

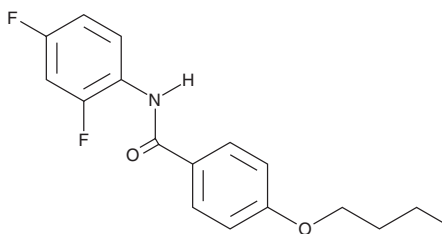
PRODUCT INFORMATION



VU0357121

Item No. 22199

CAS Registry No.: 433967-28-3
Formal Name: 4-butoxy-N-(2,4-difluorophenyl)-benzamide
MF: C₁₇H₁₇F₂NO₂
FW: 305.3
Purity: ≥98%
UV/Vis.: λ_{max}: 262 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

VU0357121 is supplied as a crystalline solid. A stock solution may be made by dissolving the VU0357121 in the solvent of choice, which should be purged with an inert gas. VU0357121 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of VU0357121 in these solvents is approximately 15 and 20 mg/ml, respectively.

VU0357121 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, VU0357121 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. VU0357121 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

VU0357121 is a positive allosteric modulator of metabotropic glutamate receptor 5 (mGluR5) that has an EC₅₀ value of 33 nM for the rat receptor expressed in HEK293 cells in a calcium flux assay.¹ It is selective for mGluR5 and does not modulate responses at mGluR2-4 or mGluR6-8 but does inhibit mGluR1.

Reference

1. Hammond, A.X., Rodriguez, A.L., Townsend, S.D., *et al.* Discovery of a novel chemical class of mGlu₅ allosteric ligands with distinct modes of pharmacology. *ACS Chem. Neurosci.* **1**(10), 702-716 (2010).

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, 11/15/2022

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