

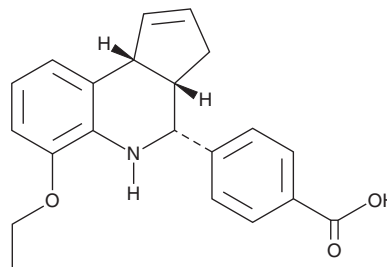
PRODUCT INFORMATION



Lin28 Inhibitor LI71

Item No. 41039

CAS Registry No.: 1357248-83-9
Formal Name: 4-[(3aS,4R,9bR)-6-ethoxy-3a,4,5,9b-tetrahydro-3H-cyclopenta[c]quinolin-4-yl]-benzoic acid
MF: C₂₁H₂₁NO₃
FW: 335.4
Purity: ≥95%
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

Lin28 inhibitor LI71 is supplied as a solid. A stock solution may be made by dissolving the Lin28 inhibitor LI71 in the solvent of choice, which should be purged with an inert gas. Lin28 inhibitor LI71 is sparingly soluble (1-10 mg/ml) in DMSO.

Description

Lin28 inhibitor LI71 is an inhibitor of the protein-microRNA precursor interaction between the RNA-binding protein Lin28 and *let-7*.¹ It binds to the cold-shock domain of Lin28 to inhibit *let-7* binding and inhibits Lin28-mediated *let-7* oligouridylation (IC₅₀ = ~27 μM). Lin28 inhibitor LI71 (100 μM) increases the levels of mature *let-7b*, *-7c*, *-7f*, *-7g*, and *-7i* as well as the levels of miR-16, which is not regulated by Lin28, in K562 leukemia cells but not the levels of miR-93, miR-99a, and miR-151, which are also not regulated by Lin28. It increases the levels of mature *let-7a-i* without increasing miR-16 levels in Lin28a/Lin28b double-knockout mouse embryonic stem cells (mESCs) expressing exogenous wild-type Lin28a. Lin28 inhibitor LI71 also inhibits the protein-protein interaction between *P. falciparum* cold-shock protein (PfCoSP) and PfDNA or Pfa/β-tubulin.² It is active against *P. falciparum* strain 3D7 and the chloroquine-resistant strain RKL-9 (IC₅₀s = 1.35 and 6.3 nM, respectively).

References

1. Wang, L., Rowe, R.G., Jaimes, A., *et al.* Small-molecule inhibitors disrupt *let-7* oligouridylation and release the selective blockade of *let-7* processing by LIN28. *Cell Rep.* **23**(10), 3091-3101 (2018).
2. Behl, A., Shoaib, R., De Leon, F., *et al.* Targeting an essential Plasmodium cold shock protein to block growth and transmission of malaria parasite. *iScience* **26**(5), 106637 (2023).

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, 06/03/2024

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