

Luminol

Key reagent for generating luminescence

Key reagent for generating luminescence. A reagent for the highly sensitive and selective detection of nitric oxide (NO) in biological samples by a luminol-H₂O₂ chemiluminescence method.

Citations: 4

[View Online »](#)

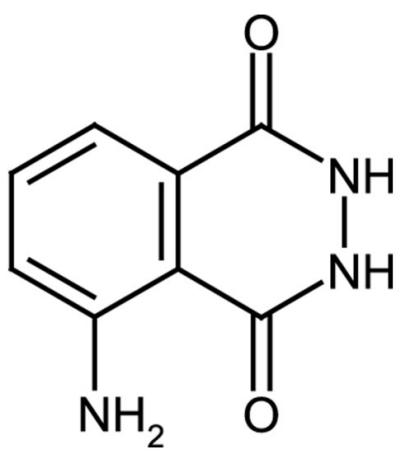
Ordering Information

[Order Online »](#)

ALX-610-002-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.

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 5-Amino-2,3-dihydro-1,4-phthalazinedione, 3-Aminophthalhydrazide

Appearance Off-white to brown solid.

CAS 521-31-3

Formula $C_8H_7N_3O_2$

MI 14: 5600

MW 177.2

Purity $\geq 97\%$ (HPLC)

RTECS TH8890060

Solubility Soluble in DMSO (50mg/ml) or in 10% ammonia (hot, 100mg/ml).

Last modified: May 29, 2024



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)