

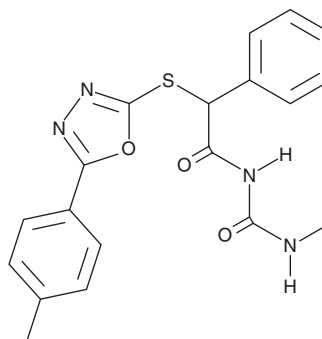
# PRODUCT INFORMATION



## STING Agonist C11

Item No. 27886

**CAS Registry No.:** 875863-22-2  
**Formal Name:** N-[(methylamino)carbonyl]- $\alpha$ -[[5-(4-methylphenyl)-1,3,4-oxadiazol-2-yl]thio]-benzeneacetamide  
**Synonym:** Stimulator of Interferon Genes Agonist C11  
**MF:** C<sub>19</sub>H<sub>18</sub>N<sub>4</sub>O<sub>3</sub>S  
**FW:** 382.4  
**Purity:**  $\geq$ 95%  
**UV/Vis.:**  $\lambda_{\text{max}}$ : 274 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:**  $\geq$ 4 years



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

### Laboratory Procedures

STING agonist C11 is supplied as a crystalline solid. A stock solution may be made by dissolving the STING agonist C11 in the solvent of choice, which should be purged with an inert gas. STING agonist C11 is soluble in the organic solvent DMSO.

### Description

STING agonist C11 is an agonist of the stimulator of interferon genes (STING) pathway.<sup>1</sup> It induces secretion of type I IFN from THF and MM6 cells when used at a concentration of 50  $\mu$ M. STING agonist C11 induces phosphorylation of IFN regulatory factor 3 (IRF3) and increases expression of IFIT1 and viperin, but not IL-1 $\beta$ , IL-6, or IL-8 in THF cells in a STING-dependent manner. It reduces viral titers of chikungunya, Venezuelan equine encephalitis, o'nyong-nyong, Mayaro, and Ross River viruses grown in THF cells ( $EC_{50}$ s = 16.44, 16.7, 18.84, 25.19, and 22.57  $\mu$ M, respectively), an effect that is dependent on the presence of STING and the IFN- $\alpha$ / $\beta$  receptor (IFNAR).

### Reference

1. Gall, B., Pryke, K., Abraham, J., *et al.* Emerging alphaviruses are sensitive to cellular states induced by a novel small-molecule agonist of the STING pathway. *J. Virol.* **92(6)**, e01913-17 (2018).

#### 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, 07/19/2023

#### 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