

Rotenone

Insecticide. Inhibitor of mitochondrial electron transport. Specifically inhibits NAD-linked substrate oxidation at the oxygen side of NADH dehydrogenase ($EC_{50}=10\text{pmol/mg}$). Inhibits mammalian cell proliferation by inhibiting microtubule assembly through tubulin binding. Can inhibit mitochondrial swelling and cytochrome c release induced by N-methyl-4-phenylpyridinium.

Citations: 19

[View Online »](#)

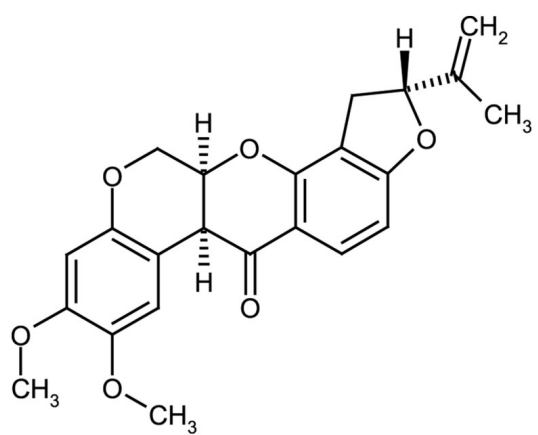
Ordering Information

[Order Online »](#)

ALX-350-360-G001	1g
ALX-350-360-G005	5g

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	White to off-white solid.
CAS	83-79-4
Couple Target	Mitochondrial complex
Couple Type	Inhibitor
Formula	$C_{23}H_{22}O_6$
MI	14: 8271
MW	394.4
Purity	≥97% (UHPLC)
RTECS	DJ2800000
Solubility	Soluble in DMSO or 100% ethanol; slightly soluble in water.
Source	Isolated from <i>Lonchocarpus</i> sp. or <i>Derris</i> sp.



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