

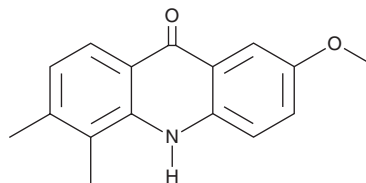
PRODUCT INFORMATION



STING Agonist 12b

Item No. 32521

CAS Registry No.: 2411100-70-2
Formal Name: 2-methoxy-5,6-dimethyl-9(10H)-acridinone
Synonym: Stimulator of Interferon Genes Agonist 12b
MF: C₁₆H₁₅NO₂
FW: 253.3
Purity: ≥98%
UV/Vis.: λ_{max}: 271 nm
Supplied as: A crystalline 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 12b is supplied as a crystalline solid. A stock solution may be made by dissolving the STING agonist 12b in the solvent of choice, which should be purged with an inert gas. STING agonist 12b is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of STING agonist 12b in these solvents is approximately 1 mg/ml.

Description

STING agonist 12b is an agonist of stimulator of interferon genes (STING).¹ It binds to STING (K_d = 26.4 μM) and induces interferon reporter gene expression in cells expressing human or mouse STING (EC₅₀s = 7.45 and 10.23 μM, respectively). STING agonist 12b (40 μM) induces expression of TNF-α, IL-6, IP-10, and IL-1β in THP-1 cells.

Reference

1. Hou, S., Lan, X.-j., Li, W., *et al.* Design, synthesis and biological evaluation of acridone analogues as novel STING receptor agonists. *Bioorg. Chem.* **95**, 103556 (2020).

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