

PRODUCT INFORMATION

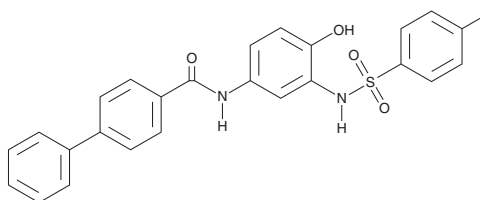


SN-011

Item No. 34890

CAS Registry No.: 2249435-90-1
Formal Name: N-[3-[[[4-(4-fluorophenyl)sulfonyl]amino]-4-hydroxyphenyl]-[1,1'-biphenyl]-4-carboxamide

Synonym: GUN35901
MF: C₂₅H₁₉FN₂O₄S
FW: 462.5
Purity: ≥98%
UV/Vis.: λ_{max}: 277 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

SN-011 is supplied as a solid. A stock solution may be made by dissolving the SN-011 in the solvent of choice, which should be purged with an inert gas. SN-011 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of SN-011 in these solvents is 3, 25, and 5 mg/ml, respectively.

Description

SN-011 is a stimulator of interferon genes (STING) antagonist.¹ It binds to the STING cyclic dinucleotide binding site ($K_d = 4.03$ nM) and inhibits 2'3'-cGAMP-induced *Irfn* expression in mouse embryonic fibroblasts (MEFs), mouse bone marrow-derived macrophages (BMDMs), and human foreskin fibroblasts (HFFs; IC_{50} s = 127.5, 107.1, and 502.8 nM, respectively). SN-011 (1 μM) impairs recruitment of IFN regulatory factor 3 (IRF3) or TANK-binding kinase 1 (TBK1) to the STING signalosome in HEK293T cells overexpressing tagged wild-type or SAVI-linked mutant STING and IRF3 or TBK1, as well as inhibits translocation of STING from the endoplasmic reticulum (ER) to the Golgi induced by herpes simplex virus 1 (HSV-1) in HFFs. *In vivo*, SN-011 (5 mg/kg) increases survival and reduces *Irfn* mRNA levels in the *Trex1*^{-/-} mouse model of Aicardi-Goutières syndrome, an autoimmune disorder characterized by constitutive activation of cGAS and IFN overproduction.

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, 11/23/2022

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