

# PRODUCT INFORMATION



**NDI-091143**

Item No. 36054

**CAS Registry No.:** 2375840-87-0  
**Formal Name:** 3-chloro-5-[[[(4,6-difluoro[1,1'-biphenyl]-3-yl)amino]sulfonyl]-4-hydroxy-benzoic acid, methyl ester

**MF:** C<sub>20</sub>H<sub>14</sub>ClF<sub>2</sub>NO<sub>5</sub>S

**FW:** 453.8

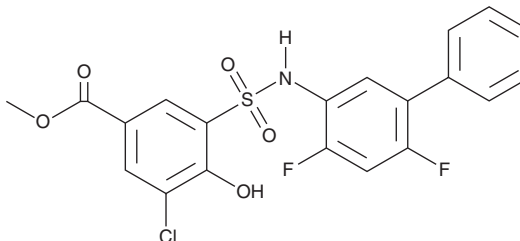
**Purity:** ≥98%

**UV/Vis.:** λ<sub>max</sub>: 292 nm

**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

NDI-091143 is supplied as a solid. A stock solution may be made by dissolving the NDI-091143 in the solvent of choice, which should be purged with an inert gas. NDI-091143 is soluble in the organic solvent DMSO at a concentration of approximately 1 mg/ml. NDI-091143 is slightly soluble in dimethyl formamide.

## Description

NDI-091143 is an allosteric inhibitor of ATP citrate lyase (ACL; K<sub>i</sub> = 7 nM).<sup>1</sup> It inhibits the production of ADP in a cell-free assay and oxaloacetate production in an enzyme-coupled assay (IC<sub>50</sub>s = 2.1 and 4.8 nM, respectively).

## Reference

1. Wei, J., Leit, S., Kuai, J., *et al.* An allosteric mechanism for potent inhibition of human ATP-citrate lyase. *Nature* **568(7753)**, 566-570 (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, 11/04/2022

## 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