

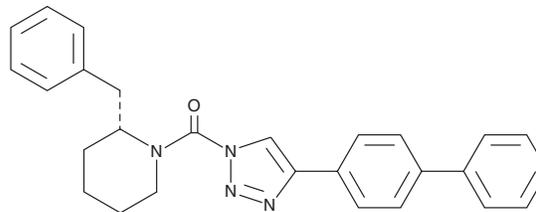
PRODUCT INFORMATION



(R)-KT109

Item No. 25682

CAS Registry No.: 2055172-60-4
Formal Name: (4-[1,1'-biphenyl]-4-yl-1H-1,2,3-triazol-1-yl)
[(2R)-2-(phenylmethyl)-1-piperidinyl]-methanone
MF: C₂₇H₂₆N₄O
FW: 422.5
Purity: ≥98%
UV/Vis.: λ_{max}: 279 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

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

Description

(R)-KT109 is the (R) isomer of the diacylglycerol lipase β (DAGL β) inhibitor KT109 (Item No. 18933).¹ (R)-KT109 is an inhibitor of DAGL β (IC₅₀ = 0.79 nM) and of DAGL α -mediated hydrolysis of 1-stearoyl-2-arachidonoyl-*sn*-glycerol (Item No. 10008650; IC₅₀ = 25.12 nM). It also inhibits α/β -hydrolase domain-containing protein 6 (ABHD6) with an IC₅₀ value of 2.51 nM. (R)-KT109 is more potent at DAGL β , DAGL α , and ABHD6 than (S)-KT109 (Item No. 25683).

Reference

- Deng, H., Kooijman, S., van den Nieuwendijk, A.M.C.H., *et al.* Triazole ureas act as diacylglycerol lipase inhibitors and prevent fasting-induced refeeding. *J. Med. Chem.* **60(1)**, 428-440 (2017).

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, 02/16/2024

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