

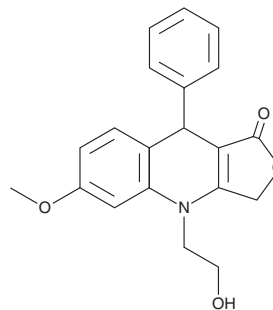
PRODUCT INFORMATION



NSC 756093

Item No. 35023

CAS Registry No.: 1629908-92-4
Formal Name: 4,9-dihydro-4-(2-hydroxyethyl)-6-methoxy-9-phenyl-furo[3,4-b]quinolin-1(3H)-one
MF: C₂₀H₁₉NO₄
FW: 337.4
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

NSC 756093 is supplied as a solid. A stock solution may be made by dissolving the NSC 756093 in the solvent of choice, which should be purged with an inert gas. NSC 756093 is soluble in the organic solvent DMSO.

Description

NSC 756093 is an inhibitor of the protein-protein interaction between guanylate binding protein 1 (GBP1) and the serine/threonine kinase Pim-1.¹ It binds to GBP1-Pim-1 ($K_d = 38$ nM) and inhibits the interaction by 65% when used at a concentration of 100 nM. NSC 756093 enhances radiation-induced cytotoxicity in FaDu hypopharyngeal carcinoma cells ($IC_{50} = 496$ nM), as well as induces apoptosis and cell cycle arrest at the G₂/M phase in the same cells.²

References

1. Andreoli, M., Persico, M., Kumar, A., *et al.* Identification of the first inhibitor of the GBP1:PIM1 interaction. Implications for the development of a new class of anticancer agents against paclitaxel resistant cancer cells. *J. Med. Chem.* **57**(19), 7916-7932 (2014).
2. Resendez, A., Tailor, D., Graves, E., *et al.* Radiosensitization of head and neck squamous cell carcinoma (HNSCC) by a podophyllotoxin. *ACS Med. Chem. Lett.* **10**(9), 1314-1321 (2019).

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, 12/02/2022

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