

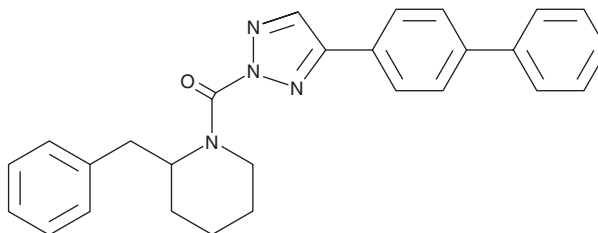
# PRODUCT INFORMATION



## KT109 N2 Regioisomer

Item No. 26127

**CAS Registry No.:** 2749638-90-0  
**Formal Name:** (4-[1,1'-biphenyl]-4-yl-2H-1,2,3-triazol-2-yl)[2-(phenylmethyl)-1-piperidiny]-methanone  
**MF:** C<sub>27</sub>H<sub>26</sub>N<sub>4</sub>O  
**FW:** 422.5  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 284 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

KT109 N2 regioisomer is supplied as a crystalline solid. A stock solution may be made by dissolving the KT109 N2 regioisomer in the solvent of choice, which should be purged with an inert gas. KT109 N2 regioisomer is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of KT109 N2 regioisomer in these solvents is approximately 10 mg/ml.

### Description

KT109 N2 regioisomer is an N2-carbamoylated regioisomer of the diacylglycerol lipase β (DAGLβ) inhibitor KT109 (Item No. 18933). It is intended for use as an internal standard for the quantification of KT109.

#### 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, 02/06/2025

#### 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](http://WWW.CAYMANCHEM.COM)