

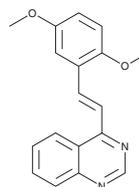
PRODUCT INFORMATION



K882

Item No. 43422

Formal Name: (E)-4-(2,5-dimethoxystyryl)quinazoline
MF: C₁₈H₁₆N₂O₂
FW: 292.3
Purity: ≥95%
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

K882 is supplied as a crystalline solid. A stock solution may be made by dissolving the K882 in the solvent of choice, which should be purged with an inert gas. K882 is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

Description

K882 is an inhibitor of Src kinase.¹ It binds to Src ($K_d = 315$ nM), decreases phosphorylation at the tyrosine in position 416 (Y416), a Src activation site, and increases phosphorylation at Y527, an inactivation site. K882 reduces the proliferation of wild-type NCI H358 non-small cell lung cancer cells and NCI H358 cells expressing a KRAS^{G12C} mutation (IC_{50} s = 2.14 and 2.03 μ M, respectively). It also inhibits migration and invasion of NCI H358 cells when used at concentrations of 2, 4, and 8 μ M, induces cell cycle arrest in the G₂/M phase at 1, 4, and 8 μ M, and induces apoptosis at 4 and 8 μ M. K882 (25, 50, and 100 mg/kg) reduces tumor volume in an NCI H358 mouse xenograft model.

Reference

1. Ma, X., Ma, Z., Qi, X., *et al.* Identification of a novel Src inhibitor K882 derived from quinazoline-based stilbenes with anti-NSCLC effect. *Bioorg. Chem.* **156**, 108185 (2025).

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, 05/14/2025

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