

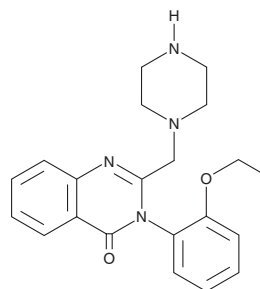
PRODUCT INFORMATION



PRLX-93936

Item No. 25631

CAS Registry No.: 903499-49-0
Formal Name: 3-(2-ethoxyphenyl)-2-(1-piperazinylmethyl)-4(3H)-quinazolinone
MF: $C_{21}H_{24}N_4O_2$
FW: 364.4
Purity: $\geq 98\%$
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

PRLX-93936 is supplied as a solid. A stock solution may be made by dissolving the PRLX-93936 in the solvent of choice which should be purged with an inert gas. PRLX-93936 is soluble in DMSO.

Description

PRLX-93936 is an analog of erastin (Item No. 17754) that has antitumor activity.¹ It inhibits the hypoxia-inducible factor 1 (HIF-1) signaling pathway under hypoxic conditions ($IC_{50} = 0.09 \mu\text{M}$ in a cell-based reporter assay).² PRLX-93936 ($1 \mu\text{M}$) also inhibits hypoxia-induced increases in HIF-1 α expression in ME-180 cervical cancer cells. It inhibits the growth of HT-1080 fibrosarcoma, OVCAR-5 ovarian cancer, BJELR tumorigenic primary fibroblast, and PANC-1 pancreatic cancer cells with IC_{50} values of less than 100 nM .¹ PRLX-93936 inhibits tumor growth in PANC-1 and HT-1080 xenograft models when administered at a dose of 50 mg/kg and induces tumor regression when administered at a dose of 100 mg/kg .

References

1. Stockwell, B.R. Erastin analogs and uses thereof. **US20070161644A1** (2007).
2. Huang, W., Huang, R., Attene Ramos, M., *et al.* Synthesis and evaluation of quinazolin-4-ones as hypoxia-inducible factor-1 α inhibitors. *Bioorg. Med. Chem. Lett.* **21(18)**, 5239-5243 (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, 10/27/2022

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