

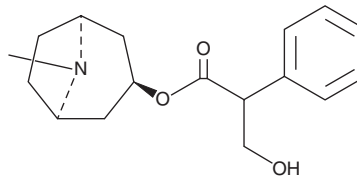
PRODUCT INFORMATION



Atropine

Item No. 12008

CAS Registry No.: 51-55-8
Formal Name: α -(hydroxymethyl)-benzeneacetic acid, (3-*endo*)-8-methyl-8-azabicyclo[3.2.1]oct-3-yl ester
Synonyms: DL-Hyoscyamine, Tropicine tropate
MF: C₁₇H₂₃NO₃
FW: 289.4
Purity: \geq 95%
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

Atropine is supplied as a crystalline solid. A stock solution may be made by dissolving the atropine in the solvent of choice, which should be purged with an inert gas. Atropine is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of atropine in these solvents is approximately 16, 10, and 2 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of atropine can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of atropine in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Atropine is a naturally occurring tropane alkaloid extracted from plants of the family Solanaceae including deadly nightshade (*A. belladonna*). It is a non-selective, competitive antagonist of the muscarinic acetylcholine receptor types M₁, M₂, M₃, M₄, and M₅ (pK_Bs range from 8.9-9.8).¹ Atropine increases firing of the sinoatrial node and conduction through the atrioventricular node of the heart, opposes the actions of the vagus nerve, blocks acetylcholine receptor sites, and decreases bronchial secretions.² It is classified as an anticholinergic (parasympatholytic) drug and commonly used to dilate the pupils, increase heart rate, reduce salivation and other secretions, and as an antidote against organophosphate poisoning.³

References

1. Caulfield, M.P. and Birdsall, N.J.M. International union of pharmacology. XVII. Classification of muscarinic acetylcholine receptors. *Pharmacol. Rev.* **50(2)**, 279-290 (1998).
2. Broadley, K.J. and Kelly, D.R. Muscarinic receptor agonists and antagonists. *Molecules* **6(3)**, 142-193 (2001).
3. Gryniewicz, G. and Gadzikowska, M. Tropane alkaloids as medicinally useful natural products and their synthetic derivatives as new drugs. *Pharmacol. Rep.* **60(4)**, 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, 07/20/2023

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