

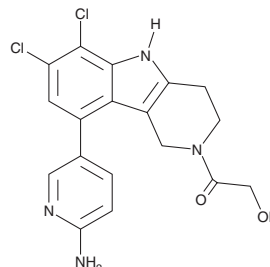
# PRODUCT INFORMATION



## G150

Item No. 28318

**CAS Registry No.:** 2369751-30-2  
**Formal Name:** 1-[9-(6-amino-3-pyridinyl)-6,7-dichloro-1,3,4,5-tetrahydro-2H-pyrido[4,3-b]indol-2-yl]-2-hydroxy-ethanone  
**MF:** C<sub>18</sub>H<sub>16</sub>Cl<sub>2</sub>N<sub>4</sub>O<sub>2</sub>  
**FW:** 391.3  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 234, 301 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

G150 is supplied as a crystalline solid. A stock solution may be made by dissolving the G150 in the solvent of choice, which should be purged with an inert gas. G150 is soluble in DMSO.

### Description

G150 is an inhibitor of cyclic GMP-AMP synthase (cGAS; IC<sub>50</sub> = 10.2 nM for the human enzyme).<sup>1</sup> It is selective for human cGAS over mouse cGAS (IC<sub>50</sub> = 25,000 nM). G150 inhibits dsDNA-induced expression of *IFNB1* in THP-1 cells and isolated human monocyte-derived macrophages (IC<sub>50</sub>s = 1.96 and 0.62 μM, respectively) but not LPS-induced expression of *TNF* or poly(I:C)-induced rRNA degradation in the same cells when used at a concentration of 10 μM.

### Reference

1. Lama, L., Adura, C., Xie, W., *et al.* Development of human cGAS-specific small-molecule inhibitors for repression of dsDNA-triggered interferon expression. *Nat. Commun.* **10**(1), 2261 (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, 12/13/2022

#### 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](http://WWW.CAYMANCHEM.COM)