

L-Norepinephrine . bitartrate . monohydrate

Adrenergic neurotransmitter

Adrenergic neurotransmitter.

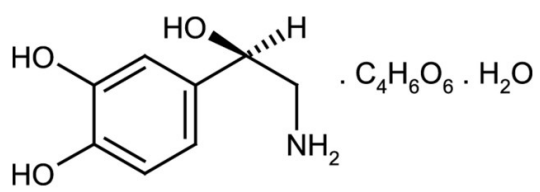
Ordering Information

[Order Online »](#)

ALX-550-076-G001	1g
------------------	----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Solutions may oxidize when exposed to air.
Handling	Protect from light.
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	L-4-(2-Amino-1-hydroxyethyl)-1,2-benzenediol . C ₄ H ₆ O ₆ . H ₂ O, L-Noradrenaline . C ₄ H ₆ O ₆ . H ₂ O
Appearance	White to off-white crystalline powder.
CAS	108341-18-0
Formula	C ₈ H ₁₁ NO ₃ . C ₄ H ₆ O ₆ . H ₂ O
MI	14: 6695
MW	169.2 . 150.1 . 18.0
Purity	≥98% (Assay)
Solubility	Soluble in water (46mg/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)