

PRODUCT INFORMATION



PF-00835231

Item No. 36047

CAS Registry No.: 870153-29-0

Formal Name: N-[(1S)-1-[[[(1S)-3-hydroxy-2-oxo-1-[[[(3S)-2-oxo-3-pyrrolidinyl]methyl]propyl]amino]carbonyl]-3-methylbutyl]-4-methoxy-1H-indole-2-carboxamide

MF: C₂₄H₃₂N₄O₆

FW: 472.5

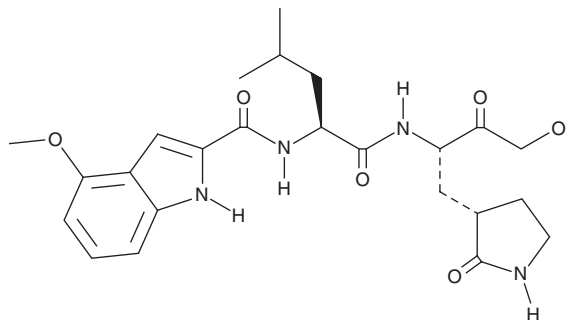
Purity: ≥98%

UV/Vis.: λ_{max}: 236, 292 nm

Supplied as: A 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

PF-00835231 is supplied as a solid. A stock solution may be made by dissolving the PF-00835231 in the solvent of choice, which should be purged with an inert gas. PF-00835231 is soluble in DMSO.

Description

PF-00835231 is an inhibitor of severe acute respiratory syndrome coronavirus (SARS-CoV) and SARS-CoV-2 main protease (M^{Pro}), also known as 3C-like protease (3CL^{Pro}; K_is = 4 and 0.27 nM, respectively).¹ It reduces the infectivity of SARS-CoV and common cold human coronavirus 229E (HCoV-229E) in Vero cells (EC₅₀s = 4.8 and 0.09 μM, respectively).

Reference

1. Hoffman, R.L., Kania, R.S., Brothers, M.A., *et al.* Discovery of ketone-based covalent inhibitors of coronavirus 3CL proteases for the potential therapeutic treatment of COVID-19. *J. Med. Chem.* **63**(21), 12725-12747 (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, 12/19/2022

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