

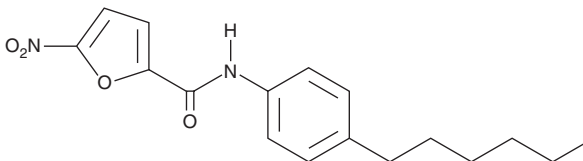
PRODUCT INFORMATION



C-171

Item No. 30159

CAS Registry No.: 2244881-69-2
Formal Name: N-(4-hexylphenyl)-5-nitro-2-furancarboxamide
MF: C₁₇H₂₀N₂O₄
FW: 316.4
Purity: ≥98%
UV/Vis.: λ_{max}: 252, 309 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

C-171 is supplied as a crystalline solid. A stock solution may be made by dissolving the C-171 in the solvent of choice, which should be purged with an inert gas. C-171 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of C-171 in these solvents is approximately 25 mg/ml.

Description

C-171 is an inhibitor of stimulator of interferon genes (STING).^{1,2} It binds to STING, inhibits its palmitoylation, and prevents the recruitment and phosphorylation of TBK1.¹ It selectively reduces human and mouse STING-, but not RIG-I or TBK1-, mediated IFN-β reporter activity in HEK293 cells when used at concentrations ranging from 0.02 to 2 μM.

References

1. Haag, S.M., Gulen, M.F., Reymond, L., *et al.* Targeting STING with covalent small-molecule inhibitors. *Nature* **559**(7713), 269-273 (2018).
2. Zhang, H., You, Q.-D., and Xu, X.-L. Targeting stimulator of interferon genes (STING): A medicinal chemistry perspective. *J. Med. Chem.* **63**(8), 3785-3816 (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, 05/07/2020

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