

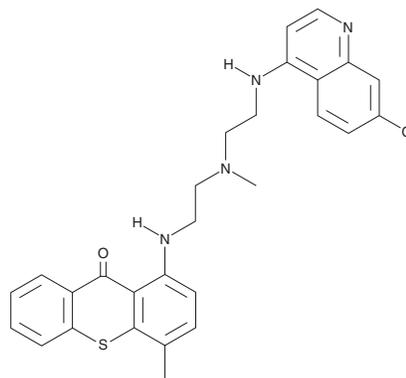
PRODUCT INFORMATION



ROC-325

Item No. 33566

CAS Registry No.: 1859141-26-6
Formal Name: 1-[[2-[[2-[(7-chloro-4-quinolinyl)amino]ethyl]methylamino]ethyl]amino]-4-methyl-9H-thioxanthen-9-one
MF: C₂₈H₂₇ClN₄OS
FW: 503.1
Purity: ≥98%
UV/Vis.: λ_{max}: 220, 257 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

ROC-325 is supplied as a solid. A stock solution may be made by dissolving the ROC-325 in the solvent of choice, which should be purged with an inert gas. ROC-325 is soluble in the organic solvent chloroform at a concentration of approximately 10 mg/ml. ROC-325 is also slightly soluble in ethanol, DMSO, and dimethyl formamide.

Description

ROC-325 is an autophagy inhibitor.¹ It inhibits autophagic flux and induces lysosomal deacidification and apoptosis in 786-O renal cell carcinoma cells when used at a concentration of 5 μM. ROC-325 decreases viability in a panel of 12 cancer cell lines, including renal, lung, pancreatic, and prostate cancer cells, with IC₅₀ values ranging from 4.9 to 11 μM. It reduces tumor growth in a 786-O mouse xenograft model when administered at a dose of 50 mg/kg. ROC-325 also decreases the cytopathic effect of severe acute respiratory syndrome coronavirus 2 (SARS-CoV-2) in infected Vero E6 cells (EC₅₀ = 3.28 μM).²

References

- Carew, J.S., Espitia, C.M., Zhao, W., *et al.* Disruption of autophagic degradation with ROC-325 antagonizes renal cell carcinoma pathogenesis. *Clin. Cancer Res.* **23(11)**, 2869-2879 (2017).
- Gorshkov, K., Chen, C.Z., Bostwick, R., *et al.* The SARS-CoV-2 cytopathic effect is blocked by lysosome alkalinizing small molecules. *ACS Infect. Dis.* (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, 12/06/2022

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