

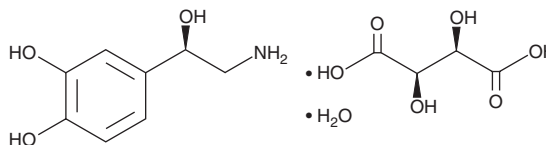
PRODUCT INFORMATION



(-)-Norepinephrine (bitartrate hydrate)

Item No. 16673

CAS Registry No.: 108341-18-0
Formal Name: 4-[(1R)-2-amino-1-hydroxyethyl]-1,2-benzenediol (2R,3R)-2,3-dihydroxybutanedioate, monohydrate
Synonyms: Arterenol, Levarterenol, NE, L-Noradrenaline
MF: C₈H₁₁NO₃ • C₄H₆O₆ • H₂O
FW: 337.3
Purity: ≥95%
UV/Vis.: λ_{max}: 224, 283 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

(-)-Norepinephrine (bitartrate hydrate) is supplied as a crystalline solid. A stock solution may be made by dissolving the (-)-norepinephrine (bitartrate hydrate) in the solvent of choice, which should be purged with an inert gas. (-)-Norepinephrine (bitartrate hydrate) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of (-)-norepinephrine (bitartrate hydrate) in these solvents is approximately 25 and 10 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of (-)-norepinephrine (bitartrate hydrate) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of (-)-norepinephrine (bitartrate hydrate) in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

(-)-Norepinephrine is the endogenous isomer of (±)-norepinephrine (Item No. 35580) that acts as a neurotransmitter and hormone and is biosynthesized from dopamine by dopamine β-hydroxylase.¹⁻³ It is an agonist of adrenergic receptors (K_i values are 330, 56, and 740 nM for α₁, α₂, and β₁ adrenoceptors, respectively).³ Through these receptors, (-)-norepinephrine regulates diverse processes, including those in neurological, immunological, and vascular systems.³⁻⁶ Formulations containing (-)-norepinephrine have been used for the restoration of blood pressure in acute hypotensive states.

References

1. von Euler, U.S. *Pharmacol. Rev.* **3(3)**, 247-277 (1951).
2. Gannon, M., Che, P., Chen, Y., et al. *Frontiers in Neuroscience* **9**, 1-12 (2015).
3. Ramos, B.P. and Arnsten, A.F.T. *Pharmacol. Ther.* **113(3)**, 523-536 (2007).
4. Lorton, D. and Bellinger, D.L. *Int. J. Mol. Sci.* **16(3)**, 5635-5665 (2015).
5. Nance, D.M. and Sanders, V.M. *Brain Behav. Immun.* **21(6)**, 736-745 (2007).
6. Raiteri, M. *Pharmacol. Rev.* **58(2)**, 162-193 (2006).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 04/10/2023

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM