

PRODUCT INFORMATION



Aposcopolamine

Item No. 36689

CAS Registry No.: 535-26-2

Formal Name: α -methylene-benzeneacetic acid,
(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

Synonym: Apohyoscine

MF: C₁₇H₁₉NO₃

FW: 285.3

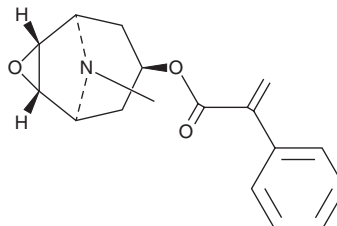
Purity: $\geq 95\%$

Supplied as: A solid

Storage: -20°C

Stability: ≥ 4 years

Item Origin: Synthetic



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

Laboratory Procedures

Aposcopolamine is supplied as a solid. A stock solution may be made by dissolving the aposcopolamine in the solvent of choice, which should be purged with an inert gas. Aposcopolamine is slightly soluble in DMSO (warmed) and methanol (sonicated).

Description

Aposcopolamine is a tropane alkaloid that has been found in *D. ferox* and is an active metabolite of the muscarinic acetylcholine receptor (mAChR) antagonist scopolamine (Item No. 40307).^{1,2} It is selective for muscarinic acetylcholine receptors (mAChRs) over nicotinic AChRs (nAChRs; IC₅₀s = 0.0192 and 188 μ M, respectively).³

References

1. Vitale, A.A., Acher, A., and Pomilio, A.B. Alkaloids of *Datura ferox* from Argentina. *J. Ethnopharmacol.* **49**(2), 81-89 (1995).
2. Chen, H., Chen, Y., Du, P., *et al.* Liquid chromatography-electrospray ionization ion trap mass spectrometry for analysis of in vivo and in vitro metabolites of scopolamine in rats. *J. Chromatogr. Sci.* **46**(1), 74-80 (2008).
3. Schmeller, T., Sporer, F., Sauerwein, M., *et al.* Binding of tropane alkaloids to nicotinic and muscarinic acetylcholine receptors. *Pharmazie* **50**(7), 493-495 (1995).

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, 03/26/2024

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