

PRODUCT INFORMATION

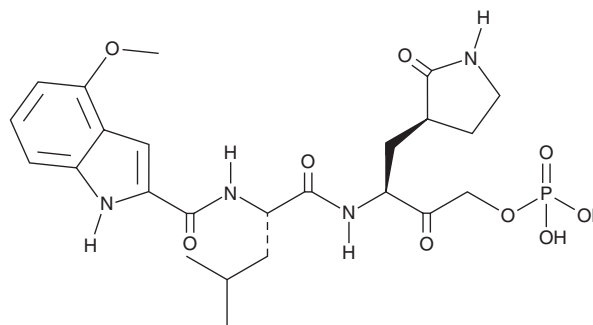


PF-07304814

Item No. 37026

CAS Registry No.: 2468015-78-1
Formal Name: 4-methoxy-N-[(1S)-3-methyl-1-[[[(1S)-2-oxo-1-[[[(3S)-2-oxo-3-pyrrolidinyl]methyl]-3-(phosphonoxy)propyl]amino]carbonyl]butyl]-1H-indole-2-carboxamide

Synonym: Lufotrelvir
MF: C₂₄H₃₃N₄O₉P
FW: 552.5
Purity: ≥98%
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-07304814 is supplied as a solid. A stock solution may be made by dissolving the PF-07304814 in the solvent of choice, which should be purged with an inert gas. PF-07304814 is soluble in DMSO.

Description

PF-07304814 is a prodrug form of the severe acute respiratory syndrome coronavirus (SARS-CoV) and SARS-CoV-2 main protease (M^{Pro}), also known as 3C-like protease (3CL^{Pro}), inhibitor PF-00835231 (Item No. 36047).¹

Reference

1. Boras, B., Jones, R.M., Anson, B.J., *et al.* Preclinical characterization of an intravenous coronavirus 3CL protease inhibitor for the potential treatment of COVID19. *Nat. Commun.* **12**(1), 6055 (2021).

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, 03/08/2024

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