

PRODUCT INFORMATION



STING18

Item No. 28319

CAS Registry No.: 2706536-26-5
Formal Name: 2-[3-chloro-4-(1,1-dimethylethyl)phenyl]-3-(2,3-dihydro-1,4-benzodioxin-6-yl)-7-fluoro-1,2,3,4-tetrahydro-1-oxo-4-isoquinolineacetic acid

Synonym: Stimulator of Interferon Genes 18

MF: $C_{29}H_{27}ClFNO_5$

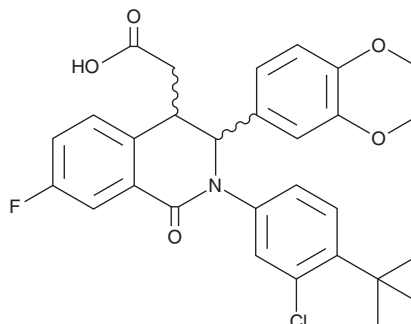
FW: 524.0

Purity: $\geq 98\%$

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

STING18 is supplied as a crystalline solid. A stock solution may be made by dissolving the STING18 in the solvent of choice, which should be purged with an inert gas. STING18 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of STING18 in these solvents is approximately 5 and 10 mg/ml, respectively.

STING18 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, STING18 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. STING18 has a solubility of approximately 0.11 mg/ml in a 1:8 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

STING18 is a competitive ligand of stimulator of interferon genes (STING; $IC_{50} = 0.068 \mu\text{M}$ in a radioligand binding assay).¹ It inhibits cGAMP-induced IFN- β production ($IC_{50} = 11 \mu\text{M}$) but does not stimulate IFN- β production ($EC_{50} = >30 \mu\text{M}$) in THP-1 cells.

Reference

1. Siu, T., Altman, M.D., Baltus, G.A., *et al.* Discovery of a novel cGAMP competitive ligand of the inactive form of STING. *Med. Chem. Lett.* **10(1)**, 92-97 (2019).

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, 03/18/2024

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