

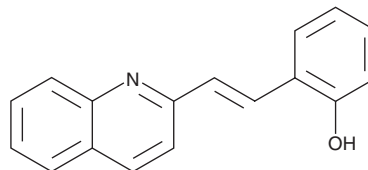
PRODUCT INFORMATION



Quininib

Item No. 19838

CAS Registry No.: 143816-42-6
Formal Name: 2-[(1E)-2-(2-quinolinyl)ethenyl]-phenol
MF: C₁₇H₁₃NO
FW: 247.3
Purity: ≥98%
UV/Vis.: λ_{max}: 225, 282, 359 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

Quininib is supplied as a crystalline solid. A stock solution may be made by dissolving the quinininib in the solvent of choice, which should be purged with an inert gas. Quininib is soluble in DMSO.

Description

Quininib is an antagonist of the cysteinyl leukotriene 1 (CysLT₁) and CysLT₂ receptors (IC₅₀s = 1.4 and 52 μM, respectively).¹ It inhibits tubule formation in HMEC-1 cells when used at concentrations of 3.16 and 10 μM and inhibits sprout formation in isolated mouse aortic rings in an *ex vivo* model of angiogenesis when used at 10 μM. Quininib (3 μM) decreases angiogenesis in an oxygen-induced retinopathy mouse model of ocular angiogenesis, preventing revascularization when used at concentrations of 0.5 and 3 μM, as well as increasing neovascularization and vascular density when used at 3 μM. It reduces tumor growth and inhibits angiogenesis in an HT-29 colorectal cancer mouse xenograft model when administered at a dose of 50 mg/kg every three days.²

References

1. Reynolds, A.L., Alvarez, Y., Sasore, T., *et al.* Phenotype-based discovery of 2-[(E)-2-(quinolin-2-yl)viny] phenol as a novel regulator of ocular angiogenesis. *J. Biol. Chem.* **291(14)**, 7242-7255 (2016).
2. Murphy, A.G., Casey, R., Maguire, A., *et al.* Preclinical validation of the small molecule drug quinininib as a novel therapeutic for colorectal cancer. *Sci. Rep.* **6**, 34523 (2016).

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, 11/21/2022

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