

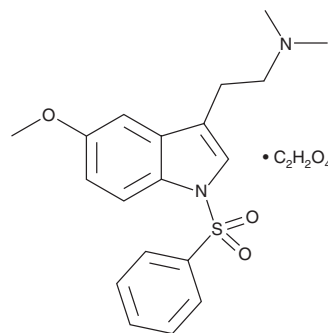
PRODUCT INFORMATION



MS 245 (oxalate)

Item No. 11936

CAS Registry No.: 275363-58-1
Formal Name: 5-methoxy-N,N-dimethyl-1-(phenylsulfonyl)-1H-indole-3-ethanamine, monoethanedioate
MF: C₁₉H₂₂N₂O₃S • C₂H₂O₄
FW: 448.5
Purity: ≥98%
UV/Vis.: λ_{max}: 212, 217, 259 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

MS 245 (oxalate) is supplied as a crystalline solid. A stock solution may be made by dissolving the MS 245 (oxalate) in the solvent of choice, which should be purged with an inert gas. MS 245 (oxalate) is soluble in the organic solvent DMSO.

Description

MS 245 is an antagonist of the serotonin (5-HT) receptor subtype 5-HT₆ (K_i = 2.3 nM for the human recombinant receptor).¹ It is selective for 5-HT₆ over 5-HT_{2A} and 5-HT_{2C} receptors (K_is = 130 and 23 nM, respectively), as well as 5-HT_{1A}, 5-HT_{1B}, 5-HT_{1E}, 5-HT₃, and 5-HT₇ receptors (K_is = 720, 9,200, 4,220, 2,390, and 600 nM, respectively). MS 245 decreases cAMP production induced by 5-HT in HEK293 cells expressing 5-HT₆ (pA₂ = 8.88 nM). It potentiates drug discrimination induced by (+)-amphetamine, but has no effect on cocaine or (-)-nicotine (Item No. 20887) drug discrimination in rats when administered at a dose of 5 mg/kg.^{2,3}

References

1. Tsai, Y., Dukat, M., Slassi, A., *et al.* N₁-(Benzenesulfonyl)tryptamines as novel 5-HT₆ antagonists. *Bioorg. Med. Chem. Lett.* **10(20)**, 2295-2299 (2000).
2. Pullagurla, M., Bondareva, T., Young, R., *et al.* Modulation of the stimulus effects of (+)amphetamine by the 5-HT₆ antagonist MS-245. *Pharmacol. Biochem. Behav.* **78(2)**, 263-268 (2004).
3. Young, R., Bondareva, T., Wesolowska, A., *et al.* Effect of the 5-HT₆ serotonin antagonist MS-245 on the actions of (-)nicotine. *Pharmacol. Biochem. Behav.* **85(1)**, 170-177 (2006).

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, 11/30/2022

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