# PRODUCT INFORMATION



H-151

Item No. 25857

CAS Registry No.: 941987-60-6

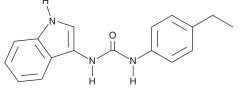
Formal Name: N-(4-ethylphenyl)-N'-1H-indol-3-yl-urea

MF:  $C_{17}H_{17}N_3O$ FW: 279.3 **Purity:** 

 $\lambda_{max}$ : 236, 258 nm UV/Vis.: 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**

H-151 is supplied as a crystalline solid. A stock solution may be made by dissolving the H-151 in the solvent of choice. H-151 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of H-151 in ethanol is approximately 10 mg/ml and approximately 30 mg/ml in DMSO and DMF.

H-151 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, H-151 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. H-151 has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

H-151 is an inhibitor of stimulator of interferon genes (STING).1 It forms an adduct with mouse STING in serum and acts in an irreversible manner. It inhibits the type I interferon (IFN) response and reduces the phosphorylation of TBK1 and palmitoylation of human STING (hsSTING) in vitro. It also reduces increases in mouse serum levels of IFN- $\beta$  and IL-6 induced by the STING agonist 10-carboxymethyl-9-acridanone (CMA).

### Reference

1. Haag, S.M., Gulen, M.F., Reymond, L., et al. Targeting STING with covalent small-molecule inhibitors. Nature 559(7713), 269-273 (2018).

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

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

uyer 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, 02/20/2020

### **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