

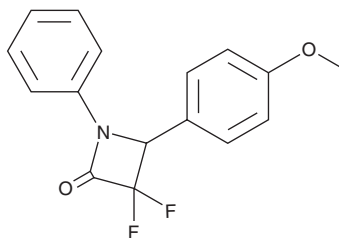
PRODUCT INFORMATION



NCRW0005-F05

Item No. 43215

CAS Registry No.: 342779-66-2
Formal Name: 3,3-difluoro-4-(4-methoxyphenyl)-1-phenyl-2-azetidinone
MF: C₁₆H₁₃F₂NO₂
FW: 289.3
Purity: ≥98%
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

NCRW0005-F05 is supplied as a solid. A stock solution may be made by dissolving the NCRW0005-F05 in the solvent of choice, which should be purged with an inert gas. NCRW0005-F05 is soluble (≥10 mg/ml) in DMSO. NCRW0005-F05 is sparingly soluble (1-10 mg/ml) in ethanol.

Description

NCRW0005-F05 is an antagonist of G protein-coupled receptor 139 (GPR139; IC₅₀ = 0.21 μM).¹

Reference

1. Wang, J., Zhu, L., Liu, Q., *et al.* High-throughput screening of antagonists for the orphan G-protein coupled receptor GPR139. *Acta Pharmacol. Sin.* **36**(7), 874-878 (2015).

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, 04/18/2025

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