

PRODUCT INFORMATION



L1Au2

Item No. 43244

CAS Registry No.: 3038386-42-1

Formal Name: [3-[[4-[(3-chloro-4-fluorophenyl)amino]-7-methoxy-6-quinazolinyl]oxy]-1-propyn-1-yl] (1,3,5-triaza-7-phosphatricyclo[3.3.1.1^{3,7}]decane-κP⁷)-gold

Synonyms: Alkynyl-Au(I) Gefitinib, N-(3-Chloro-4-fluorophenyl)-7-methoxy-6-(prop-2-yn-1-yloxy)-quinazolin-4-amine(1,3,5-triaza-7-phosphaadamantane)gold(I), TrxR/EGFR-IN-1

MF: C₂₄H₂₄AuClFN₆O₂P

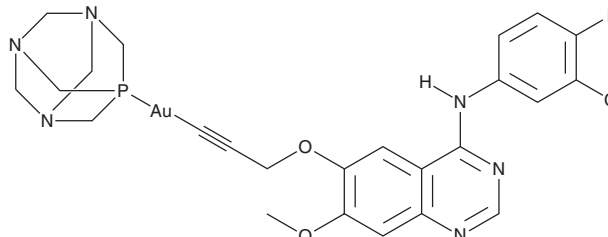
FW: 710.9

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

L1Au2 is supplied as a solid. A stock solution may be made by dissolving the L1Au2 in the solvent of choice, which should be purged with an inert gas. L1Au2 is slightly soluble (0.1-1 mg/ml) in DMSO.

Description

L1Au2 is a dual inhibitor of EGFR and thioredoxin reductase (TrxR; IC₅₀ = 0.99 μM).¹ It inhibits the phosphorylation of EGFR, Akt, and ERK in PC-9 lung cancer cells, which express EGFR containing the exon 19 deletion, when used at concentrations of 0.01 or 0.05 μM. L1Au2 inhibits the proliferation of PC-9 and gefitinib-resistant PC-9 cells (IC₅₀s = 0.012 and 1.76 μM, respectively) and induces ferroptosis in gefitinib-resistant PC-9 cells. *In vivo*, L1Au2 (10 mg/kg) decreases tumor weight and intratumoral glutathione peroxidase 4 (Gpx4) levels in a gefitinib-resistant PC-9 mouse xenograft model.

Reference

1. Wang, L., Ma, X., Zhou, L., et al. Dual-targeting TrxR-EGFR alkynyl-Au(I) gefitinib complex induces ferroptosis in gefitinib-resistant lung cancer via degradation of GPX4. *J. Med. Chem.* **68**(5), 5275-5291 (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, 04/28/2025

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