

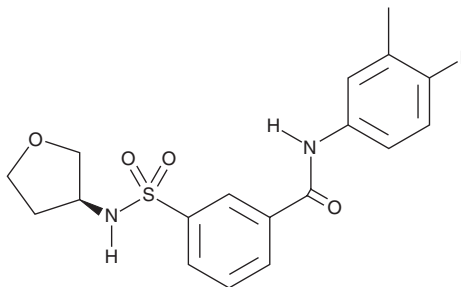
PRODUCT INFORMATION



JNJ-632

Item No. 26325

CAS Registry No.: 1572510-42-9
Formal Name: N-(4-fluoro-3-methylphenyl)-3-[[[(3S)-tetrahydro-3-furanyl]amino]sulfonyl]-benzamide
MF: C₁₈H₁₉FN₂O₄S
FW: 378.4
Purity: ≥98%
Supplied as: A crystalline 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

JNJ-632 is supplied as a crystalline solid. A stock solution may be made by dissolving the JNJ-632 in the solvent of choice, which should be purged with an inert gas. JNJ-632 is soluble in the organic solvent DMSO at a concentration of approximately 50 mg/ml.

Description

JNJ-632 is a modulator of hepatitis B virus (HBV) capsid assembly.^{1,2} It reduces viral DNA load in HBV-infected HepG2.2.15 cells (EC₅₀ = 121 nM) without affecting cell growth.² It also reduces viral DNA load in primary human hepatocytes infected with the HBV genotypes Ae, Bj, C, and D (EC₅₀s = 101, 240, 119, and 200 nM, respectively). JNJ-632 prevents formation of covalently closed circular DNA in HBV-infected primary human hepatocytes in a concentration-dependent manner when administered in combination with the HBV viral inoculum. *In vivo*, JNJ-632 (200 mg/kg per day) reduces plasma HBV DNA content in HBV-infected humanized mice.¹

References

1. Vandyck, K., Rombouts, G., Stoops, B., *et al.* Synthesis and evaluation of N-phenyl-3-sulfamoyl-benzamide derivatives as capsid assembly modulators inhibiting hepatitis B virus (HBV). *J. Med. Chem.* **61**(14), 6247-6260 (2018).
2. Berke, J.M., Dehertogh, P., Vergauwen, K., *et al.* Capsid assembly modulators have a dual mechanism of action in primary human hepatocytes infected with hepatitis B virus. *Antimicrob. Agents Chemother.* **61**(8), pii: e00560-176 (2017).

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, 12/20/2022

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