

# PRODUCT INFORMATION



## Dimethyl-W84 (dibromide)

Item No. 16927

**CAS Registry No.:** 402475-33-6  
**Formal Name:** N<sup>1</sup>,N<sup>6</sup>-bis[3-(1,3-dihydro-5-methyl-1,3-dioxo-2H-isoindol-2-yl)propyl]-N<sup>1</sup>,N<sup>1</sup>,N<sup>6</sup>,N<sup>6</sup>-tetramethyl-1,6-hexanediaminium, dibromide

**MF:** C<sub>34</sub>H<sub>48</sub>N<sub>4</sub>O<sub>4</sub> • 2Br

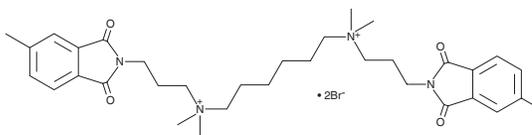
**FW:** 736.6

**Purity:** ≥98%

**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

Dimethyl-W84 (dibromide) is supplied as a solid. A stock solution may be made by dissolving the dimethyl-W84 (dibromide) in the solvent of choice. Dimethyl-W84 (dibromide) is soluble in organic solvents such as DMSO which should be purged with an inert gas. Dimethyl-W84 (dibromide) is also soluble in water. We do not recommend storing the aqueous solution for more than one day.

### Description

Dimethyl-W84 is a selective allosteric modulator of the M<sub>2</sub> muscarinic acetylcholine receptor.<sup>1,2</sup> It hinders the dissociation of the orthosteric antagonist N-methylscopolamine from the M<sub>2</sub> receptor with an EC<sub>50</sub> value of 3 nM.<sup>1,2</sup>

### References

1. Tränkle, C., Mies-Klomfass, E., Cid, M.H.B., et al. Identification of a [<sup>3</sup>H]Ligand for the common allosteric site of muscarinic acetylcholine M<sub>2</sub> receptors. *Mol. Pharmacol.* **54(1)**, 139-145 (1998).
2. Tränkle, C., Weyand, O., Voigtländer, U., et al. Interactions of orthosteric and allosteric ligands with [<sup>3</sup>H] dimethyl-W84 at the common allosteric site of muscarinic M<sub>2</sub> receptors. *Mol. Pharmacol.* **64(1)**, 180-190 (2003).

#### 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