

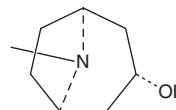
PRODUCT INFORMATION



Pseudotropine

Item No. 36166

CAS Registry No.: 135-97-7
Formal Name: (3-exo)-8-methyl-8-azabicyclo[3.2.1]octan-3-ol
Synonyms: NSC 43871, Pseudotropanol, 3 β -Tropanol, Ψ -Tropane
MF: C₈H₁₅NO
FW: 141.2
Purity: \geq 98%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years
Item Origin: Synthetic



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

Laboratory Procedures

Pseudotropine is supplied as a solid. A stock solution may be made by dissolving the pseudotropine in the solvent of choice, which should be purged with an inert gas. Pseudotropine is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of pseudotropine in these solvents is approximately 10, 2, and 1 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 pseudotropine can be prepared by directly dissolving the solid in aqueous buffers. The solubility of pseudotropine in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Pseudotropine is an alkaloid and a plant metabolite of the tropane tropinone (Item No. 34287) that has been found in *A. belladonna*.¹ It is biosynthesized from tropinone by tropinone reductase II. Pseudotropine has been found as a contaminant in foodstuffs due to coharvesting of pseudotropine-containing weeds in agriculture.²

References

1. 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).
2. de Nijs, M., Crews, C., Dorgelo, F., *et al.* Emerging issues on tropane alkaloid contamination of food in Europe. *Toxins (Basel)* **15**(2), 98 (2023).

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, 06/12/2023

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