

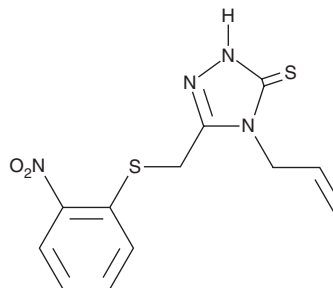
PRODUCT INFORMATION



SSYA10-001

Item No. 38324

CAS Registry No.: 675104-49-1
Formal Name: 2,4-dihydro-5-[[[(2-nitrophenyl)thio]methyl]-4-(2-propen-1-yl)-3H-1,2,4-triazole-3-thione
MF: C₁₂H₁₂N₄O₂S₂
FW: 308.4
Purity: ≥98%
UV/Vis.: λ_{max}: 252 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

SSYA10-001 is supplied as a solid. A stock solution may be made by dissolving the SSYA10-001 in the solvent of choice, which should be purged with an inert gas. SSYA10-001 is soluble in DMSO.

Description

SSYA10-001 is an inhibitor of severe acute respiratory syndrome coronavirus (SARS-CoV) helicase nsp13 (IC₅₀s = 5.3 and 5.7 μM for dsDNA and dsRNA, respectively) and the protein-protein interaction between human papillomavirus (HPV) E6 and E6-associated proteins (E6-AP; IC₅₀ = 10 μM).¹⁻³ It is selective for SARS-CoV nsp13 over hepatitis C virus (HCV) and dengue virus helicases at 2 μM.¹ SSYA10-001 inhibits the replication of SARS-CoV in HEK293T cells, Middle East respiratory syndrome CoV (MERS-CoV) in Vero E6 cells, and mouse hepatitis virus (MHV) in mouse WM-L2 fibroblasts (EC₅₀s = 8.95, 25, and 12 μM, respectively).^{1,2} It decreases HPV E6-induced p53 degradation in SiHa and HeLa cells when used at concentrations of 1, 5, and 10 μM.³

References

1. Adedeji, A.O., Singh, K., Calcaterra, N.E., *et al.* Severe acute respiratory syndrome coronavirus replication inhibitor that interferes with the nucleic acid unwinding of the viral helicase. *Antimicrob Agents Chemother.* **56**(9), 4718-4728 (2012).
2. Adedeji, A.O., Singh, K., Kassim, A., *et al.* Evaluation of SSYA10-001 as a replication inhibitor of severe acute respiratory syndrome, mouse hepatitis, and Middle East respiratory syndrome coronaviruses. *Antimicrob Agents Chemother.* **58**(8), 4894-4898 (2014).
3. Malechka, K.A., Fera, D., Schultz, D.C., *et al.* Identification and characterization of small molecule human papillomavirus E6 inhibitors. *ACS Chem. Biol.* **9**(7), 1603-1612 (2014).

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, 05/09/2023

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