

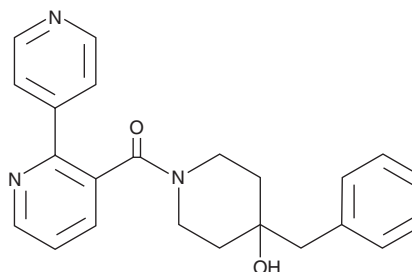
PRODUCT INFORMATION



Soticlestat

Item No. 34928

CAS Registry No.: 1429505-03-2
Formal Name: [2,4'-bipyridin]-3-yl[4-hydroxy-4-(phenylmethyl)-1-piperidinyl]-methanone
Synonyms: OV935, TAK-935
MF: C₂₃H₂₃N₃O₂
FW: 373.4
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

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

Description

Soticlestat is an inhibitor of cholesterol 24-hydroxylase/CYP46A1 ($IC_{50} = 0.0074 \mu\text{M}$).¹ It is selective for cholesterol 24-hydroxylase/CYP46A1 over the cytochrome P450 (CYP) isoforms CYP2C8, -2C9, -2D6, -3A4, -1A2, and 2C19 (IC_{50} s = 62, 19, >100, 66, >100, and 14 μM , respectively). Soticlestat (0.02% in the diet) reduces seizure frequency and severity and improves survival in a model of Dravet syndrome induced by hyperthermia in mice with heterozygous deletion of *Scn1a* (*Scn1a*^{+/-}), the gene encoding voltage-gated sodium channel 1.1 ($Na_v1.1$).²

References

1. Koike, T., Yoshikawa, M., Ando, H.K., *et al.* Discovery of soticlestat, a potent and selective inhibitor for cholesterol 24-hydroxylase (CH24H). *J. Med. Chem.* **64**(16), 12228-12244 (2021).
2. Hawkins, N.A., Jurado, M., Thaxton, T.T., *et al.* Soticlestat, a novel cholesterol 24-hydroxylase inhibitor, reduces seizures and premature death in Dravet syndrome mice. *Epilepsia* **62**(11), 2845-2857 (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, 12/14/2022

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