

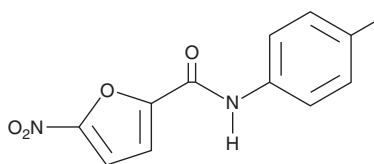
PRODUCT INFORMATION



C-176

Item No. 25859

CAS Registry No.: 314054-00-7
Formal Name: N-(4-iodophenyl)-5-nitro-2-furancarboxamide
MF: C₁₁H₇IN₂O₄
FW: 358.1
Purity: ≥98%
UV/Vis.: λ_{max}: 258, 320 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

C-176 is supplied as a crystalline solid. A stock solution may be made by dissolving the C-176 in the solvent of choice, which should be purged with an inert gas. C-176 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of C-176 in these solvents is approximately 30 mg/ml.

C-176 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, C-176 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. C-176 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

C-176 is a covalent inhibitor of stimulator of interferon genes (STING).¹ It selectively reduces STING-, but not RIG-I or TBK1-, mediated IFN-β reporter activity in HEK293 cells when used at concentrations ranging from 0.01 to 5 μM. C-176 reduces serum levels of type-I interferons (IFNs) and cardiac inflammation in the *Trex^{-/-}* mouse model of multi-organ inflammation.

Reference

1. Haag, S.M., Gulen, M.F., Reymond, L., *et al.* Targeting STING with covalent small-molecule inhibitors. *Nature* **559(7713)**, 269-273 (2018).

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, 11/30/2022

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