

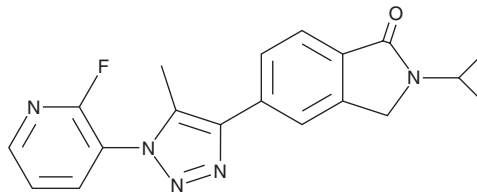
PRODUCT INFORMATION



CFMTI

Item No. 37546

CAS Registry No.: 864864-17-5
Formal Name: 2-cyclopropyl-5-[1-(2-fluoro-3-pyridinyl)-5-methyl-1H-1,2,3-triazol-4-yl]-2,3-dihydro-1H-isoindol-1-one
MF: C₁₉H₁₆FN₅O
FW: 349.4
Purity: ≥90%
UV/Vis.: λ_{max}: 272 nm
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

CFMTI is supplied as a solid. A stock solution may be made by dissolving the CFMTI in the solvent of choice, which should be purged with an inert gas. CFMTI is soluble in organic solvents such as chloroform, DMSO, and dimethyl formamide (DMF). The solubility of CFMTI in chloroform and DMF is approximately 1 and 10 mg/ml, respectively. CFMTI is slightly soluble in DMSO.

Description

CFMTI is a metabotropic glutamate receptor 1 (mGluR1) allosteric antagonist (IC₅₀ = 2.6 nM in CHO cells expressing the human receptor).^{1,2} It is selective for mGluR1 over mGluR5 by greater than 2,000-fold, as well as over mGluR2, -4, -6, -7, or -8 at 10 μM in CHO cells expressing the human receptors.² CFMTI reduces methamphetamine-induced hyperlocomotion in mice when administered orally at doses ranging from 1 to 10 mg/kg.

References

1. Ito, S., Hirata, Y., Nagatomi, Y., *et al.* Discovery and biological profile of isoindolinone derivatives as novel metabotropic glutamate receptor 1 antagonists: A potential treatment for psychotic disorders. *Bioorg. Med. Chem. Lett.* **19(18)**, 5310-5313 (2009).
2. Satow, A., Suzuki, G., Maehara, S., *et al.* Unique antipsychotic activities of the selective metabotropic glutamate receptor 1 allosteric antagonist 2-cyclopropyl-5-[1-(2-fluoro-3-pyridinyl)-5-methyl-1H-1,2,3-triazol-4-yl]-2,3-dihydro-1H-isoindol-1-one. *J. Pharmacol. Exp. Ther.* **330(1)**, 179-190 (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, 02/27/2023

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