

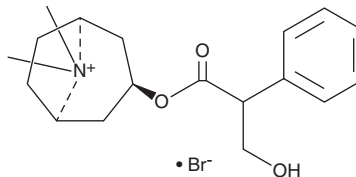
# PRODUCT INFORMATION



## Methylatropine (bromide)

Item No. 33433

**CAS Registry No.:** 2870-71-5  
**Formal Name:** 3-(3-hydroxy-1-oxo-2-phenylpropoxy)-8,8-dimethyl-8-azoniabicyclo[3.2.1]octane, monobromide  
**Synonym:** NSC 61810  
**MF:** C<sub>18</sub>H<sub>26</sub>NO<sub>3</sub> • Br  
**FW:** 384.3  
**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

Methylatropine (bromide) is supplied as a crystalline solid. A stock solution may be made by dissolving the methylatropine (bromide) in water. The solubility of methylatropine (bromide) in water is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Methylatropine is an antagonist of muscarinic acetylcholine receptors (IC<sub>50</sub> = <0.1 nM in a radioligand binding assay using isolated porcine brain membranes) and a derivative of atropine (Item No. 12008).<sup>1,2</sup> It reduces acetylcholine-induced decreases in blood pressure in rats when administered intravenously with an ED<sub>50</sub> value of 5.5 µg/kg.<sup>2</sup> Methylatropine reduces salivation, induces mydriasis, and increases heart rate in dogs.<sup>3</sup>

### References

- Schmeller, T., Sporer, F., Sauerwein, M., *et al.* Binding of tropane alkaloids to nicotinic and muscarinic acetylcholine receptors. *Pharmazie* **50(7)**, 493-495 (1995).
- Brezenoff, H.E., Xiao, Y.-F., and Vargas, H. A comparison of the central and peripheral antimuscarinic effects of atropine and methylatropine injected systemically and into the cerebral ventricles. *Life Sci.* **42(8)**, 905-911 (1988).
- Albanus, L. Central and peripheral effects of anticholinergic compounds. *Acta Pharmacol. Toxicol. (Copenh)* **28(4)**, 305-326 (1970).

#### 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, 11/04/2022

#### 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