

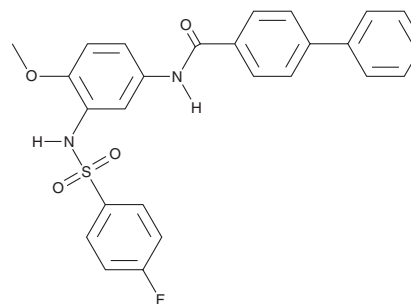
PRODUCT INFORMATION



SN-001

Item No. 38899

CAS Registry No.: 727699-84-5
Formal Name: N-[3-[[[4-(4-fluorophenyl)sulfonyl]amino]-4-methoxyphenyl]-[1,1'-biphenyl]-4-carboxamide
Synonyms: DUN99845, Stimulator of Interferon Genes Inhibitor 1, STING Inhibitor 1
MF: C₂₆H₂₁FN₂O₄S
FW: 476.5
Purity: ≥95%
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

SN-001 is supplied as a solid. A stock solution may be made by dissolving the SN-001 in the solvent of choice, which should be purged with an inert gas. SN-001 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of SN-001 in ethanol is approximately 5 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Description

SN-001 is an inhibitor of stimulator of interferon genes (STING).¹ It inhibits herring testis DNA-induced *lfnb* mRNA expression in L-929 cells (IC₅₀ = 3.82 μM).

Reference

1. Hong, Z., Mei, J., Li, C., *et al.* STING inhibitors target the cyclic dinucleotide binding pocket. *Proc. Natl. Acad. Sci. USA* **118**(24), e2105465118 (2021).

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, 07/06/2023

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