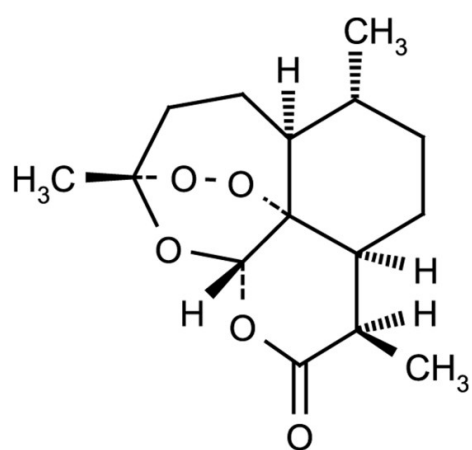
Ordering Information

[Order Online »](#)

ALX-350-219-M100	100mg
ALX-350-219-G001	1g

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	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Qinghaosu
Appearance	White to off-white crystalline solid.
CAS	63968-64-9
Couple Target	NOS
Couple Type	Inhibitor
Formula	C ₁₅ H ₂₂ O ₅
Identity	Determined by IR.
MI	14: 817
MW	282.4
Purity	≥98% (TLC)
RTECS	KD4170000
Solubility	Soluble in methanol, 100% ethanol, dimethyl formamide, DMSO or acetone; almost insoluble in water.
Source	Isolated from the traditional Chinese anti-malarial herb <i>Artemisia annua</i> L.



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)