

# 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.

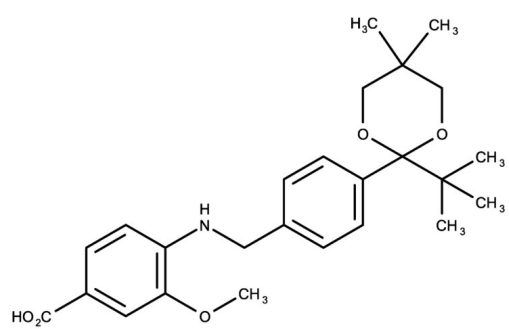
### Ordering Information

[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**  $C_{25}H_{33}NO_5$

**Identity** Determined by NMR.

**MW** 427.54

**Purity** ≥98% (HPLC)

**Solubility** Soluble in DMSO (40 mg/ml).



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)