

# PRODUCT INFORMATION

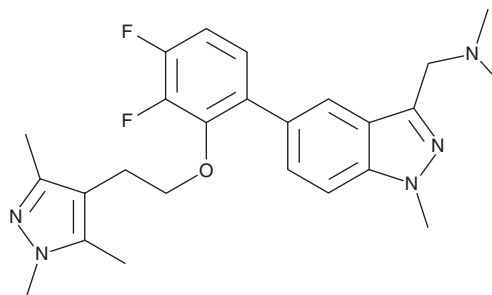


## IMP-1088

Item No. 25366

**CAS Registry No.:** 2059148-82-0  
**Formal Name:** 5-[3,4-difluoro-2-[2-(1,3,5-trimethyl-1H-pyrazol-4-yl)ethoxy]phenyl]-N,N,1-trimethyl-1H-indazole-3-methanamine

**MF:** C<sub>25</sub>H<sub>29</sub>F<sub>2</sub>N<sub>5</sub>O  
**FW:** 453.5  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 217, 244, 305 nm  
**Supplied as:** A solution in methyl acetate  
**Storage:** -20°C  
**Stability:** ≥2 years



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

### Description

IMP-1088 is an inhibitor of N-myristoyltransferase 1 (NMT1) and NMT2 (IC<sub>50</sub> = <1 nM for both human enzymes).<sup>1</sup> It inhibits replication of rhinovirus (RV), the virus most commonly responsible for the common cold, by inhibiting myristoylation of the RV capsid protein VP0 from strain RV-A16 in HeLa cells. IMP-1088 (500 nM) reduces RV-A16, RV-A1, RV-B14, and RV-A28 viral titers in HeLa cells without affecting cell growth (IC<sub>50</sub> = >1,000 nM). It also blocks production of infectious RV-16 virus in primary human bronchial epithelial cells (hBECs) when added 3 hours post infection *via* inhibition of viral capsid assembly.

### Reference

1. Mousnier, A., Bell, A.S., Swieboda, D.P., *et al.* Fragment-derived inhibitors of human N-myristoyltransferase block capsid assembly and replication of the common cold virus. *Nat. Chem.* **10**(6), 599-606 (2018).

#### 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, 01/14/2021

#### 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