

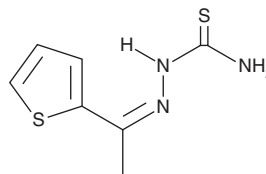
PRODUCT INFORMATION



SSAA09E1

Item No. 32582

CAS Registry No.: 433212-75-0
Formal Name: (2Z)-2-[1-(2-thienyl)ethylidene]-hydrazinecarbothioamide
MF: C₇H₉N₃S₂
FW: 199.3
Purity: ≥98%
UV/Vis.: λ_{max}: 257, 327 nm
Supplied as: A crystalline 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

SSAA09E1 is supplied as a crystalline solid. A stock solution may be made by dissolving the SSAA09E1 in the solvent of choice, which should be purged with an inert gas. SSAA09E1 is soluble in the organic solvent ethanol.

Description

SSAA09E1 is an inhibitor of severe acute respiratory syndrome coronavirus (SARS-CoV) viral entry.¹ It reduces infection of HEK293T cells transiently transfected with angiotensin-converting enzyme 2 (ACE2) by an HIV-based virus system pseudotyped with SARS-CoV spike glycoprotein (EC₅₀ = 6.7 μM). SSAA09E1 also inhibits the proteolytic activity of cathepsin L (IC₅₀ = 5.33 μM) but not cathepsin B when used at a concentration of 20 μM.

Reference

1. Adedeji, A.O., Severson, W., Jonsson, C., *et al.* Novel inhibitors of severe acute respiratory syndrome coronavirus entry that act by three distinct mechanisms. *J. Virol.* **87**(14), 8017-8028 (2013).

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, 11/22/2022

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