

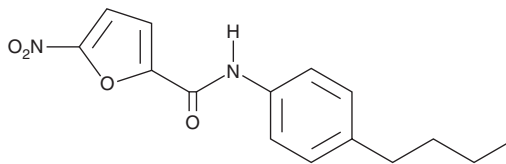
# PRODUCT INFORMATION



## C-170

Item No. 30157

CAS Registry No.: 346691-38-1  
Formal Name: N-(4-butylphenyl)-5-nitro-2-furancarboxamide  
MF:  $C_{15}H_{16}N_2O_4$   
FW: 288.3  
Purity:  $\geq 98\%$   
UV/Vis.:  $\lambda_{\max}$ : 254, 304 nm  
Supplied as: A crystalline solid  
Storage:  $-20^{\circ}\text{C}$   
Stability:  $\geq 4$  years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

C-170 is supplied as a crystalline solid. A stock solution may be made by dissolving the C-170 in the solvent of choice, which should be purged with an inert gas. C-170 is soluble in the organic solvent DMSO.

### Description

C-170 is an inhibitor of stimulator of interferon genes (STING).<sup>1,2</sup> It binds to STING, inhibits its palmitoylation, and prevents the recruitment and phosphorylation of TBK1.<sup>1</sup> It selectively reduces human and mouse STING-, but not RIG-I- or TBK1-, mediated IFN- $\beta$  reporter activity in HEK293 cells when used at concentrations ranging from 0.02 to 2  $\mu\text{M}$ .

### 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, 03/21/2024

#### 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