

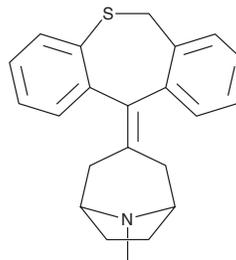
PRODUCT INFORMATION



Tropatepine

Item No. 43415

CAS Registry No.: 27574-24-9
Formal Name: 3-dibenzo[b,e]thiepin-11(6H)-ylidene-8-methyl-8-azabicyclo[3.2.1]octane
MF: C₂₂H₂₃NS
FW: 333.5
Purity: ≥95%
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

Tropatepine is supplied as a solid. A stock solution may be made by dissolving the tropatepine in the solvent of choice, which should be purged with an inert gas. Tropatepine is soluble in methanol and DMSO.

Description

Tropatepine is an anticholinergic agent.^{1,2} It protects against sinistrotorsion syndrome induced by oxotremorine (Item No. 35707) in guinea pigs (ED₅₀ = 0.2-0.5 mg/kg).² Formulations containing tropatepine have been used in the treatment of neuroleptic-induced parkinsonian symptoms.

References

1. Nigam, M., Leu-Semenescu, S., and Arnulf, I. Successful treatment of drug-resistant cataplexy with the anticholinergic drug tropatepine. *J. Clin. Sleep Med.* **17(4)**, 849-851 (2021).
2. Boissier, J.R., and Ratouis, R. New tropane derivatives and their salts. **PV170767** (1970).

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, 05/13/2025

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