

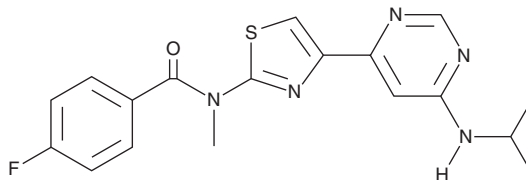
PRODUCT INFORMATION



FITM

Item No. 29219

CAS Registry No.: 932737-65-0
Formal Name: 4-fluoro-N-methyl-N-[4-[6-[(1-methylethyl)amino]-4-pyrimidinyl]-2-thiazolyl]-benzamide
MF: C₁₈H₁₈FN₅OS
FW: 371.4
Purity: ≥98%
UV/Vis.: λ_{max}: 226, 260 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

FITM is supplied as a crystalline solid. A stock solution may be made by dissolving the FITM in the solvent of choice, which should be purged with an inert gas. FITM is soluble in organic solvents such as ethanol and DMSO. The solubility of FITM in these solvents is approximately 20 and 100 mM, respectively.

Description

FITM is an antagonist of metabotropic glutamate receptor 1 (mGluR1; IC₅₀ = 5.1 nM for the human receptor).¹ It is selective for mGluR1 over mGluR2, mGluR5, and mGluR8 (IC₅₀s = >10, 7, and >10 μM, respectively). FITM (0.3 mg/kg) decreases methamphetamine-induced hyperlocomotion in mice and reverses disruptions in prepulse inhibition induced by methamphetamine and ketamine in rats when administered at a dose of 1 mg/kg.

Reference

1. Satoh, A., Nagatomi, Y., Hirata, Y., *et al.* Discovery and in vitro and in vivo profiles of 4-fluoro-N-[4-[6-(isopropylamino)pyrimidin-4-yl]-1,3-thiazol-2-yl]-N-methylbenzamide as novel class of an orally active metabotropic glutamate receptor 1 (mGluR1) antagonist. *Bioorg. Med. Chem. Lett.* **19(18)**, 5464-5468 (2009).

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, 11/08/2022

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