

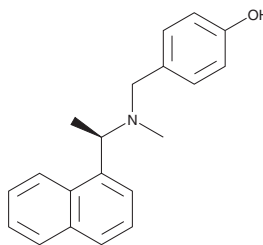
PRODUCT INFORMATION



Jun9-72-2

Item No. 37195

CAS Registry No.: 2657659-42-0
Formal Name: 4-[[methyl[(1R)-1-(1-naphthalenyl)ethyl]amino]methyl]-phenol
MF: C₂₀H₂₁NO
FW: 291.4
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

Jun9-72-2 is supplied as a solid. A stock solution may be made by dissolving the Jun9-72-2 in the solvent of choice, which should be purged with an inert gas. Jun9-72-2 is soluble in DMSO.

Description

Jun9-72-2 is an inhibitor of the severe acute respiratory syndrome coronavirus 2 (SARS-CoV-2) papain-like protease (PL^{pro}; IC₅₀ = 0.67 μM).¹ It inhibits viral replication of SARS-CoV-2 in Vero E6 and Caco2-hACE2 cells (EC₅₀s = 8.31 and 7.9 μM, respectively).

Reference

1. Ma, C., Sacco, M.D., Xia, Z., *et al.* Discovery of SARS-CoV-2 papain-like protease inhibitors through a combination of high-throughput screening and a FlipGFP-based reporter assay. *ACS Cent. Sci.* **7**(7), 1245-1260 (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, 10/17/2022

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