

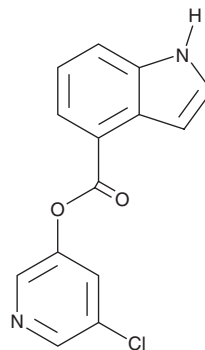
PRODUCT INFORMATION



GRL-0496

Item No. 36114

CAS Registry No.: 1087243-14-8
Formal Name: 1H-indole-4-carboxylic acid,
5-chloro-3-pyridinyl ester
MF: C₁₄H₉ClN₂O₂
FW: 272.7
Purity: ≥98%
UV/Vis.: λ_{max}: 212, 230 nm
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

GRL-0496 is supplied as a solid. A stock solution may be made by dissolving the GRL-0496 in the solvent of choice, which should be purged with an inert gas. GRL-0496 is soluble in organic solvents such as ethanol and DMSO. The solubility of GRL-0496 in these solvents is approximately 25 and 30 mg/ml, respectively.

Description

GRL-0496 is an inhibitor of severe acute respiratory syndrome coronavirus (SARS-CoV) and SARS-CoV-2 main protease (M^{Pro}), also known as 3C-like protease (3CL^{Pro}; IC₅₀s = 0.03 and 0.038 μM, respectively).^{1,2} It reduces viral infectivity in SARS-CoV-infected Vero cells (EC₅₀ = 6.9 μM).¹

References

1. Ghosh, A.K., Gong, G., Grum-Tokars, V., *et al.* Design, synthesis and antiviral efficacy of a series of potent chloropyridyl ester-derived SARS-CoV 3CL^{Pro} inhibitors. *Bioorg. Med. Chem. Lett.* **18(20)**, 5684-5688 (2008).
2. Cao, W., Cho, C.-C.D., Geng, Z.Z., *et al.* Evaluation of SARS-CoV-2 main protease inhibitors using a novel cell-based assay. *ACS Cent. Sci.* **8(2)**, 192-204 (2022).

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, 12/19/2022

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