

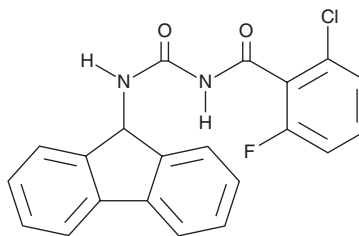
# PRODUCT INFORMATION



## TMN 355

Item No. 31155

CAS Registry No.: 1186372-20-2  
Formal Name: 2-chloro-N-[(9H-fluoren-9-ylamino) carbonyl]-6-fluoro-benzamide  
MF:  $C_{21}H_{14}ClFN_2O_2$   
FW: 380.8  
Purity:  $\geq 98\%$   
Supplied as: A solid  
Storage:  $-20^{\circ}\text{C}$   
Stability:  $\geq 4$  years



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

### Laboratory Procedures

TMN 355 is supplied as a solid. A stock solution may be made by dissolving the TMN 355 in the solvent of choice, which should be purged with an inert gas. TMN 355 is soluble in the organic solvent DMSO at a concentration of approximately 40 mM.

### Description

TMN 355 is an inhibitor of cyclophilin A ( $IC_{50} = 1.52$  nM).<sup>1</sup> It inhibits HIV-1 pseudoparticle infection of HEK293/T-17 cells when used at concentrations ranging from 0.1 to 1  $\mu\text{M}$ .<sup>2</sup>

### References

1. Ni, S., Yuan, Y., Huang, J., *et al.* Discovering potent small molecule inhibitors of cyclophilin A using de novo drug design approach. *J. Med. Chem.* **52**(17), 5295-5298 (2009).
2. Gale, T.V., Horton, T.M., Hoffmann, A.R., *et al.* Host proteins identified in extracellular viral particles as targets for broad-spectrum antiviral inhibitors. *J. Proteome Res.* **18**(1), 7-17 (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, 10/24/2022

#### 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