

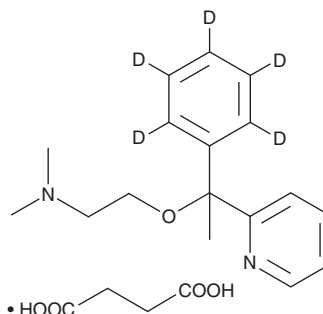
PRODUCT INFORMATION



Doxylamine-d₅ (succinate)

Item No. 40674

CAS Registry No.: 1216840-94-6
Formal Name: butanedioic acid, compd. with N,N-dimethyl-2-[1-phenyl-d₅-1-(2-pyridinyl)ethoxy]ethanamine
MF: C₁₇H₁₇D₅N₂O • C₄H₆O₄
FW: 393.5
Chemical Purity: ≥95% (Doxylamine)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₅); ≤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

Doxylamine-d₅ (succinate) is intended for use as an internal standard for the quantification of doxylamine 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).

Description

Doxylamine is an antihistamine.^{1,2} It selectively inhibits contractions induced by histamine (Item No. 33828) over acetylcholine (ACh; Item No. 23829) in isolated guinea pig ileum (IC₅₀s = 0.01 and 0.58 μg/ml, respectively).³ Doxylamine (2-4 mg/kg) protects rabbits against anaphylaxis induced by beef serum antigen.² It reduces ethanol-induced deficits in behavior performance in rats when administered at a dose of 60 mg/kg.⁴ Formulations containing doxylamine have been used in the treatment of nausea and vomiting in pregnant women and insomnia.

References

1. Brown, B.B., Werner, H.W., and Peters, E.L. The pharmacologic properties of 2-[α-(2-dimethylaminoethoxy)-α-methylbenzyl]-pyridine succinate, a new antihistaminic agent. *J. Lab. Clin. Med.* **33**(3), 325-331 (1948).
2. Brown, B.B. and Werner, H.W. The effects of decapryl succinate, a new antihistamine agent, in some natural and acquired hypersensitivities in animals; 2-[α-dimethylaminoethoxy)-α-methyl-benzyl] pyridine succinate. *Ann. Allergy* **6**(2), 122-130 (1948).
3. Werle, E. and Lorenz, W. The antikinin action of some antihistaminic drugs on the isolated guinea-pig ileum, rat uterus and blood pressure of the anesthetized dog. *Bradykinin and Related Kinins Cardiovascular, Biochemical, and Neural Actions. Advances in Experimental Medicine and Biology*. Scuteri, F., Rocha e Silva, M., and Back, N., editors, Plenum Press, New York, New York (1970).
4. Allen, L.E., Ferguson, H.C., and McKinney, G.R. A survey of selected drugs on behavioral performance in ethanol-treated rats. *Eur. J. Pharmacol.* **15**(3), 371-374 (1971).

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, 03/21/2024

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