

Lucigenin

ROS dye

Chemiluminescent probe for the detection of superoxide in biological systems.

Citations: 12

[View Online »](#)

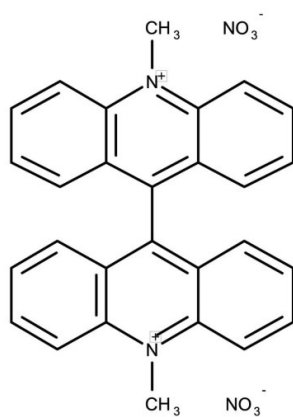
Ordering Information

[Order Online »](#)

ALX-620-061-M050	50mg
------------------	------

Manuals, SDS & CofA

[View Online »](#)



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 bis-N-Methylacridinium nitrate, N,N'-Dimethyl-9,9'-biacridinium dinitrate

Appearance Off-white to brown powder.

CAS 2315-97-1

Formula $C_{28}H_{22}N_4O_6$

MW 510.5

Purity ≥97% (HPLC)

Solubility Soluble in DMSO (~12mg/ml), 100% ethanol (~0.5mg/ml) or acetic acid (~10mg/ml, heat up solution).

Last modified: May 29, 2024



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