

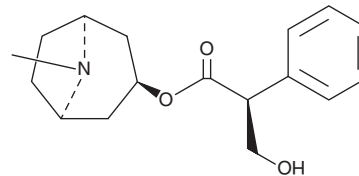
# PRODUCT INFORMATION



## (-)-Hyoscyamine

Item No. 25644

**CAS Registry No.:** 101-31-5  
**Formal Name:** (αS)-α-(hydroxymethyl)-benzeneacetic acid, (3-endo)-8-methyl-8-azabicyclo[3.2.1]oct-3-yl ester  
**Synonyms:** L-Hyoscyamine, (S)-Hyoscyamine  
**MF:** C<sub>17</sub>H<sub>23</sub>NO<sub>3</sub>  
**FW:** 289.4  
**Purity:** ≥98%  
**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

(-)-Hyoscyamine is supplied as a crystalline solid. A stock solution may be made by dissolving the (-)-hyoscyamine in the solvent of choice. (-)-Hyoscyamine is soluble in the organic solvent chloroform, which should be purged with an inert gas, at a concentration of approximately 50 mg/ml. (-)-Hyoscyamine is also soluble in ethanol and methanol.

### Description

(-)-Hyoscyamine is a tropane alkaloid, a muscarinic receptor antagonist, and an isomer of atropine (Item No. 12008).<sup>1,2</sup> (-)-Hyoscyamine inhibits the arterial depressor response induced by the muscarinic agonist methacholine in dogs.<sup>1</sup> It is more effective than atropine at inhibiting EEG arousal in dogs following sciatic nerve stimulation when administered at doses of 0.5 and 1 mg/kg. (-)-Hyoscyamine induces convulsions in and is toxic to mice (LD<sub>50</sub> = 95 mg/kg).<sup>2</sup>

### References

1. Domino, E.F. and Hudson, R.D. Observations on the pharmacological actions of the isomers of atropine. *J. Pharmacol. Exp. Ther.* **127(4)**, 305-312 (1959).
2. European Food Safety Authority. Scientific Opinion on Tropane alkaloids in food and feed EFSA Panel on contaminants in the food chain (CONTAM). *EFSA J.* **11(10)**, 3386 (2013).

#### 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, 11/10/2022

#### 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