

# PRODUCT INFORMATION

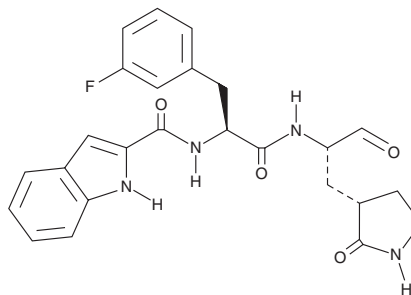


## M<sup>PRO</sup> Inhibitor 11b

Item No. 31345

**CAS Registry No.:** 2413716-71-7  
**Formal Name:** N-[(1S)-1-[(3-fluorophenyl)methyl]-2-[[[(1S)-1-formyl-2-[(3S)-2-oxo-3-pyrrolidinyl]ethyl]amino]-2-oxoethyl]-1H-indole-2-carboxamide

**MF:** C<sub>25</sub>H<sub>25</sub>FN<sub>4</sub>O<sub>4</sub>  
**FW:** 464.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 297 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

M<sup>PRO</sup> inhibitor 11b is supplied as a crystalline solid. A stock solution may be made by dissolving the M<sup>PRO</sup> inhibitor 11b in the solvent of choice, which should be purged with an inert gas. M<sup>PRO</sup> inhibitor 11b is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of M<sup>PRO</sup> inhibitor 11b in these solvents is approximately 30 mg/ml.

### Description

M<sup>PRO</sup> inhibitor 11b is an inhibitor of severe acute respiratory syndrome coronavirus 2 (SARS-CoV-2) main protease (M<sup>PRO</sup>; IC<sub>50</sub> = 0.04 μM in a TR-FRET assay).<sup>1</sup> It reduces viral yield in the culture supernatant of SARS-CoV-2-infected Vero E6 cells (EC<sub>50</sub> = 0.72 μM). M<sup>PRO</sup> inhibitor 11b also reduces viral RNA copy numbers in the same model when used at concentrations ranging from 1.85 to 50 μM.

### Reference

1. Dai, W., Zhang, B., Jiang, X.-M., *et al.* Structure-based design of antiviral drug candidates targeting the SARS-CoV-2 main protease. *Science* **368**(6497), 1331-1335 (2020).

#### 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, 03/21/2024

#### 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