

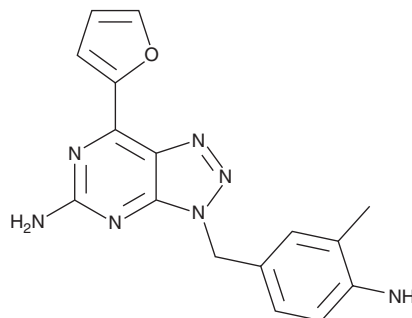
PRODUCT INFORMATION



Vipadenant

Item No. 20239

CAS Registry No.: 442908-10-3
Formal Name: 3-[(4-amino-3-methylphenyl)methyl]-7-(2-furanyl)-3H-1,2,3-triazolo[4,5-d]pyrimidin-5-amine
Synonym: BIIB-014
MF: C₁₆H₁₅N₇O
FW: 321.3
Purity: ≥98%
UV/Vis.: λ_{max}: 231, 301, 349 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

Vipadenant is supplied as a crystalline solid. A stock solution may be made by dissolving the vipadenant in the solvent of choice, which should be purged with an inert gas. Vipadenant is soluble in the organic solvent DMSO at a concentration of approximately 43 mg/ml.

Description

Vipadenant is a selective adenosine 2A (A_{2A}) receptor antagonist with K_i values of 1.3, 68, 63, and 1,005 nM for A_{2A}, A₁, A_{2B}, and A₃ receptors, respectively, in a radioligand binding assay.¹ It acts as a functional antagonist of human recombinant A_{2A} (K_A = 0.58 nM) in a calcium mobilization-based FLIPR assay. Vipadenant also reverses haloperidol-induced hypolocomotion in mice and rats.

Reference

1. Gillespie, R.J., Bamford, S.J., Botting, R., *et al.* Antagonists of the human A_{2A} adenosine receptor. 4. Design, synthesis, and preclinical evaluation of 7-aryltriazolo[4,5-d]pyrimidines. *J. Med. Chem.* **52**(1), 33-47 (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, 12/19/2022

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