

PRODUCT INFORMATION

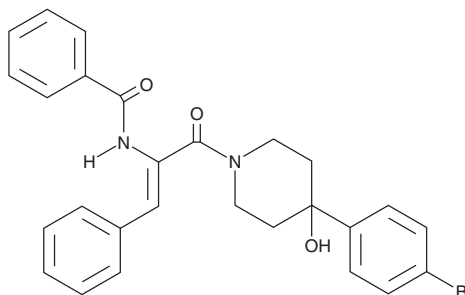


K22

Item No. 31578

CAS Registry No.: 2141978-86-9
Formal Name: N-[(1Z)-1-[[4-(4-bromophenyl)-4-hydroxy-1-piperidinyl]carbonyl]-2-phenylethenyl]-benzamide

MF: C₂₇H₂₅BrN₂O₃
FW: 505.4
Purity: ≥98%
UV/Vis.: λ_{max}: 221, 282 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

K22 is supplied as a crystalline solid. A stock solution may be made by dissolving the K22 in the solvent of choice, which should be purged with an inert gas. K22 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of K22 in these solvents is approximately 30 mg/ml.

Description

K22 is an antiviral agent.¹ It inhibits the replication of severe acute respiratory syndrome coronavirus (SARS-CoV), Middle East respiratory syndrome coronavirus (MERS-CoV), human coronavirus 229E (HCoV-229E), feline coronavirus (FCoV), and mouse hepatitis virus (MHV) in Vero cells when used at a concentration of 40 μM. K22 inhibits the formation of double membrane vesicles, a hallmark of coronavirus replication, and inhibits viral RNA synthesis in HCoV-229E-infected MRC-5 cells. It also reduces viral titers in cells infected with various *Arteriviridae* or *Torovirinae* viruses.²

References

1. Lundin, A., Dijkman, R., Bergström, T., *et al.* Targeting membrane-bound viral RNA synthesis reveals potent inhibition of diverse coronaviruses including the Middle East respiratory syndrome virus. *PLoS Pathog.* **10(5)**, e1004166 (2014).
2. Rappe, J.C.F., de Wilde, A., Di, H., *et al.* Antiviral activity of K22 against members of the order Nidovirales. *Virus Res.* **246**, 28-34 (2018).

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, 11/03/2020

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