

PRODUCT INFORMATION

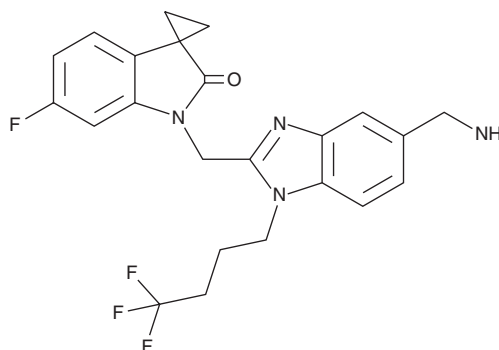


Sisunatovir

Item No. 42124

CAS Registry No.: 1903763-82-5
Formal Name: 1'-[[5-(aminomethyl)-1-(4,4,4-trifluorobutyl)-1H-benzimidazol-2-yl]methyl]-6'-fluoro-spiro[cyclopropane-1,3'-[3H]indol]-2'(1'H)-one

Synonym: RV521
MF: C₂₃H₂₂F₄N₄O
FW: 446.4
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

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

Description

Sisunatovir is an inhibitor of respiratory syncytial virus (RSV) fusion (IC₅₀ = 0.9 nM).¹ It reduces RSV A2 replication in a plaque reduction assay *in vitro* (EC₅₀ = 1.3 nM). Sisunatovir (1, 10, or 50 mg/kg) decreases lung viral titers in a mouse model of RSV infection.

Reference

1. Cockerill, G.S., Angell, R.M., Bedernjak, A., *et al.* Discovery of sisunatovir (RV521), an inhibitor of respiratory syncytial virus fusion. *J. Med. Chem.* **64(7)**, 3658-3676 (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, 11/04/2024

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