

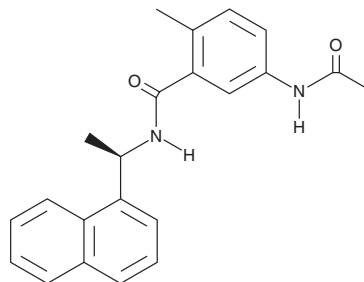
PRODUCT INFORMATION



HY-17542

Item No. 33354

CAS Registry No.: 1093070-14-4
Formal Name: 5-(acetylamino)-2-methyl-N-[(1R)-1-(1-naphthalenyl)ethyl]-benzamide
Synonym: PL^{pro} Inhibitor 6
MF: C₂₂H₂₂N₂O₂
FW: 346.4
Purity: ≥98%
UV/Vis.: λ_{max}: 224 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

HY-17542 is supplied as a crystalline solid. A stock solution may be made by dissolving the HY-17542 in the solvent of choice, which should be purged with an inert gas. HY-17542 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of HY-17542 in these solvents is approximately 30 mg/ml.

Description

HY-17542 is an inhibitor of the severe acute respiratory syndrome coronavirus (SARS-CoV) papain-like protease (PL^{pro}; IC₅₀ = 2.6 μM).¹ It increases the viability of SARS-CoV-infected Vero E6 cells (EC₅₀ = 13.1 μM).

Reference

1. Ratia, K., Pegan, S., Takayama, J., *et al.* A noncovalent class of papain-like protease/deubiquitinase inhibitors blocks SARS virus replication. *Proc. Natl. Acad. Sci. USA* **105**(42), 16119-16124 (2008).

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, 10/12/2022

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