

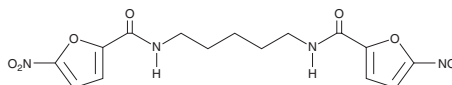
# PRODUCT INFORMATION



## STING Agonist 23

Item No. 38345

**Formal Name:** N,N'-(pentane-1,5-diyl)bis(5-nitrofur-2-carboxamide)  
**Synonym:** Stimulator of Interferon Genes Agonist 23  
**MF:** C<sub>15</sub>H<sub>16</sub>N<sub>4</sub>O<sub>8</sub>  
**FW:** 380.3  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 215, 308 nm  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

STING agonist 23 is supplied as a solid. A stock solution may be made by dissolving the STING agonist 23 in the solvent of choice, which should be purged with an inert gas. STING agonist 23 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of STING agonist 23 in these solvents is approximately 14 and 16 mg/ml, respectively.

### Description

STING agonist 23 is an agonist of stimulator of interferon genes (STING).<sup>1</sup> It increases the phosphorylation of TANK-binding kinase 1 (TBK1) in wild-type, but not STING-deficient, HT-1080 cells when used at a concentration of 20 μM. STING agonist 23 induces dimerization of STING in THP-1 cells and mouse bone marrow-derived macrophages (BMDMs) in a concentration-dependent manner. It inhibits the replication of herpes simplex virus (HSV), rhesus rotavirus (RRV), and severe acute respiratory syndrome coronavirus 2 (SARS-CoV-2) *in vitro* (EC<sub>50</sub>s = 6.15, 6.8, and 10.71 μM, respectively). *In vivo*, STING agonist 23 increases serum levels of IFN-β and TNF-α in mice.

### Reference

1. Zang, R., Xue, L., Zhang, M., *et al.* Design and syntheses of a bimolecular STING agonist based on the covalent STING antagonist. *Eur. J. Med. Chem.* **250**, 115184 (2023).

#### 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, 05/02/2023

#### 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