

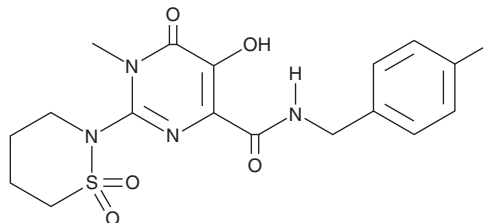
PRODUCT INFORMATION



HIV-1 Integrase Inhibitor 1

Item No. 36807

CAS Registry No.: 729607-74-3
Formal Name: N-[(4-fluorophenyl)methyl]-1,6-dihydro-5-hydroxy-1-methyl-6-oxo-2-(tetrahydro-1,1-dioxido-2H-1,2-thiazin-2-yl)-4-pyrimidinecarboxamide
Synonym: BMS-707035
MF: C₁₇H₁₉FN₄O₅S
FW: 410.4
Purity: ≥98%
UV/Vis.: λ_{max}: 253, 302 nm
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

HIV-1 integrase inhibitor 1 is supplied as a solid. A stock solution may be made by dissolving the HIV-1 integrase inhibitor 1 in the solvent of choice, which should be purged with an inert gas. HIV-1 integrase inhibitor 1 is slightly soluble in DMSO.

Description

HIV-1 integrase inhibitor 1 is an inhibitor of HIV-1 integrase DNA strand transfer (IC₅₀ = 14 nM).¹ It is also an inhibitor of severe acute respiratory syndrome coronavirus 2 (SARS-CoV-2) main protease (M^{pro}), also known as 3C-like protease (3CL^{pro}), when used at a concentration of 10 μM in a fluorescence-based assay.² HIV-1 integrase inhibitor 1 decreases SARS-CoV-2 replication (EC₅₀ = 0.27 μM) and reduces particle-forming units (PFUs) in infected Huh7.5 cells in a concentration-dependent manner.

References

- Langley, D.R., Samanta, H.K., Lin, Z., *et al.* The terminal (catalytic) adenosine of the HIV LTR controls the kinetics of binding and dissociation of HIV integrase strand transfer inhibitors. *Biochemistry* **47(51)**, 13481-13488 (2008).
- Narayanan, A., Narwal, M., Majowicz, S.A., *et al.* Identification of SARS-CoV-2 inhibitors targeting Mpro and PLpro using in-cell-protease assay. *Commun. Biol.* **5(1)**, 169 (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, 06/19/2023

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