

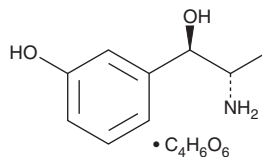
PRODUCT INFORMATION



Metaraminol (tartrate)

Item No. 23938

CAS Registry No.: 33402-03-8
Formal Name: αR-[(1S)-1-aminoethyl]-3-hydroxy-benzenemethanol, 2R,3R-dihydroxybutanedioate)
MF: C₉H₁₃NO₂ • C₄H₆O₆
FW: 317.3
Purity: ≥98%
Supplied as: A 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

Metaraminol (tartrate) is supplied as a solid. A stock solution may be made by dissolving the metaraminol (tartrate) in the solvent of choice, which should be purged with an inert gas. Metaraminol (tartrate) is slightly soluble in DMSO.

Description

Metaraminol is a vasoconstrictor.^{1,2} *In vivo*, metaraminol (0.05 mg/kg) increases cardiac output, mean arterial pressure, and coronary blood flow and decreases pulse rate in a canine model of cardiogenic shock.¹ It increases survival in a canine model of hemorrhagic hypotension.² Metaraminol also reverses acetoxycycloheximide-induced transient amnesia in a shock avoidance test in rats.³

References

1. Sarnoff, S.J., Case, R.B., Berglund, E., *et al.* Ventricular function. V. The circulatory effects of aramine; mechanism of action of vasopressor drugs in cardiogenic shock. *Circulation* **10(1)**, 84-93 (1954).
2. Sarnoff, S.J. and Kaufman, H.E. Effect of aramine on the survival of dogs subjected to hemorrhagic hypotension. *Circ. Res.* **2(5)**, 420-423 (1954).
3. Serota, R.G., Roberts, R.B., and Flexner, L.B. Acetoxycycloheximide-induced transient amnesia: Protective effects of adrenergic stimulants. *Proc. Natl. Acad. Sci. U.S.A.* **69(2)**, 340-342 (1972).

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, 11/17/2022

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