

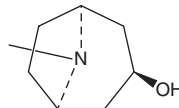
PRODUCT INFORMATION



Tropine

Item No. 18627

CAS Registry No.: 120-29-6
Formal Name: (3-endo)-8-methyl-8-Azabicyclo[3.2.1]octan-3-ol
Synonyms: NSC 43870, 3-Tropanol
MF: C₈H₁₅NO
FW: 141.2
Purity: ≥95%
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.

Laboratory Procedures

Tropine is supplied as a crystalline solid. A stock solution may be made by dissolving the tropine in the solvent of choice, which should be purged with an inert gas. Tropine is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of tropine in these solvents is approximately 30 mg/ml.

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 tropine can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of tropine in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Tropine is a naturally-occurring tropane alkaloid extracted primarily from plants of the *Solanaceae* family.¹ It serves as an intermediate in the synthesis of a variety of bioactive alkaloids, including atropine (Item No. 12008), benztropine (Item No. 18214), and scopolamine, many of which have potent neurological actions.^{1,2}

References

1. Gryniewicz, G. and Gadzikowska, M. Tropane alkaloids as medicinally useful natural products and their synthetic derivatives as new drugs. *Pharmacol. Rep.* **60**, 439-463 (2008).
2. Richter, U., Rothe, G., Fabian, A.-K., *et al.* Overexpression of tropinone reductases alters alkaloid composition in *Atropa belladonna* root cultures. *J. Exp. Bot.* **56(412)**, 645-652 (2005).

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, 12/05/2022

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