NVS-STG2

Molecular glue STING agonist

NVS-STG2 is a STING agonist that activates human STING by binding in a pocket between the transmembrane domains of neighboring STING dimers in a molecular glue-like manner. Inactive with mouse STING. NVS-STG2 slowed tumor growth significantly in MC38 tumors inoculated into hSTING knock-in mice.

Ordering Information

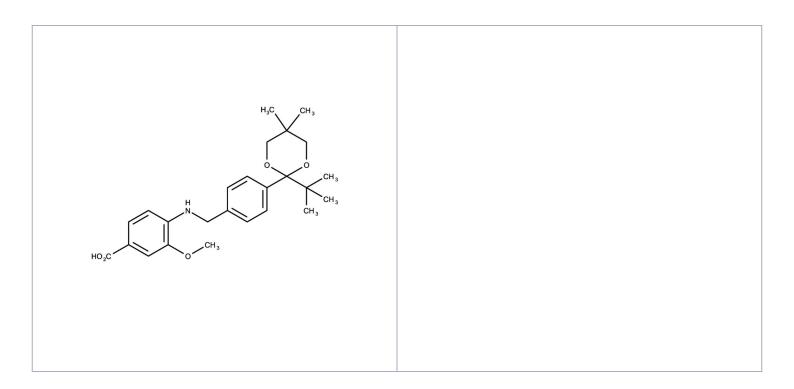
Order Online »

ENZ-CHM450-0025

25mg

Manuals, SDS & CofA

View Online »



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 4-((4-(2-(tert-Butyl)-5,5-dimethyl-1,3-dioxan-2-

yl)benzyl)amino)-3-methoxybenzoic acid

Appearance White solid.

CAS 3030588-01-0

Couple Type Activator

Formula ${\rm C}_{25}{\rm H}_{33}{\rm NO}_5$

Identity Determined by NMR.

MW 427.54

Purity ≥98% (HPLC)

Solubility Soluble in DMSO (40 mg/ml).

Last modified: January 24, 2025