

# PRODUCT INFORMATION



NC03

Item No. 45389

CAS Registry No.: 568558-66-7

Formal Name: N-(2,3-dihydro-1,4-benzodioxin-6-yl)-2-[[5-(3,4,5-trimethoxyphenyl)-1,3,4-oxadiazol-2-yl]thio]-acetamide

MF: C<sub>21</sub>H<sub>21</sub>N<sub>3</sub>O<sub>7</sub>S

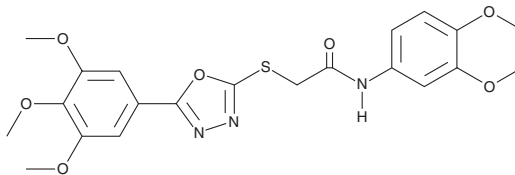
FW: 459.5

Purity: ≥98%

Supplied as: A 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

NC03 is supplied as a solid. A stock solution may be made by dissolving the NC03 in the solvent of choice, which should be purged with an inert gas. NC03 is sparingly soluble (1-10 mg/ml) in DMSO.

## Description

NC03 is an inhibitor of PI4KII $\alpha$ .<sup>1</sup> It reduces the levels of the PI4KII $\alpha$  product phosphatidylinositol 4-phosphate (PI4P) in the Golgi apparatus and endosomes of COS-7 cells when used at a concentration of 10  $\mu$ M. NC03 also reduces PI4P production by 40-50% in the Rab7<sup>+</sup> late endosomes of HEK-AT1 cells that express the rat angiotensin II type 1a (AT<sub>1a</sub>) receptor.

## Reference

1. Sengupta, N., Jović, M., Barnaeva, E., *et al.* A large scale high-throughput screen identifies chemical inhibitors of phosphatidylinositol 4-kinase type II alpha. *J. Lipid Res.* **60**(3), 683-693 (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, 04/02/2026

## 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