

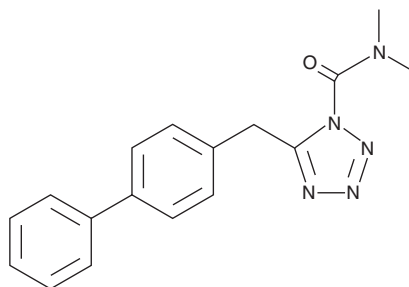
PRODUCT INFORMATION



LY2183240

Item No. 10008663

CAS Registry No.: 874902-19-9
Formal Name: 5-([1,1'-biphenyl]-4-ylmethyl)-N,N-dimethyl-1H-tetrazole-1-carboxamide
MF: C₁₇H₁₇N₅O
FW: 307.4
Purity: ≥98%
UV/Vis.: λ_{max}: 253 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

LY2183240 is supplied as a crystalline solid. A stock solution may be made by dissolving the LY2183240 in the solvent of choice. LY2183240 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of LY2183240 in these solvents is approximately 10, 2, and 3 mg/ml, respectively.

LY2183240 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, LY2183240 should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. LY2183240 has a solubility of approximately 0.1 mg/ml in a 1:5 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

LY2183240 is a potent, competitive inhibitor of anandamide uptake (IC₅₀ = 270 pM; K_i = 540 pM) and hydrolysis.¹⁻³ It increases anandamide levels in rat cerebellum (ED₅₀ = 1.37 mg/kg) and displays dose-dependent efficacy (3-30 mg/kg) in several rodent models of persistent pain.¹

References

1. Moore, S.A., Nomikos, G.G., Dickason-Chesterfield, A.K., *et al.* Identification of a high-affinity binding site involved in the transport of endocannabinoids. *Proc. Natl. Acad. Sci. USA* **102**(49), 17852-17857 (2005).
2. Ortar, G., Cascio, M.G., Moriello, A.S., *et al.* Carbamoyl tetrazoles as inhibitors of endocannabinoid inactivation: A critical re-visitation. *Eur. J. Med. Chem.* **43**, 62-72 (2008).
3. Di Marzo, V. Targeting the endocannabinoid system: To enhance or reduce? *Nat. Rev. Drug Discov.* **7**, 438-455 (2008).

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, 07/12/2017

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