

PRODUCT INFORMATION



Lenacapavir

Item No. 41830

CAS Registry No.: 2189684-44-2
Formal Name: (3bS,4aR)-N-[(1S)-1-[3-[4-chloro-3-[(methylsulfonyl)amino]-1-(2,2,2-trifluoroethyl)-1H-indazol-7-yl]-6-[3-methyl-3-(methylsulfonyl)-1-butyn-1-yl]-2-pyridinyl]-2-(3,5-difluorophenyl)ethyl]-5,5-difluoro-3b,4,4a,5-tetrahydro-3-(trifluoromethyl)-1H-cyclopropa[3,4]cyclopenta[1,2-c]pyrazole-1-acetamide
Synonyms: GS-6207, GS-714207, GS-CA2

MF: C₃₉H₃₂ClF₁₀N₇O₅S₂

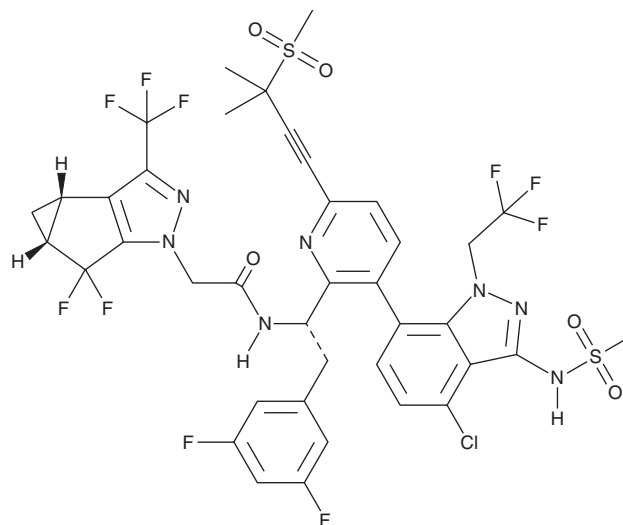
FW: 968.3

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

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

Description

Lenacapavir is an inhibitor of the protein-protein interaction between HIV capsid protein and various HIV proteins involved in viral replication.¹ It inhibits viral replication in HIV-1-infected primary human CD4⁺ T cells and macrophages (mean IC₅₀s = 32 and 56 pM, respectively), two HIV-2 clinical isolates (mean IC₅₀ = 885 pM), and 23 HIV-1 clinical isolates (mean IC₅₀ = 50 pM). Lenacapavir (25 mg/kg, s.c.) prevents simian-tropic HIV-1 (stHIV-A19) infection in pigtail macaques.² Formulations containing lenacapavir have been used in the treatment of HIV-1.

References

1. Link, J.O., Rhee, M.S., Tse, W.C., *et al.* Clinical targeting of HIV capsid protein with a long-acting small molecule. *Nature* **584**(7822), 614-618 (2020).
2. Swanstrom, A.E., Gorelick, R.J., Welker, J.L., *et al.* Long-acting lenacapavir protects macaques against intravenous challenge with simian-tropic HIV. *EBioMedicine* **95**, 104764 (2023).

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, 12/16/2024

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