

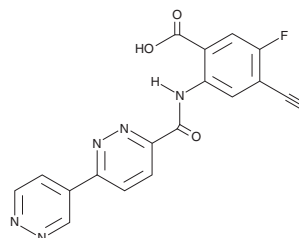
PRODUCT INFORMATION



STING Agonist 12L

Item No. 38344

CAS Registry No.: 2913152-30-2
Formal Name: 2-([3,4'-bipyridazine]-6-carboxamido)-4-ethynyl-5-fluorobenzoic acid
Synonym: Stimulator of Interferon Genes Agonist 12L
MF: C₁₈H₁₀FN₅O₃
FW: 363.3
Purity: ≥98%
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 12L is supplied as a solid. A stock solution may be made by dissolving the STING agonist 12L in the solvent of choice, which should be purged with an inert gas. STING agonist 12L is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of STING agonist 12L in ethanol is approximately 5 mg/ml and approximately 10 mg/ml in DMSO and DMF.

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 12L can be prepared by directly dissolving the solid in aqueous buffers. The solubility of STING agonist 12L in PBS (pH 7.2) is approximately 2 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

STING agonist 12L is an agonist of stimulator of interferon genes (STING).¹ It binds to wild-type STING (IC₅₀ = 1.15 μM), as well as the STING R232, AQ, and Q variants (IC₅₀s = 1.06, 0.61, and 1.12 μM, respectively) and induces reporter gene expression in THP-1 and RAW 264.7 cells (EC₅₀s = 0.38 and 12.94 μM, respectively). STING agonist 12L (5 μM) induces *IFNB1*, *CXCL10*, and *IL6* mRNA expression in THP-1 cells. *In vivo*, STING agonist 12L (10 mg/kg) increases plasma IFN-β levels and reduces tumor volume and the number of lung metastases in a B16/F10 murine melanoma model.

Reference

1. Shan, B., Hou, H., Zhang, K., *et al.* Design, synthesis, and biological evaluation of bipyridazine derivatives as stimulator of interferon genes (STING) receptor agonists. *J. Med. Chem.* **66**(5), 3327-3347 (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, 07/21/2023

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