

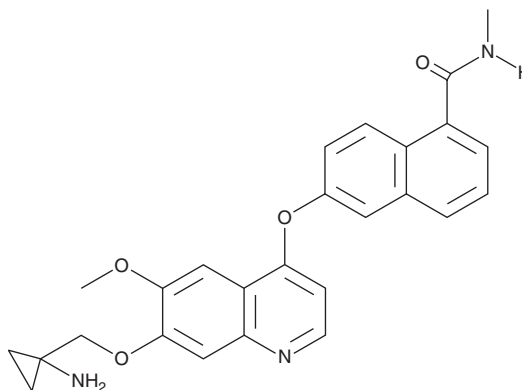
PRODUCT INFORMATION



Lucitanib

Item No. 41711

CAS Registry No.: 1058137-23-7
Formal Name: 6-[[7-[(1-aminocyclopropyl)methoxy]-6-methoxy-4-quinolinyl]oxy]-N-methyl-1-naphthalenecarboxamide
Synonyms: AL 3810, E-3810
MF: C₂₆H₂₅N₃O₄
FW: 443.5
Purity: ≥98%
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

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

Description

Lucitanib is a dual inhibitor of VEGFR and FGFR that inhibits VEGFR1, VEGFR2, and VEGFR3 (IC₅₀s = 7, 25, and 10 nM, respectively) and FGFR1 and FGFR2 (IC₅₀s = 17.5 and 82.5 nM, respectively).¹ It is selective for these receptors over FGFR3 and FGFR4 (IC₅₀s = 237.5 and >1,000 nM, respectively), as well as the PDGFR family members c-Kit, PDGFRα, and PDGFRβ (IC₅₀s = 456, 175, and 525 nM, respectively), but does not inhibit FMS, also known as CSF-1 receptor tyrosine kinase (IC₅₀ = 5 nM). Lucitanib inhibits the proliferation of human umbilical vein endothelial cells (HUVECs) induced by VEGF or basic FGF (bFGF; IC₅₀s = 40 and 50 nM, respectively) but inhibits the proliferation of A2780 ovarian, A498 renal, and SK-N-SH neuroblastoma cells only at higher concentrations (IC₅₀s = 27.7, 19.6, and 12.79 μM, respectively). It reduces tumor growth in HT-29 colon and A2780 ovarian cancer mouse xenograft models, as well as in RXF 393, A498, and SK-N-SH neuroblastoma mouse xenograft models in a dose-dependent manner. It also completely inhibits bFGF-induced angiogenesis in a Matrigel™ plug assay in mice when administered at a dose of 20 mg/kg.

Reference

1. Bello, E., Colella, G., Scarlato, V., et al. E-3810 is a potent dual inhibitor of VEGFR and FGFR that exerts antitumor activity in multiple preclinical models. *Cancer Res.* **71**(4), 1396-1405 (2011).

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, 09/11/2024

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