

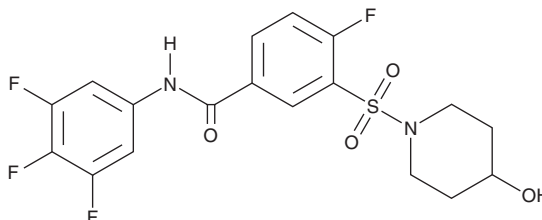
PRODUCT INFORMATION



NVR 3-778

Item No. 42725

CAS Registry No.: 1445790-55-5
Formal Name: 4-fluoro-3-[(4-hydroxy-1-piperidinyl)sulfonyl]-N-(3,4,5-trifluorophenyl)-benzamide
Synonyms: AL 3778, JNJ-63595948
MF: C₁₈H₁₆F₄N₂O₄S
FW: 432.4
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

NVR 3-778 is supplied as a solid. A stock solution may be made by dissolving the NVR 3-778 in the solvent of choice, which should be purged with an inert gas. NVR 3-778 is soluble (≥10 mg/ml) in ethanol and DMSO.

Description

NVR 3-778 is a hepatitis B virus (HBV) capsid assembly modulator.¹ It inhibits viral replication of HBV genotypes A-H in HepG2 2.2.15 cells (mean EC₅₀s = 0.2-0.58 μM). NVR 3-778 inhibits the production of HBV surface antigen (HBsAg), HBV e antigen (HBeAg), and HBV DNA in a primary human hepatocyte model of *de novo* HBV infection (EC₅₀s = 4.8, 4.5, and 0.81 μM, respectively). *In vivo*, NVR 3-778 (40 mg/kg twice per day) reduces serum levels of HBV DNA in HBV-infected mice with humanized livers.²

References

1. Lam, A.M., Espiritu, C., Vogel, R., *et al.* Preclinical characterization of NVR 3-778, a first-in-class capsid assembly modulator against hepatitis B virus. *Antimicrob. Agents Chemother.* **63**(1), e01734-18 (2018).
2. Klumpp, K., Shimada, T., Allweiss, L., *et al.* Efficacy of NVR 3-778, alone and In combination With pegylated interferon, vs entecavir in uPA/SCID mice With humanized livers and HBV infection. *Gastroenterology* **154**(3), 652-662.e8 (2018).

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, 02/12/2025

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