

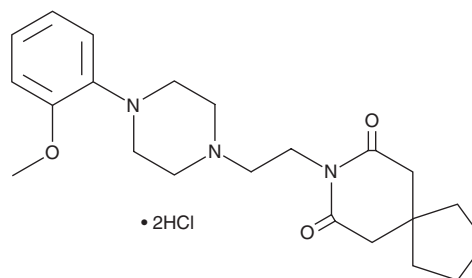
PRODUCT INFORMATION



BMY 7378

Item No. 21369

CAS Registry No.: 21102-95-4
Formal Name: 8-[2-[4-(2-methoxyphenyl)-1-piperazinyl]ethyl]-8-azaspiro[4.5]decane-7,9-dione, dihydrochloride
MF: C₂₂H₃₁N₃O₃ • 2HCl
FW: 458.4
Purity: ≥98%
UV/Vis.: λ_{max}: 208, 240, 279 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

BMY 7378 is supplied as a crystalline solid. A stock solution may be made by dissolving the BMY 7378 in the solvent of choice, which should be purged with an inert gas. BMY 7378 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of BMY 7378 in these solvents is approximately 0.2, 25, and 10 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 BMY 7378 can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of BMY 7378 in PBS (pH 7.2) is approximately 0.5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

BMY 7378 is an antagonist of the α_{2C}-adrenoceptor and α_{1D}-adrenoceptor with pK_i values of 6.54 and 8.2, respectively, that has been used in functional systems to assess the contribution of these receptors in vascular smooth muscle contraction.^{1,2} It also acts as a partial agonist at the serotonin 5-HT_{1A} receptor with a pK_i value of 8.3.¹

References

1. Goetz, A.S., King, H.K., Ward, S.D., *et al.* BMY 7378 is a selective antagonist of the D subtype of alpha 1-adrenoceptors. *Eur. J. Pharmacol.* **272(2-3)**, R5-R6 (1995).
2. Cleary, L., Murad, K., Bexis, S., *et al.* The alpha (1D)-adrenoceptor antagonist BMY 7378 is also an alpha (2C)-adrenoceptor antagonist. *Auton. Autacoid Pharmacol.* **25(4)**, 135-141 (2005).

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