

PRODUCT INFORMATION



NVS-STG2

Item No. 40159

CAS Registry No.: 3030588-01-0

Formal Name: 4-[[[4-[2-(1,1-dimethylethyl)-5,5-dimethyl-1,3-dioxan-2-yl]phenyl]methyl]amino]-3-methoxy-benzoic acid

MF: $C_{25}H_{33}NO_5$

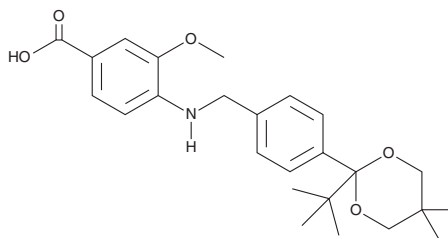
FW: 427.5

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

NVS-STG2 is supplied as a solid. A stock solution may be made by dissolving the NVS-STG2 in the solvent of choice, which should be purged with an inert gas. NVS-STG2 is slightly soluble (0.1-1 mg/ml) in DMSO.

Description

NVS-STG2 is an agonist of stimulator of interferon genes (STING).¹ It binds to STING and induces STING activation in a reporter assay ($IC_{50} = 5.2 \mu\text{M}$). Intratumoral injection of NVS-STG2 (400 or 800 $\mu\text{g}/\text{animal}$) reduces tumor volume and increases intratumoral IFN- γ levels in a B16.SIY murine melanoma model.

Reference

1. Li, J., Canham, S.M., Wu, H., *et al.* Activation of human STING by a molecular glue-like compound. *Nat. Chem. Biol.* **20**(3), 365-372 (2024).

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, 10/22/2025

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