

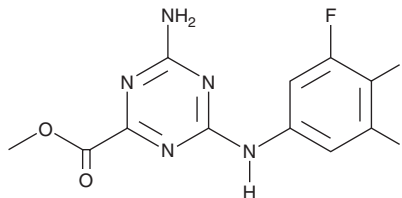
PRODUCT INFORMATION



CU-76

Item No. 31030

CAS Registry No.: 2400954-58-5
Formal Name: methyl 4-amino-6-((3,5-difluoro-4-iodophenyl)amino)-1,3,5-triazine-2-carboxylate
MF: C₁₁H₈F₂IN₅O₂
FW: 407.1
Purity: ≥95%
UV/Vis.: λ_{max}: 264 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

CU-76 is supplied as a crystalline solid. A stock solution may be made by dissolving the CU-76 in the solvent of choice, which should be purged with an inert gas. CU-76 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of CU-76 in these solvents is approximately 30 mg/ml.

CU-76 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CU-76 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. CU-76 has a solubility of approximately 0.1 mg/ml in a 1:5 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

CU-76 is an inhibitor of cyclic GMP-AMP synthase (cGAS; IC₅₀ = 0.24 μM).¹ It reduces DNA-, but not Sendai virus-, induced dimerization of IFN regulatory factor 3 (IRF3; Item No. 22811) in THP-1 cells, indicating selectivity for the cGAS DNA sensing pathway over the RIG-I-MAVS RNA sensing pathway. CU-76 decreases IFN-stimulatory DNA-induced production of IFN-β in THP-1 cells when used at concentrations of 10, 30, and 100 μM.

Reference

1. Padilla-Salinas, R., Sun, L., Anderson, R., *et al.* Discovery of small-molecule cyclic GMP-AMP synthase inhibitors. *J. Org. Chem.* **85**(3), 1579-1600 (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, 06/25/2020

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