

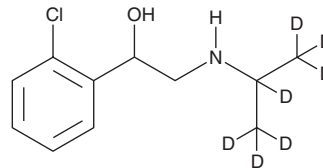
PRODUCT INFORMATION



(±)-Clorprenaline-d₇

Item No. 27614

CAS Registry No.: 2748354-86-9
Formal Name: 1-(2-chlorophenyl)-2-((propan-2-yl-d₇)amino)ethan-1-ol
Synonym: Isoprophenamine-d₇
MF: C₁₁H₉ClD₇NO
FW: 220.7
Chemical Purity: ≥98% (Clorprenaline)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₇); ≤1% d₀
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

(±)-Clorprenaline-d₇ is intended for use as an internal standard for the quantification of clorprenaline by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

(±)-Clorprenaline-d₇ is supplied as a solid. A stock solution may be made by dissolving the (±)-clorprenaline-d₇ in the solvent of choice, which should be purged with an inert gas. (±)-Clorprenaline-d₇ is soluble in the organic solvent chloroform.

Description

Clorprenaline is a β_2 -adrenergic receptor agonist.¹ It reverses pilocarpine-induced constriction of isolated guinea pig trachea, rabbit lung, and bronchi of pithed dogs.² It has been detected in pork, beef, mutton, and chicken muscle food samples by UPLC-Q-orbitrap-HRMS.¹ Formulations containing clorprenaline have been used as partitioning agents and growth promoters in livestock.

References

1. Guo, C., Shi, F., Gong, L., *et al.* Ultra-trace analysis of 12 β_2 -agonists in pork, beef, mutton and chicken by ultrahigh-performance liquid-chromatography-quadrupole-orbitrap tandem mass spectrometry. *J. Pharm. Biomed. Anal.* **107**, 526-534 (2015).
2. Powell, C.E., Gibson, W.R., and Swanson, E.E. The pharmacological action of 1-o-chlorophenyl-2-isopropylaminoethanol (isoprophenamine); A bronchodilator. *J. Am. Pharm. Assoc.* **45(12)**, 785-787 (1956).

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, 02/10/2023

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