

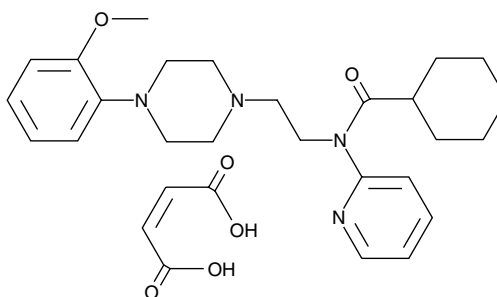
Product Information



WAY-100635 (maleate)

Item No. 14599

CAS Registry No.: 1092679-51-0
Formal Name: N-[2-[4-(2-methoxyphenyl)-1-piperazinyl]ethyl]-N-2-pyridinyl-cyclohexanecarboxamide, 2Z-butenedioate
MF: C₂₅H₃₄N₄O₂ • C₄H₄O₄
FW: 538.6
Purity: ≥95%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid
UV/Vis.: λ_{max}: 271 nm



Laboratory Procedures

For long term storage, we suggest that WAY-100635 (maleate) be stored as supplied at -20°C. It should be stable for at least two years.

WAY-100635 (maleate) is supplied as a crystalline solid. A stock solution may be made by dissolving the WAY-100635 (maleate) in the solvent of choice. WAY-100635 (maleate) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of WAY-100635 (maleate) in these solvents is approximately 10, 15, and 25 mg/ml, respectively.

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 WAY-100635 (maleate) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of WAY-100635 (maleate) in PBS, pH 7.2, is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

WAY-100635 is a potent, silent antagonist of serotonin 5-HT_{1A} receptors with an IC₅₀ value of 2.2 nM (K_i = 0.8 nM) for inhibiting 5-HT_{1A} receptors in rat hippocampal membranes.^{1,2} Because it displays 100-fold selectivity for 5-HT_{1A} over other 5-HT subtypes, WAY-100635 is classically used to examine the distribution and function of 5-HT_{1A} receptors.³ However, WAY-100635 has also been shown to exhibit agonist activity at dopamine D₄ receptors (K_d = 2.4 nM).⁴

References

1. Zhuang, Z.-P., Kung, M.-P., and Kung, H.F. Synthesis and evaluation of 4-(2'-methoxyphenyl)-1-[2'-[N-(2''-pyridinyl)-p-iodobenzamido]ethyl]piperazine (p-MPPI): A new iodinated 5-HT_{1A} ligand. *J. Med. Chem.* **37**(10), 1406-1407 (1994).
2. Mensonides-Harsema, M.M., Liao, Y., Böttcher, H., *et al.* Synthesis and *in vitro* and *in vivo* functional studies of ortho-substituted phenylpiperazine and N-substituted 4-N-(o-methoxyphenyl)aminopiperidine analogues of WAY100635. *J. Med. Chem.* **43**(3), 432-439 (2000).
3. Forster, E.A., Cliffe, I.A., Bill, D.J., *et al.* A pharmacological profile of the selective silent 5-HT_{1A} receptor antagonist, WAY-100635. *Eur. J. Pharmacol.* **281**(1), 81-88 (1995).
4. Chemel, B.R., Roth, B.L., Armbruster, B., *et al.* WAY-100635 is a potent dopamine D₄ receptor agonist. *Psychopharmacology (Berl)* **188**(2), 244-251 (2006).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/14599

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY. NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. 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

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery.**

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy located on our website and in our catalog.**

Copyright Cayman Chemical Company, 02/10/2015

Cayman Chemical

Mailing address
1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone
(800) 364-9897
(734) 971-3335

Fax
(734) 971-3640

E-Mail
custserv@caymanchem.com

Web
www.caymanchem.com