

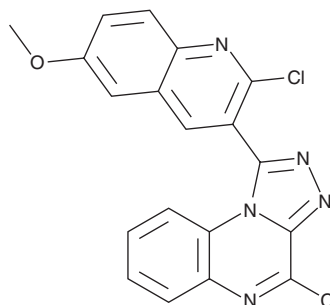
PRODUCT INFORMATION



STING Agonist 1a

Item No. 32520

CAS Registry No.: 652142-94-4
Formal Name: 4-chloro-1-(2-chloro-6-methoxy-3-quinolinyl)-[1,2,4]triazolo[4,3-a]quinoxaline
Synonym: Stimulator of Interferon Genes Agonist 1a
MF: C₁₉H₁₁Cl₂N₅O
FW: 396.2
Purity: ≥98%
UV/Vis.: λ_{max}: 218, 250 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

STING agonist 1a is supplied as a crystalline solid. A stock solution may be made by dissolving the STING agonist 1a in the solvent of choice, which should be purged with an inert gas. STING agonist 1a is slightly soluble in DMSO and dimethyl formamide.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of STING agonist 1a can be prepared by directly dissolving the crystalline solid in aqueous buffers. STING agonist 1a is slightly soluble in PBS, pH 7.2. We do not recommend storing the aqueous solution for more than one day.

Description

STING agonist 1a is an agonist of stimulator of interferon genes (STING).¹ It induces expression of an IRF-inducible SEAP reporter gene in a cell-based assay (EC₅₀ = 16.77 μM). STING agonist 1a (12.5-100 μM) induces expression of IFN-β, IL-6, and chemokine (C-X-C motif) ligand 10 (CXCL10) in THP-1 cells, an effect that can be reversed by STING knockout or the STING inhibitor H-151 (Item No. 25857).

Reference

1. Hou, H., Yang, R., Liu, X., *et al.* Discovery of triazoloquinoxaline as novel STING agonists via structure-based virtual screening. *Bioorg. Chem.* **100**, 103958 (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, 03/21/2024

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