

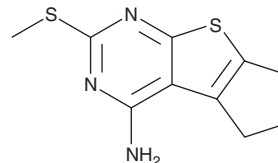
PRODUCT INFORMATION



VU0029251

Item No. 21987

CAS Registry No.: 330819-85-7
Formal Name: 6,7-dihydro-2-(methylthio)-5H-cyclopenta[4,5]thieno[2,3-d]pyrimidin-4-amine
MF: C₁₀H₁₁N₃S₂
FW: 237.3
Purity: ≥98%
UV/Vis.: λ_{max}: 241, 287 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

VU0029251 is supplied as a crystalline solid. A stock solution may be made by dissolving the VU0029251 in the solvent of choice, which should be purged with an inert gas. VU0029251 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of VU0029251 in these solvents is approximately 20 and 30 mg/ml, respectively.

VU0029251 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, VU0029251 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. VU0029251 has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

VU0029251 is a partial antagonist of metabotropic glutamate receptor 5 (mGluR5; K_i = 1.07 μM).¹ It inhibits calcium mobilization induced by glutamate in HEK293 cell membranes expressing rat mGluR5 (IC₅₀ = 1.7 μM).

Reference

1. Rodriguez, A.L., Grier, M.D., Jones, C.K., *et al.* Discovery of novel allosteric modulators of metabotropic glutamate receptor subtype 5 reveals chemical and functional diversity and in vivo activity in rat behavioral models of anxiolytic and antipsychotic activity. *Mol. Pharmacol.* **78**(6), 1105-1123 (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, 12/16/2022

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