

Dimethyl-W84 dibromide

M2 receptor modulator

Very potent allosteric modulator of muscarinic receptors, selective for M2-receptors. Retards the dissociation of the antagonist N-methylscopolamine. Affinity for the free receptor is about 10-fold higher than with W84 (Prod. No. ALX-550-399). $pK_A=7.08$, $p(\alpha K_A)=6.75$, $pEC_{50,diss}=6.77$.

Citations: 8

[View Online »](#)

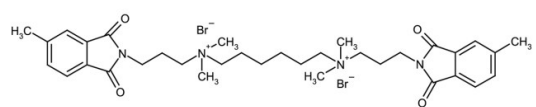
Ordering Information

[Order Online »](#)

ALX-550-400-M005	5mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Handling Avoid freeze/thaw cycles.

Long Term Storage +4°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name N,N'-Bis[3-(1,3-dihydro-5-methyl-1,3-dioxo-2H-isoindol-2-yl)propyl]-N,N,N',N'-tetramethyl-1,6-hexanediamminium dibromide

Appearance White to off-white solid.

CAS 402475-33-6

Couple Target Acetylcholine receptor

Couple Type Ligand

Formula $C_{34}H_{48}Br_2N_4O_4$

MW 736.6

Purity ≥98% (HPLC)

Solubility Soluble in DMSO, water or aqueous buffer.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)