

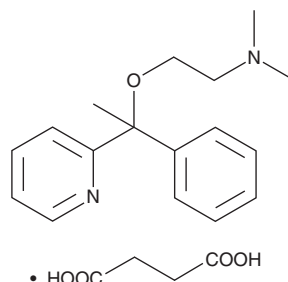
PRODUCT INFORMATION



Doxylamine (succinate)

Item No. 21826

CAS Registry No.: 562-10-7
Formal Name: butanedioic acid, compd. with
N,N-dimethyl-2-[1-phenyl-1-(2-
pyridinyl)ethoxy]ethanamine
Synonym: NSC 74772
MF: $C_{17}H_{22}N_2O \cdot C_4H_6O_4$
FW: 388.5
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 261 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.

Description

Doxylamine (succinate) (Item No. 21826) is an analytical reference standard categorized as an antihistamine and hypnotic.¹ Overdose of doxylamine (succinate) is common and has been associated with rhabdomyolysis.^{1,2} This product is intended for research and forensic applications.

References

1. Soto, L.F., Miller, C.H., and Ognibene, A.J. Severe rhabdomyolysis after doxylamine overdose. *Postgrad. Med.* **93(8)**, 227-229, 232 (1993).
2. Leybushkis, B., Fasseas, P., and Ryan, K.F. Doxylamine overdose as a potential cause of rhabdomyolysis. *Am. J. Med. Sci* **322(1)**, 48-49 (2001).

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, 03/25/2024

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