

H-151

STING inhibitor

Catalog code: inh-h151

<https://www.invivogen.com/h151>

For research use only

Version 19G19-MM

PRODUCT INFORMATION

Contents

- 10 mg H-151

Storage and stability

- H-151 is shipped at room temperature. Upon receipt, store at -20°C.
- Upon resuspension, prepare aliquots and store at -20°C. Resuspended product is stable for at least 3 months when properly stored. Avoid repeated freeze-thaw cycles.

Quality control

- Purity ≥ 95% (UHPLC)
- The absence of bacterial contamination (e.g. lipoproteins and endotoxins) has been confirmed using HEK-Blue™ TLR2 and HEK-Blue™ TLR4 cells.
- The inhibitory activity has been validated using cellular assays.

BACKGROUND

STING (stimulator of interferon genes) has become a focal point in immunology research and drug discovery^{1,2}. In a healthy individual, STING functions as a signaling hub, orchestrating immune responses to pathogenic, tumoral, or self-DNA detected in the cytoplasm². Upon activation, STING induces type I interferon (IFN) production through TANK-binding-kinase-1 (TBK1)-mediated IFN regulatory factor (IRF3) signaling². STING activation also leads to NF-κB-dependent inflammatory cytokine production². In some autoimmune diseases such as STING-associated vasculopathy with onset in infancy (SAVI), STING is constitutively activated resulting in high IFN production^{3,4}. The discovery of a mechanism to pharmacologically inhibit STING should lead to new treatments for such diseases.

PRODUCT DESCRIPTION

H-151 is a potent, irreversible and selective small molecule inhibitor of STING¹. This synthetic indole-derivative exerts its inhibitory action by covalently binding to STING at the transmembrane cysteine residue at position 91. H-151 blocks STING palmitoylation and clustering, two essential steps for STING signaling. Of note, H-151 potently inhibits both human and murine STING, *in vitro* and *in vivo* mouse models. Indeed, in models of autoinflammatory disease, H-151 blocks STING-induced expression of pro-inflammatory cytokines and reduces inflammation¹. Notably, H-151 is effective against all the STING variants tested, including constitutively active disease-associated mutants such as S154 (N154S) and M155 (V155M; see validation data sheet available on our website).

1. Haag S.M. *et al.*, 2018. Targeting STING with covalent small-molecule inhibitors. *Nature* 559:269-73. 2. Ishikawa H. & Barber G.N. 2008. STING is an endoplasmic reticulum adaptor that facilitates innate immune signalling. *Nature* 455:674-8. 3. Liu Y. *et al.*, 2014. Activated STING in a vascular and pulmonary syndrome. *N Engl J Med.* 371:507-18. 4. Jeremiah N. *et al.*, 2013. Inherited STING-activating mutation underlies a familial inflammatory syndrome with lupus-like manifestations. *J Clin Invest.* 124:5516-20.

TECHNICAL SUPPORT

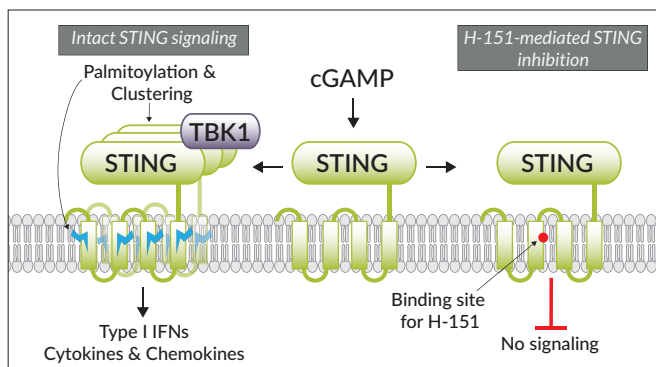
InvivoGen USA (Toll-Free): 888-457-5873

InvivoGen USA (International): +1 (858) 457-5873

InvivoGen Europe: +33 (0) 5-62-71-69-39

InvivoGen Hong Kong: +852 3622-3480

E-mail: info@invivogen.com



Mechanism of action of H-151-mediated STING inhibition.

CHEMICAL PROPERTIES

CAS number: 941987-60-6

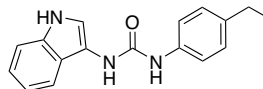
Synonym: N-(4-Ethylphenyl)-N'-1H-indol-3-yl-urea

Solubility: 20 mg/ml (71.60 mM) in DMSO

Formula: C₁₇H₁₇N₃O

Molecular weight: 279.34 g/mol

Structure:



METHODS

Preparation of 10 mg/ml (35.8 mM) stock solution

1. Add 1 ml of DMSO to 10 mg of H-151. Mix by vortexing.
2. Use immediately or store aliquots at -20°C.
3. Prepare a 1:10 dilution with DMSO to obtain a 1 mg/ml solution.
4. Further dilutions of the 1 mg/ml solution can be prepared using culture medium, such as RPMI or DMEM, containing 10% fetal calf serum.

Note: Dilutions in water or PBS may cause the product to precipitate.

Working concentration range: 4 ng/ml (15 nM) to 4 µg/ml (15 µM) for cellular assays

STING inhibition assay

On the next page is a protocol to study STING inhibition in THP1-Dual™ cells. These cells derive from the human monocytic cell line THP-1, by stable integration of two inducible reporters allowing the simultaneous study of the IRF pathway, by assessing the activity of the secreted luciferase Lucia, and the NF-κB pathway, by monitoring the activity of SEAP. These cells have been shown to express STING and respond to STING agonists.

For more information, visit <https://www.invivogen.com/thp1-dual>.

Protocol for STING inhibition in THP1-Dual™ cells

1. Add 20 µl of H-151 (final concentration 4 ng/ml to 4 µg/ml) per well of a flat-bottom 96-well plate.
2. Add 160 µl of cell suspension (~100,000 cells) per well.
3. Incubate for 2 hours at 37 °C in a 5% CO₂ incubator.
4. Add 20 µl of a test sample or a STING agonist, such as 2'3'-cGAMP (final concentration 20 µg/ml) per well.
5. Incubate the plate for 18-24 hours at 37 °C in a 5% CO₂ incubator.
6. Monitor IRF and NF-κB activation by measuring the levels of Lucia luciferase and SEAP using QUANTI-Luc™ and QUANTI-Blue™ Solution, respectively.

STING inhibition assays can also be performed in other cells lines from InvivoGen's STING reporter cell collection. InvivoGen has developed stable reporter cells where the wild-type STING gene has been replaced by a STING variant using knock-in technology (KI-STING). These cells feature IRF-inducible Lucia luciferase and an NF-κB SEAP (secreted embryonic alkaline phosphatase) secreted reporter proteins as convenient read-outs.

For more information, visit <https://www.invivogen.com/sting-reporter-cells>.

RELATED PRODUCTS

Product	Cat. Code
STING Agonists	
2'3'-cGAMP	tlrl-nacga23
3'3'-cGAMP Fluorinated	tlrl-nacgaf
2'3'-c-di-AM(PS)2 (Rp,Rp)	tlrl-nacda2r
Detection Reagents	
QUANTI-Blue™ Solution	rep-qbs
QUANTI-Luc™	rep-qlc1
Mouse Macrophage Reporter Cells Lines	
RAW-Lucia™ ISG Cells	rawl-isg
RAW-Lucia™ ISG-KO-STING cells	rawl-kostg
Human Monocytic Reporter Cells Lines	
THP1-Dual™ Cells	thpd-nfis
THP1-Dual™ KO-STING Cells	thpd-kostg
THP1-Dual™ KI-hSTING-A162 Cells	thpd-a162
THP1-Dual™ KI-hSTING-H232 Cells	thpd-h232
THP1-Dual™ KI-hSTING-M155 Cells	thpd-m155
THP1-Dual™ KI-hSTING-R232 Cells	thpd-r232
THP1-Dual™ KI-hSTING-S154 Cells	thpd-s154

TECHNICAL SUPPORT

InvivoGen USA (Toll-Free): 888-457-5873

InvivoGen USA (International): +1 (858) 457-5873

InvivoGen Europe: +33 (0) 5-62-71-69-39

InvivoGen Hong Kong: +852 3622-3480

E-mail: info@invivogen.com