

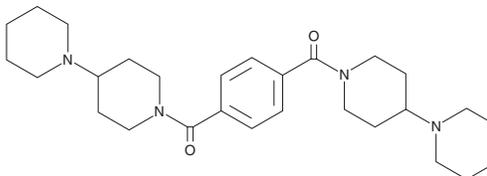
PRODUCT INFORMATION



UNC1079

Item No. 20566

CAS Registry No.: 1418741-86-2
Formal Name: 1,1'-(1,4-phenylene)bis[1-[1,4'-bipiperidin]-1'-yl]-methanone
MF: C₂₈H₄₂N₄O₂
FW: 466.7
Purity: ≥95%
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

UNC1079 is supplied as a crystalline solid. A stock solution may be made by dissolving the UNC1079 in the solvent of choice, which should be purged with an inert gas. UNC1079 is soluble in the organic solvent ethanol at concentration of approximately 1 mg/ml.

Description

UNC1079 is an analog of UNC1215 (Item No. 13968), the selective L3MBTL3 domain inhibitor. UNC1079 is a 1,000-fold weaker than UNC1215 as a L3MBTL3 domain inhibitor.¹ This compound is intended for use as a negative control in cellular studies.¹ See the Structural Genomics Consortium (SGC) website for more information.

Reference

1. James, L.I., Barsyte-Lovejoy, D., Zhong, N., *et al.* Discovery of a chemical probe for the L3MBTL3 methyllysine reader domain. *Nat. Chem. Biol.* **9**(3), 184-191 (2013).

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, 12/16/2022

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