

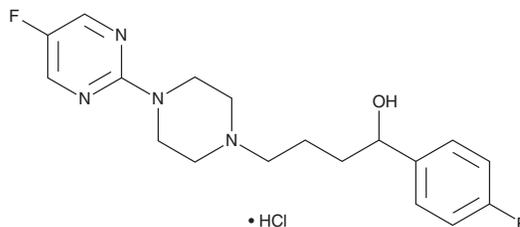
PRODUCT INFORMATION



BMY 14802 (hydrochloride)

Item No. 42264

CAS Registry No.: 105565-55-7
Formal Name: α -(4-fluorophenyl)-4-(5-fluoro-2-pyrimidinyl)-1-piperazinebutanol, monohydrochloride
Synonym: BMS 181100
MF: $C_{18}H_{22}F_2N_4O \cdot HCl$
FW: 384.9
Purity: $\geq 98\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

BMY 14802 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the BMY 14802 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. BMY 14802 (hydrochloride) is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in ethanol.

Description

BMY 14802 is an atypical antipsychotic.¹ It selectively binds to σ receptors over α_1 -adrenergic receptors (α_1 -ARs) and dopamine D_2 receptors (IC_{50} s = 112, 520, and 6,340 nM, respectively). *In vivo*, BMY 14802 inhibits the conditioned avoidance response and apomorphine-induced stereotypy in rats (ED_{50} s = 26.4 and 33 mg/kg, respectively).

Reference

1. Yevich, J.P., New, J.S., Lobeck, W.G., *et al.* Synthesis and biological characterization of α -(4-fluorophenyl)-4-(5-fluoro-2-pyrimidinyl)-1-piperazinebutanol and analogues as potential atypical antipsychotic agents. *J. Med. Chem.* **35**(24), 4516-4525 (1992).

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, 11/15/2024

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