

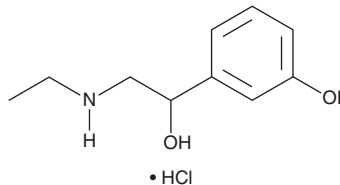
PRODUCT INFORMATION



Etilefrine (hydrochloride)

Item No. 33249

CAS Registry No.: 943-17-9
Formal Name: α -[(ethylamino)methyl]-3-hydroxy-benzenemethanol, monohydrochloride
Synonyms: Circupon, Ethylephrine, Ethylphenylephrine
MF: $C_{10}H_{15}NO_2 \cdot HCl$
FW: 217.7
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 217 nm
Supplied as: A crystalline solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Laboratory Procedures

Etilefrine (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the etilefrine (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Etilefrine (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of etilefrine (hydrochloride) in these solvents is approximately 5 mg/ml in ethanol and approximately 30 mg/ml in DMSO and DMF.

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 etilefrine (hydrochloride) can be prepared by directly dissolving the etilefrine (hydrochloride) in aqueous buffers. The solubility of etilefrine (hydrochloride) in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Etilefrine is a sympathomimetic amine.¹ It induces vasoconstriction in isolated perfused rat tail artery segments in a concentration-dependent manner. Etilefrine (45 μ g/kg, i.v.) inhibits anesthesia-induced decreases in mean arterial pressure and cardiac output in dogs.² Formulations containing etilefrine have been used in the treatment of orthostatic hypotension.

References

1. Frost, B.R., Frewin, D.B., and Gerke, D.C. The effects of etilefrine on blood vessels in the rat tail. *J. Pharm. Pharmacol.* **29(5)**, 272-275 (1977).
2. Tarnow, J., Brückner, J.B., Eberlein, H.J., et al. Haemodynamic responses to ethylphenylephrine (Effortil) during halothane-induced myocardial depression in the dog. *Acta Anaesthesiol. Scand.* **17(3)**, 184-189 (1973).

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, 10/06/2022

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