

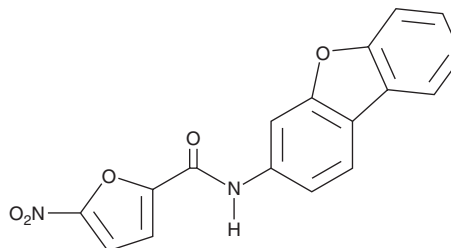
PRODUCT INFORMATION



C-178

Item No. 25860

CAS Registry No.: 329198-87-0
Formal Name: N-3-dibenzofuranyl-5-nitro-2-furancarboxamide
MF: C₁₇H₁₀N₂O₅
FW: 322.3
Purity: ≥98%
UV/Vis.: λ_{max}: 212, 262, 300 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-178 is supplied as a crystalline solid. A stock solution may be made by dissolving the C-178 in the solvent of choice, which should be purged with an inert gas. C-178 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of C-178 in these solvents is approximately 0.25, 20, and 14.2 mg/ml, respectively.

Description

C-178 is a covalent inhibitor of stimulator of interferon genes (STING).¹ It binds to Cys91 on STING to block its palmitoylation and prevents recruitment and phosphorylation of TBK1 in HEK293T cells. C-178 (0.01-1.25 μM) selectively reduces STING-, but not RIG-I- or TBK1-, mediated IFN-β reporter activity in HEK293 cells. It also prevents increases in *Irfn1* expression in bone marrow-derived macrophages (BMDMs) induced by cyclic di-GMP (Item No. 17144), double-stranded DNA, and LPS when used at a concentration of 0.5 μM.

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, 12/21/2022

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