

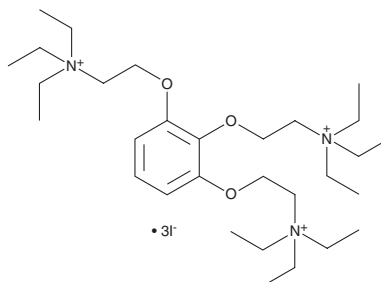
# PRODUCT INFORMATION



## Gallamine (iodide)

Item No. 26086

**CAS Registry No.:** 65-29-2  
**Formal Name:** 2,2',2''-[1,2,3-benzenetriyl tris(oxy)]tris[N,N,N-triethyl-ethanaminium, triiodide  
**MF:** C<sub>30</sub>H<sub>60</sub>N<sub>3</sub>O<sub>3</sub> • 3I  
**FW:** 891.5  
**Purity:** ≥95%  
**Supplied as:** A 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

Gallamine (iodide) is supplied as a solid. A stock solution may be made by dissolving the gallamine (iodide) in the solvent of choice, which should be purged with an inert gas. Gallamine (iodide) is slightly soluble in methanol and DMSO.

Gallamine (iodide) is slightly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

### Description

Gallamine is an allosteric modulator of muscarinic acetylcholine receptors.<sup>1</sup> It slows the dissociation of the muscarinic antagonist [<sup>3</sup>H]NMS from M<sub>1</sub>-M<sub>5</sub> muscarinic receptors, with the highest potency for M<sub>2</sub> receptors. Gallamine is cardioselective, inhibiting [<sup>3</sup>H]NMS binding in rat heart tissue homogenates more potently than in other tissues (IC<sub>50</sub>s = 0.8 and 3-30 μM, respectively).<sup>2</sup> Formulations containing gallamine have previously been used as neuromuscular blockers during surgery.

### References

1. Ellis, J., Huyler, J., and Brann, M.R. Allosteric regulation of cloned m1-m5 muscarinic receptor subtypes. *Biochem. Pharmacol.* **42**(10), 1927-1932 (1991).
2. Stockton, J.M., Birdsall, J.M., Burgen, A.S.V., et al. Modification of the binding properties of muscarinic receptors by gallamine. *Mol. Pharmacol.* **23**(3), 551-557 (1983).

#### 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, 10/28/2022

#### 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](http://WWW.CAYMANCHEM.COM)