

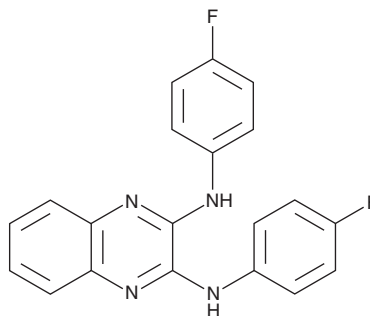
PRODUCT INFORMATION



LQZ-7I

Item No. 36602

CAS Registry No.: 195822-23-2
Formal Name: N²,N³-bis(4-fluorophenyl)-2,3-
quinoxalinediamine
MF: C₂₀H₁₄F₂N₄
FW: 348.4
Purity: ≥98%
UV/Vis.: λ_{max}: 244, 270 nm
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

LQZ-7I is supplied as a solid. A stock solution may be made by dissolving the LQZ-7I in the solvent of choice, which should be purged with an inert gas. LQZ-7I is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of LQZ-7I in these solvents is approximately 1, 2, and 5 mg/ml, respectively.

Description

LQZ-7I is an inhibitor of survivin.¹ It inhibits survivin dimerization in a secreted alkaline phosphatase (SEAP) assay using DU145 prostate cancer cells expressing survivin and induces survivin degradation in C4-2 and PC3 prostate cancer cells in a proteasome-dependent manner. LQZ-7I is cytotoxic to C4-2 and PC3 cells (IC₅₀s = 3.1 and 4.8 μM, respectively). It reduces tumor growth in a PC3 mouse xenograft model when administered at a dose of 100 mg/kg.

Reference

1. Peery, R., Kyei-Baffour, K., Dong, Z., *et al.* Synthesis and identification of a novel lead targeting survivin dimerization for proteasome-dependent degradation. *J. Med. Chem.* **63**(13), 7243-7251 (2020).

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, 12/20/2022

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