

Monocrotaline

Pulmonary fibrosis inducer

Monocrotaline is a naturally occurring pyrrolizidine alkaloid used to create an animal model of pulmonary arterial hypertension (PAH). It can mimic several important aspects of human PAH including vascular remodeling, proliferation of smooth muscle cells, endothelial dysfunction, production of inflammatory cytokines and right ventricle failure. Induces endoplasmic reticulum stress in a rat model of PAH. Induces sinusoidal obstruction syndrome, a form of drug-induced liver injury. May be used to establish a mouse model of pulmonary fibrosis.

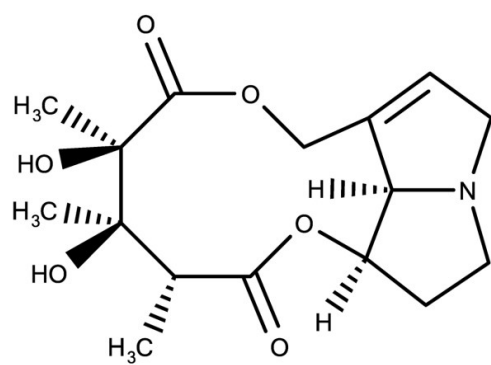
Ordering Information

[Order Online »](#)

ENZ-CHM469-0250	250mg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name (3R,4R,5R)-4,5,8,10,12,13,13aR,13bR-octahydro-4,5-dihydroxy-3,4,5-trimethyl-2H-[1,6]dioxacycloundecino[2,3,4-gh]pyrrolizine-2,6(3H)-dione

Appearance White solid.

CAS 315-22-0

Formula $C_{16}H_{23}NO_6$

Identity Determined by NMR.

MW 325.36

Purity ≥98% (HPLC)

Solubility Soluble in DMSO (50 mg/ml), ethanol (10 mg/ml) or organic solvents such as chloroform (50 mg/ml).

Last modified: February 27, 2025



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)