

diABZI

diABZI (compound 3) trihydrochloride; non-CDN STING agonist

Catalog code: tlr-diabzi

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

For research use only

Version 22K25-MM

PRODUCT INFORMATION

Contents

- 1 mg diABZI
- 1.5 ml endotoxin-free water

Storage and stability

- diABZI is provided as a lyophilized product and shipped at room temperature. Upon receipt, store product at -20°C.
- Resuspended product is stable for up to 3 months when properly stored at -20°C.
- Avoid repeated freeze-thaw cycles.

Quality control

- Purity: ≥95% (UHPLC)
- Activation of STING has been confirmed using cellular assays.
- The absence of bacterial contamination (e.g. lipoproteins and endotoxins) has been confirmed using HEK-Blue™ cellular assays.

PRODUCT DESCRIPTION

diABZI, also known as diABZI (compound 3) trihydrochloride, is a small molecule, non-cyclic dinucleotide, that potently activates STING. Notably, diABZI (compound 3) trihydrochloride is water soluble.

Development of a non-CDN STING agonist

A series of small molecule amidobenzimidazoles (ABZI) were identified to effectively compete with the classical STING agonist, 2'3'-cGAMP. To take advantage of the symmetrical nature of STING, two molecules of the lead compound were joined to create a single dimeric ligand (diABZI; compound 2). Importantly, this afforded a 1000x increase in the STING binding affinity, when compared to 2'3'-cGAMP. This dimeric ligand was then further optimized into 'diABZI (compound 3)'¹.

Activation of STING by diABZI

In contrast to classical CDNs, diABZI activates STING while maintaining its open conformation¹. Further research is needed to understand the implications of this difference. However, similar to 2'3'-cGAMP, diABZI induces STING-dependent activation of type-I interferon and pro-inflammatory cytokines *in vitro* and *in vivo*¹.

diABZI has therapeutic potential in the treatment of cancer with significant inhibition of tumor growth observed in a syngeneic mouse model of colorectal cancer¹. Furthermore, diABZI has been shown to suppress infection by diverse strains of SARS-CoV-2, including variants of concern (e.g. B.1.351), through the induction of a effective IFN response².

1. Ramanjulu, J.M. et al. 2018. Design of amidobenzimidazole STING receptor agonists with systemic activity. *Nature* 564, 439-443. **2. Li, M. et al. 2021.** Pharmacological activation of STING blocks SARS-CoV-2 infection. *Science* 6(59), eabi9007.

CHEMICAL PROPERTIES

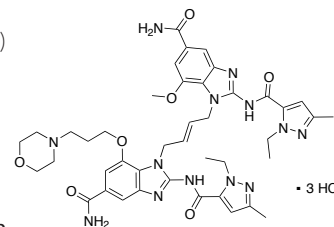
CAS Number: 2138299-34-8

Synonyms: diABZI (Compound 3)

Formula: C₄₂H₅₁N₁₃O₇•3HCl

Molecular weight: 959.33 g/mol

Solubility: 2mg/ml H₂O



METHODS

Preparation of a stock solution (1 mg/ml; 1.04 mM)

1. Before opening the vial, centrifuge briefly and open the lid carefully to avoid any loss of product.
2. Add 1 ml of endotoxin-free water (provided) to the vial of diABZI to obtain a stock solution at 1 mg/ml.
3. Vortex until completely dissolved.

Working concentration: 0.01 - 30 µM

Activation of STING in THP1-Dual™ cells

Below is a protocol for monitoring the activation of STING by diABZI using InvivoGen's THP1-Dual™ cells. These cells allow the simultaneous study of the NF-κB pathway, by monitoring the activity of SEAP, and the IRF (interferon regulatory factor) pathway, by assessing the activity of the secreted Lucia luciferase. For more information, please visit <https://www.invivogen.com/thp1-dual>.

1. Add 20 µl of diABZI (10x final concentration) per well of a flat-bottom 96-well plate.
2. Add 20 µl of a positive control (i.e. 2'3'-cGAMP) to another well.
3. Prepare a suspension of THP1-Dual™ cells (~500,000 cells per ml) as detailed in the cell line data sheet.
4. Add 180 µl of cell suspension (~100,000 cells) per well.
5. Incubate the plate at 37°C in a 5% CO₂ incubator for 18-24 hours.
6. Prepare QUANTI-Luc™ (IRF assessment) and/or QUANTI-Blue™ Solution (NF-κB assessment) and carry out the measurements following the instructions on the data sheet.

RELATED PRODUCTS

Product	Description	Cat. Code
2'3'-cGAMP	STING ligand	tlrl-nacga23
cAIM(PS)2 Difluor (Rp/Sp)	STING ligand	tlrl-nacairs
THP1-Dual™ cells	Reporter monocytes	thpd-nfis
QUANTI-Blue™ Solution	SEAP detection reagent	rep-qbs
QUANTI-Luc™	Luciferase detection reagent	rep-qlc1

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 Asia: +852 3622-3480

E-mail: info@invivogen.com