

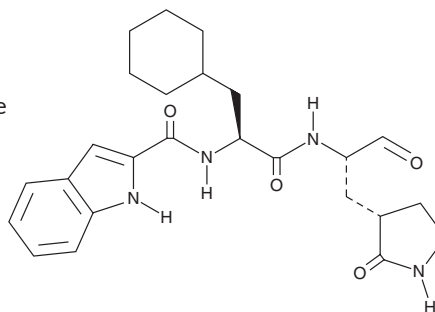
PRODUCT INFORMATION



Bofutrelvir

Item No. 31344

CAS Registry No.: 2103278-86-8
Formal Name: N-[(1S)-1-(cyclohexylmethyl)-2-[[[(1S)-1-formyl-2-[(3S)-2-oxo-3-pyrrolidinyl]ethyl]amino]-2-oxoethyl]-1H-indole-2-carboxamide
Synonym: FB2001
MF: C₂₅H₃₂N₄O₄
FW: 452.6
Purity: ≥95%
UV/Vis.: λ_{max}: 218, 296 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

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

Description

Bofutrelvir is an inhibitor of severe acute respiratory syndrome coronavirus 2 (SARS-CoV-2) main protease (M^{pro}; IC₅₀ = 0.053 μM in a TR-FRET assay).¹ It reduces viral yield in the culture supernatant of SARS-CoV-2-infected Vero E6 cells (EC₅₀ = 0.53 μM). Bofutrelvir 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