

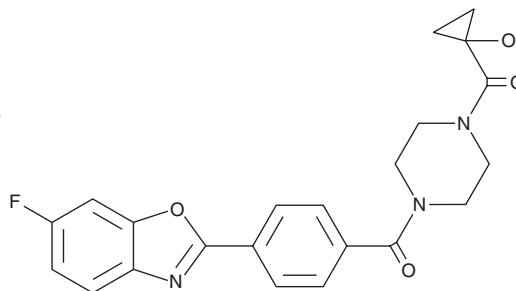
PRODUCT INFORMATION



FT113

Item No. 37545

CAS Registry No.: 1630808-89-7
Formal Name: [4-[4-(6-fluoro-2-benzoxazolyl)benzoyl]-1-piperazinyl](1-hydroxycyclopropyl)-methanone
MF: C₂₂H₂₀FN₃O₄
FW: 409.4
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

FT113 is supplied as a solid. A stock solution may be made by dissolving the FT113 in the solvent of choice, which should be purged with an inert gas. FT113 is soluble in ethanol and methanol.

Description

FT113 is an inhibitor of fatty acid synthase (FASN; IC₅₀ = 213 nM).¹ It inhibits the proliferation of PC3 human prostate cancer and MV4-11 human acute myeloid leukemia cells (IC₅₀s = 47 and 26 nM, respectively) and inhibits FASN activity in a ¹⁴C-acetate incorporation assay using BT474 human breast cancer cells (IC₅₀ = 90 nM). FT113 (25 and 50 mg/kg) reduces tumor growth and increases intratumoral levels of the FASN substrate malonyl-CoA in an MV4-11 mouse xenograft model. It also inhibits infection of HEK293T cells expressing human angiotensin-converting enzyme 2 (HEK293T-hACE2 cells) with an infectious-clone-derived mNeonGreen severe acute respiratory syndrome coronavirus 2 reporter virus (SARS-CoV-2-mNG) with an EC₅₀ value of 17 nM.²

References

1. Martin, M.W., Lancia, D.R., Jr., Li, H., *et al.* Discovery and optimization of novel piperazines as potent inhibitors of fatty acid synthase (FASN). *Bioorg. Med. Chem. Lett.* **29(8)**, 1001-1006 (2019).
2. Chu, J., Xing, C., Du, Y., *et al.* Pharmacological inhibition of fatty acid synthesis blocks SARS-CoV-2 replication. *Nat. Metab.* **3(11)**, 1466-1475 (2021).

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, 11/02/2023

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