

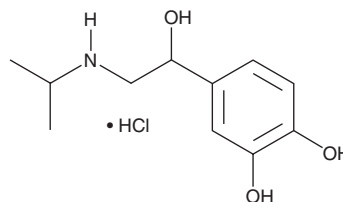
PRODUCT INFORMATION



Isoproterenol (hydrochloride)

Item No. 15592

CAS Registry No.: 51-30-9
Formal Name: 4-[1-hydroxy-2-[(1-methylethyl)amino]ethyl]-1,2-benzenediol, monohydrochloride
Synonyms: Isoprenaline, NSC 37745, NSC 89747
MF: C₁₁H₁₇NO₃ • HCl
FW: 247.7
Purity: ≥98%
UV/Vis.: λ_{max}: 225, 282 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

Isoproterenol (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the isoproterenol (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Isoproterenol (hydrochloride) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of isoproterenol (hydrochloride) in these solvents is approximately 10 and 2 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 isoproterenol (hydrochloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of isoproterenol (hydrochloride) in PBS (pH 7.2) is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Isoproterenol is an agonist of β_1 - and β_2 -adrenergic receptors (β_1 - and β_2 -ARs; K_i s = 224 and 458 nM, respectively).¹ It is selective for β_1 - and β_2 -ARs over β_3 -ARs (K_i = 1,570 nM). Isoproterenol inhibits contractions in isolated field-stimulated rat vas deferens (EC_{50} = 45.6 nM).² *In vivo*, isoproterenol (0.33 mg/kg) decreases blood pressure and increases water intake in nephrectomized rats.³ It reduces blood pressure and increases heart rate in renal hypertensive rabbits.⁴ Isoproterenol inhibits histamine-induced bronchospasms in anesthetized dogs.⁵ Formulations containing isoproterenol have been used in the treatment of bradycardias and to improve breathing during anesthesia.

References

1. Hoffmann, C., Leitz, M.R., Oberdorf-Maass, S., et al. *Naunyn Schmiedebergs Arch. Pharmacol.* **369**(2), 151-159 (2004).
2. Lotti, V.J., Cerino, D., and Kling, P. *J. Auton. Pharmacol.* **2**(3), 169-174 (1982).
3. Hosutt, J.A., Rowland, N., and Stricker, E.M. *Physiol. Behav.* **21**(4), 593-598 (1978).
4. van Boom, M. and Saxena, P.R. *Clin. Exp. Pharmacol. Physiol.* **8**(3), 227-239 (1981).
5. Wasserman, M.A. and Levy, B. *J. Pharmacol. Exp. Ther.* **189**(2), 445-455 (1974).

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