

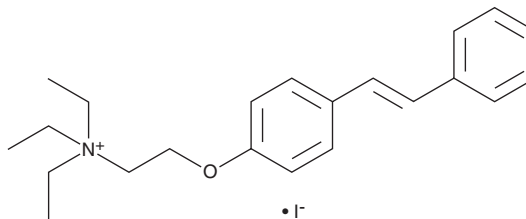
# PRODUCT INFORMATION



## MG624

Item No. 24298

**CAS Registry No.:** 77257-42-2  
**Formal Name:** N,N,N-triethyl-2-[4-[(1E)-2-phenylethenyl]phenoxy]-ethanaminium, monoiodide  
**MF:** C<sub>22</sub>H<sub>30</sub>NO • I  
**FW:** 451.4  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 204, 224, 302 nm  
**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

MG624 is supplied as a crystalline solid. A stock solution may be made by dissolving the MG624 in the solvent of choice, which should be purged with an inert gas. MG624 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of MG624 in DMSO and DMF is approximately 20 mg/ml and approximately 10 mM in ethanol.

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 MG624 can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of MG624 in PBS (pH 7.2) is approximately 0.3 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

MG624 is an antagonist of neuronal nicotinic acetylcholine receptors (nAChRs; K<sub>i</sub> = 0.055 μM in neuronal chick optic lobe membranes).<sup>1</sup> It is selective for neuronal nAChRs over muscle-type AChRs (K<sub>i</sub> = 70 μM in TE671 cells that express muscle-type AChRs). MG624 inhibits currents evoked by acetylcholine in *Xenopus* oocytes expressing chick α7 subunit-containing nAChRs (IC<sub>50</sub> = 94 nM). It is selective for α7 subunit-containing nAChRs over α6, β2, and β4 subunit-containing nAChRs (K<sub>i</sub>s = 4.52, 12, and 9.2 μM, respectively).<sup>2</sup> It selectively decreases vagus nerve stimulation-induced contractions of isolated guinea pig vagus nerve-stomach preparations (EC<sub>50</sub> = 49.4 μM) over isolated rat phrenic nerve-hemidiaphragm preparations (EC<sub>50</sub> = 486 μM). MG624 decreases protein levels of early growth response gene 1 (Egr-1) induced by nicotine in human microvascular endothelial cells of the lung (HMEC-Ls).<sup>3</sup> It also decreases proliferation of HMEC-Ls and inhibits angiogenesis *in vitro* and *ex vivo* as well as in an H69 small cell lung cancer (SCLC) mouse xenograft model concomitant with a reduction in tumor growth when administered in the diet at a dose of approximately 10 mg/kg per day.

### References

1. Gotti, C., Balestra, B., Morretti, M., *et al.* 4-Oxystilbene compounds are selective ligands for neuronal nicotinic αBungarotoxin receptors. *Br. J. Pharmacol.* **124**(6), 1197-1206 (1998).
2. Vailati, S., Moretti, M., Longhi, R., *et al.* Developmental expression of heteromeric nicotinic receptor subtypes in chick retina. *Mol. Pharmacol.* **63**(6), 1329-1337 (2003).
3. Brown, K.C., Lau, J.K., Dom, A.M., *et al.* MG624, an α7-nAChR antagonist, inhibits angiogenesis via the Egr-1/FGF2 pathway. *Angiogenesis* **15**(1), 99-114 (2012).

#### 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, 11/29/2022

#### 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