

PRODUCT INFORMATION



NSC 668394

Item No. 16863

CAS Registry No.: 382605-72-3

Formal Name: 7-[[2-(3,5-dibromo-4-hydroxyphenyl)ethyl]amino]-5,8-quinolinedione

Synonym: UA 668394

MF: C₁₇H₁₂Br₂N₂O₃

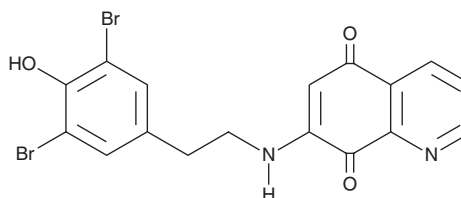
FW: 452.1

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 668394 is supplied as a solid. A stock solution may be made by dissolving the NSC 668394 in the solvent of choice, which should be purged with an inert gas. NSC 668394 is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

NSC 668394 is a binder of ezrin ($K_d = 10.6 \mu\text{M}$).¹ It inhibits PKC α -induced phosphorylation of ezrin in a cell-free assay ($\text{IC}_{50} = 8.1 \mu\text{M}$).² NSC 668394 (10 μM) inhibits the interaction between ezrin and actin in K7M2 osteosarcoma cells. It inhibits the migration and invasion of K7M2 cancer cells when used at a concentration of 2 μM .¹ NSC 668394 (10 μM) reduces bacterial invasion but is not bactericidal in HEp-2 cells infected with group A *Streptococcus* (GAS).³ It induces developmental delays and cyclopia in *Xenopus* larvae when used at a concentration of 10 μM and induces toxicity at 40 μM .²

References

1. Paige, M., Kosturko, G., Bulut, G., *et al.* Design, synthesis and biological evaluation of ezrin inhibitors targeting metastatic osteosarcoma. *Bioorg. Med. Chem.* **22**(1), 478-487 (2014).
2. Bulut, G., Hong, S.-H., Chen, K., *et al.* Small molecule inhibitors of ezrin inhibit the invasive phenotype of osteosarcoma cells. *Oncogene* **31**(3), 269-281 (2021).
3. Rox, K., Rohde, M., Chhatwal, G.S., *et al.* Disorazoles block group A streptococcal invasion into epithelial cells via interference with the host factor ezrin. *Cell Chem. Biol.* **24**(2), 159-170 (2017).

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, 09/30/2024

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