

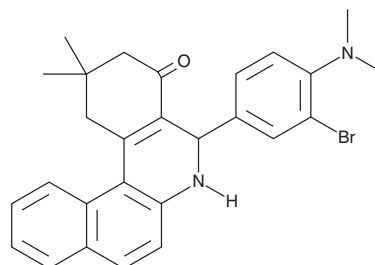
PRODUCT INFORMATION



Glutaminase Inhibitor Compound 968

Item No. 17199

CAS Registry No.: 311795-38-7
Formal Name: 5-[3-bromo-4-(dimethylamino)phenyl]-2,3,5,6-tetrahydro-2,2-dimethyl-benzo[a]phenanthridin-4(1H)-one
MF: C₂₇H₂₇BrN₂O
FW: 475.4
Purity: ≥98%
UV/Vis.: λ_{max}: 216, 232, 281, 292, 339, 371 nm
Supplied as: A crystalline 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

Glutaminase inhibitor compound 968 is supplied as a crystalline solid. A stock solution may be made by dissolving the glutaminase inhibitor compound 968 in the solvent of choice, which should be purged with an inert gas. Glutaminase inhibitor compound 968 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of glutaminase inhibitor compound 968 in these solvents is approximately 0.5, 14, and 16 mg/ml, respectively.

Glutaminase inhibitor compound 968 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, glutaminase inhibitor compound 968 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Glutaminase inhibitor compound 968 has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Glutaminase inhibitor compound 968 is a cell-permeable, reversible inhibitor of a mitochondrial glutaminase splice variant, glutaminase C (IC₅₀ = ~2.5 μM), which is commonly found in cancer cells.^{1,2} Through this action, glutaminase inhibitor compound 968 blocks transformation induced by Rho GTPases and inhibits the proliferation of non-invasive epithelial (T-47D and MDA-MB-361) and invasive mesenchymal (MDA-MB-231 and Hs578T) cancer cells, without affecting normal cells.²⁻⁴ Glutaminase inhibitor compound 968 induces apoptosis and alters both histone methylation and gene expression in cancer cells.⁴

References

1. Stalneck, C. A., Ulrich, S. M., Li, Y., *et al.* Mechanism by which a recently discovered allosteric inhibitor blocks glutamine metabolism in transformed cells. *Proc. Natl. Acad. Sci. USA* **112**(2), 394-399 (2015).
2. Wang, J. B., Erickson, J. W., Fuji, R., *et al.* Targeting mitochondrial glutaminase activity inhibits oncogenic transformation. *Cancer Cell* **18**(3), 2007-219 (2010).
3. Katt, W. P., Ramachandran, S., Erickson, J. W., *et al.* Dibenzophenanthridines as inhibitors of glutaminase C and cancer cell proliferation. *Mol. Cancer Ther.* **11**(6), 1269-1278 (2012).
4. Simpson, N. E., Tryndyak, V. P., Pogribna, M., *et al.* Modifying metabolically sensitive histone marks by inhibiting glutamine metabolism affects gene expression and alters cancer cell phenotype. *Epigenetics* **7**(12), 1413-1420 (2012).

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, 10/13/2022

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