

PRODUCT INFORMATION



BI-99179

Item No. 41307

CAS Registry No.: 1291779-76-4

Formal Name: N-[4-(2-benzoxazolyl)phenyl]-N-methyl-3S-[(1R-oxopropyl)amino]-cyclopentanecarboxamide

MF: C₂₃H₂₅N₃O₃

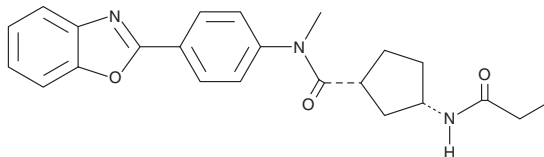
FW: 391.5

Purity: ≥98%

Supplied as: A 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

BI-99179 is supplied as a solid. A stock solution may be made by dissolving the BI-99179 in the solvent of choice, which should be purged with an inert gas. BI-99179 is slightly soluble (0.1-1 mg/ml) in acetonitrile and sparingly soluble (1-10 mg/ml) in DMSO.

Description

BI-99179 is an inhibitor of fatty acid synthase type I (FAS-I; IC₅₀ = 79 nM), a multifunctional enzyme that catalyzes the conversion of acetyl-coenzyme A (acetyl-CoA; Item No. 16160) and malonyl-CoA (Item No. 16455) to long-chain fatty acids.¹ It is selective for FAS-I over 30 other enzymes, receptors, and ion channels at 10 μM. BI-99179 inhibits the incorporation of ¹⁴C-acetate in mouse N-42 hypothalamus cells (IC₅₀ = 570 nM). It reduces the proliferation of GaMG human glioma cells when used at a concentration of 10 μM.²

References

1. Kley, J.T., Mack, J.B.C., Hamilton, B., *et al.* Discovery of BI 99179, a potent and selective inhibitor of type I fatty acid synthase with central exposure. *Bioorg. Med. Chem. Lett.* **21(19)**, 5924-5927 (2011).
2. Singha, P.K., Mäklin, K., Vihavainen, T., *et al.* Evaluation of FASN inhibitors by a versatile toolkit reveals differences in pharmacology between human and rodent FASN preparations and in antiproliferative efficacy *in vitro* vs. *in situ* in human cancer cells. *Eur. J. Pharm. Sci.* **149**, 105321 (2020).

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, 07/15/2024

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