

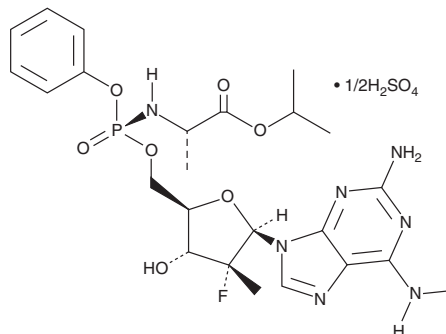
PRODUCT INFORMATION



AT-511 (hemisulfate)

Item No. 33870

CAS Registry No.: 2241337-84-6
Formal Name: N-[[P(S),2'R]-2-amino-2'-deoxy-2'-fluoro-N,2'-dimethyl-P-phenyl-5'-adenylyl]-L-alanine, 1-methylethyl ester, hemisulfate
Synonyms: AT-527 (hemisulfate), RO7496998
MF: C₂₄H₃₃FN₇O₇P • 1/2H₂SO₄
FW: 630.6
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

AT-511 (hemisulfate) is supplied as a solid. A stock solution may be made by dissolving the AT-511 (hemisulfate) in the solvent of choice, which should be purged with an inert gas. AT-511 (hemisulfate) is soluble in the organic solvent DMSO.

Description

AT-511 is a prodrug form of the antiviral nucleoside analog AT-9010.¹ It reduces the infectious virus yield in cells infected with Middle East respiratory syndrome coronavirus (MERS-CoV), severe acute respiratory syndrome CoV (SARS-CoV), SARS-CoV-2, human CoV 229E (HCoV-229E), and HCoV-OC43 (EC₅₀s = 37, 0.47, 2.8, 1.2, and 0.5 μM, respectively).

Reference

1. Good, S.S., Westover, J., Jung, K.H., *et al.* AT-527, a double prodrug of a guanosine nucleotide analog, is a potent inhibitor of SARS-CoV-2 *in vitro* and a promising oral antiviral for treatment of COVID-19. *Antimicrob. Agents Chemother.* **65**(4), e02479-20 (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, 12/01/2023

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