

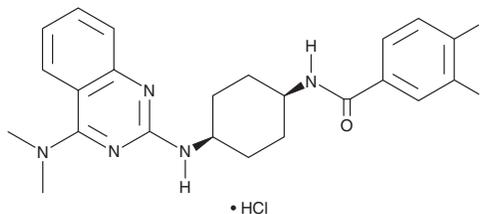
PRODUCT INFORMATION



ATC 0175

Item No. 42376

CAS Registry No.: 510733-97-8
Formal Name: N-[cis-4-[[4-(dimethylamino)-2-quinazoliny]amino]cyclohexyl]-3,4-difluoro-benzamide, monohydrochloride
MF: C₂₃H₂₅F₂N₅O • HCl
FW: 461.9
Purity: ≥95%
Supplied as: A 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

ATC 0175 is supplied as a solid. A stock solution may be made by dissolving the ATC 0175 in the solvent of choice, which should be purged with an inert gas. ATC 0175 is slightly soluble (0.1-1 mg/ml) in ethanol and DMSO.

Description

ATC 0175 is a melanocortin receptor 1 (MC1R) antagonist (IC₅₀ = 7.23 nM).¹ It is selective for MC1R over MC2R (IC₅₀ = >10,000 nM) but does not antagonize the serotonin (5-HT) receptor subtypes 5-HT_{2B} and 5-HT_{1A} (IC₅₀s = 9.66 and 16.9 nM, respectively). ATC 0175 (3 nM) inhibits [³⁵S]GTPγS binding and calcium influx induced by melanin-concentrating hormone (MCH; Item No. 24462) in HEK293 cells expressing human MC1R. *In vivo*, ATC 0175 (1, 3, and 10 mg/kg) reduces stress-induced decreases in the time spent in the open arms of the elevated plus maze in mice. It also decreases immobility time in the forced swim test in mice.

Reference

1. Chaki, S., Yamaguchi, J.-i., Yamada, H., *et al.* ATC0175: An orally active melanin-concentrating hormone receptor 1 antagonist for the potential treatment of depression and anxiety. *CNS Drug Rev.* **11(4)**, 341-152 (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, 12/05/2024

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