

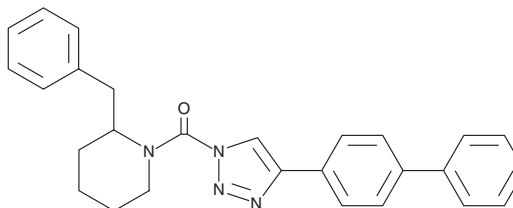
PRODUCT INFORMATION



KT109

Item No. 18933

CAS Registry No.: 1402612-55-8
Formal Name: (4-[1,1'-biphenyl]-4-yl-1H-1,2,3-triazol-1-yl)[2-(phenylmethyl)-1-piperidiny]-methanone
MF: C₂₇H₂₆N₄O
FW: 422.5
Purity: ≥98%
UV/Vis.: λ_{max}: 278 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

KT109 is supplied as a crystalline solid. A stock solution may be made by dissolving the KT109 in the solvent of choice. KT109 is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of KT109 in these solvents is approximately 10 mg/ml.

KT109 is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

In humans, two forms of diacylglycerol lipase, DAGLα and DAGLβ, generate the endocannabinoid 2-arachidonoyl glycerol (2-AG; Item No. 62160) by attacking DAG at the *sn*-1 position. KT109 is a selective inhibitor of DAGLβ with an IC₅₀ value of 42 nM.¹ It demonstrates ~60-fold selectivity for DAGLβ over DAGLα, and negligible activity against other key enzymes involved in endocannabinoid signaling, including FAAH, MAGL, and ABHD11.¹ KT109 has been shown to disrupt the lipid network involved in macrophage inflammatory responses, lowering 2-AG, as well as arachidonic acid (Item No. 90010) and eicosanoids, in mouse peritoneal macrophages.¹

Reference

1. Hsu, K.-L., Tsuboi, K., Adibekian, A., *et al.* DAGLβ inhibition perturbs a lipid network involved in macrophage inflammatory responses. *Nat. Chem. Biol.* **8**(12), 999-1007 (2012).

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, 06/12/2018

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