

PRODUCT INFORMATION

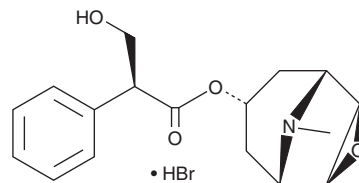


Scopolamine (hydrobromide)

Item No. 14108

CAS Registry No.: 114-49-8
Formal Name: (1 α ,2 β ,4 β ,5 α ,7 β)-9-methyl-3-oxa-9-azatricyclo[3.3.1.0^{2,4}]non-7-yl ester α S-(hydroxymethyl)-benzeneacetic acid, monohydrobromide

Synonym: Hyoscine
MF: C₁₇H₂₁NO₄ • HBr
FW: 384.3
Purity: \geq 95%
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 5 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

Scopolamine is a tropane alkaloid that can be found in many plants of the *Solanaceae* (nightshade) family.¹ It is a muscarinic receptor antagonist that can be used to induce memory impairment in animals.²⁻⁴ Scopolamine prevents motion sickness, nausea, and vomiting in animals.^{3,5}

References

1. Ziegler, J. and Facchini, P.J. *Annu. Rev. Plant Biol.* **59**, 735-769 (2008).
2. Woods, S., Clarke, N.N., Layfield, R., et al. *Brit. J. Pharmacol.* **167**, 436-449 (2012).
3. Soto, E. and Vega, R. *Curr. Neuropharmacol.* **8(1)**, 26-40 (2010).
4. Shannon, H.E., Hart, J.C., Bymaster, F.P., et al. *J. Pharmacol. Exp. Ther.* **290(2)**, 901-907 (1999).
5. Gryniewicz, G. and Gadzikowska, M. *Pharmacol. Rep.* **60**, 439-463 (2008).

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, 06/15/2023

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