

PRODUCT INFORMATION



SSAA09E3

Item No. 31386

CAS Registry No.: 52869-18-8

Formal Name: N-(9,10-dihydro-9,10-dioxo-2-anthracenyl)-benzamide

MF: $C_{21}H_{13}NO_3$

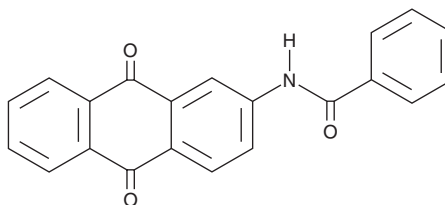
FW: 327.3

Purity: $\geq 98\%$

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

SSAA09E3 is supplied as a solid. A stock solution may be made by dissolving the SSAA09E3 in the solvent of choice, which should be purged with an inert gas. SSAA09E3 is soluble (≥ 10 mg/ml) in DMSO.

Description

SSAA09E3 is an inhibitor of severe acute respiratory syndrome coronavirus (SARS-CoV) viral membrane fusion with the host cell membrane.¹ It inhibits SARS/HIV pseudotyped viral entry into HEK293T cells ($EC_{50} = 9.7 \mu\text{M}$) and inhibits SARS-CoV infection of Vero cells ($EC_{50} = 0.15 \mu\text{M}$).

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, 11/25/2025

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