

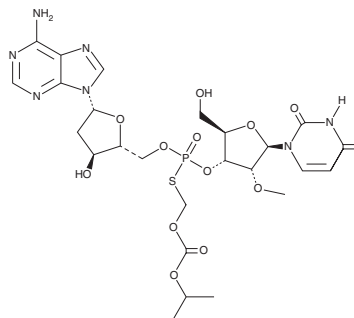
PRODUCT INFORMATION



SB-9200

Item No. 42317

CAS Registry No.: 942123-43-5
Formal Name: 2'-O-methyl-P(S)-[[[(1-methylethoxy)carbonyl]oxy]methyl]-P-thiouridylyl-(3'→5')-2'-deoxy-adenosine
Synonyms: GS-9992, Inarigivir Soproxil
MF: C₂₅H₃₄N₇O₁₃PS
FW: 703.6
Purity: ≥95%
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

SB-9200 is supplied as a solid. A stock solution may be made by dissolving the SB-9200 in the solvent of choice, which should be purged with an inert gas. SB-9200 is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of SB-9200 can be prepared by directly dissolving the solid in aqueous buffers. SB-9200 is sparingly soluble (1-10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

SB-9200 is an antiviral agent and a prodrug form of SB-9000.¹ It reduces replication of wild-type, lamivudine-resistant, or adefovir dipivoxil-resistant hepatitis B virus (HBV) in HepG2/2.15 cells (EC₅₀s = 2.5, 2.1, and 2.8 μM, respectively). SB-9200 inhibits replication in genotype 1a or 1b hepatitis C virus (HCV) replicon cells (EC₅₀s = 2.2 and 1 μM).² It also decreases viability, migration and invasion, and the levels of STAT3 and phosphorylated STAT3 and increases the levels of DEAD-box helicase 58 (DDX58) in HCT116 and SW620 colon cancer cells when used at a concentration of 80 μM.³ SB-9200 (100 mg/kg) reduces hepatic levels of HBV DNA in HBV-infected mice.¹

References

1. Coughlin, J.E., Padmanabhan, S., Zhang, G., *et al.* Orally bioavailable anti-HBV dinucleotide acyloxyalkyl prodrugs. *Bioorg. Med. Chem. Lett.* **20**(5), 1783-1786 (2010).
2. Jones, M., Cunningham, M.E., Wing, P., *et al.* SB 9200, a novel agonist of innate immunity, shows potent antiviral activity against resistant HCV variants. *J. Med. Virol.* **89**(9), 1620-1628 (2017).
3. Deng, Y., Fu, H., Han, X., *et al.* Activation of DDX58/RIG-I suppresses the growth of tumor cells by inhibiting STAT3/CSE signaling in colon cancer. *Int. J. Oncol.* **61**(4), 120 (2022).

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, 12/04/2024

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