

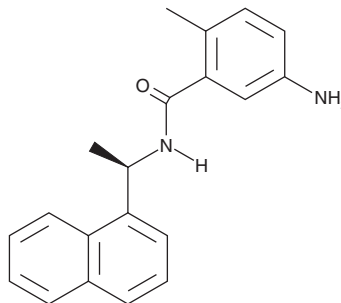
PRODUCT INFORMATION



GRL-0617

Item No. 33348

CAS Registry No.: 1093070-16-6
Formal Name: 5-amino-2-methyl-N-[(1R)-1-(1-naphthalenyl)ethyl]-benzamide
MF: C₂₀H₂₀N₂O
FW: 304.4
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

GRL-0617 is supplied as a crystalline solid. A stock solution may be made by dissolving the GRL-0617 in the solvent of choice, which should be purged with an inert gas. GRL-0617 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of GRL-0617 in these solvents is approximately 30 mg/ml.

Description

GRL-0617 is an inhibitor of severe acute respiratory syndrome coronavirus (SARS-CoV) and SARS-CoV-2 papain-like proteases (PL^{Pro}s; IC₅₀s = 0.6 and 1.5 μM, respectively).^{1,2} It increases the viability of SARS-CoV-infected Vero E6 cells (EC₅₀ = 14.5 μM).¹ GRL-0617 (6.25-50 μM) reduces viral replication in SARS-CoV-2-infected Caco-2 cells.²

References

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).
2. Shin, D., Mukherjee, R., Grewe, D., *et al.* Papain-like protease regulates SARS-CoV-2 viral spread and innate immunity. *Nature* **587**(7835), 657-662 (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, 10/26/2022

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