

PRODUCT INFORMATION



TVB-2640

Item No. 35703

CAS Registry No.: 1399177-37-7
Formal Name: 4-[1-[4-cyclobutyl-2-methyl-5-(3-methyl-1H-1,2,4-triazol-5-yl)benzoyl]-4-piperidinyl]-benzonitrile

Synonyms: Denifanstat,
Fatty Acid Synthase Inhibitor 2,
FASN-IN-2

MF: C₂₇H₂₉N₅O

FW: 439.6

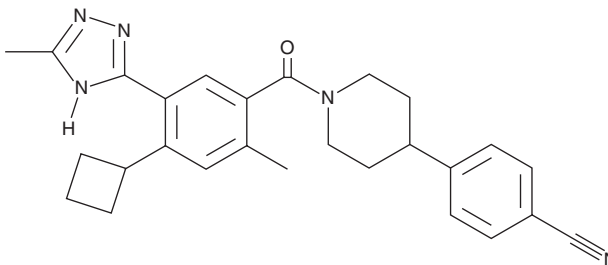
Purity: ≥98%

UV/Vis.: λ_{max}: 231 nm

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

TVB-2640 is supplied as a solid. A stock solution may be made by dissolving the TVB-2640 in the solvent of choice, which should be purged with an inert gas. TVB-2640 is soluble in DMSO.

Description

TVB-2640 is an inhibitor of fatty acid synthase (FASN; IC₅₀ = 0.5 μM).¹ It inhibits infection of an infectious-clone-derived severe acute respiratory syndrome coronavirus 2 (SARS-CoV-2) reporter virus in HEK293T cells expressing human angiotensin-converting enzyme 2 (ACE2; EC₅₀ = 4 nM).²

References

1. Syed-Abdul, M.M., Parks, E.J., Gaballah, A.H., *et al.* Fatty acid synthase inhibitor TVB-2640 reduces hepatic *de novo* lipogenesis in males with metabolic abnormalities. *Hepatology* **72(1)**, 103-118 (2020).
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, 10/24/2022

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