

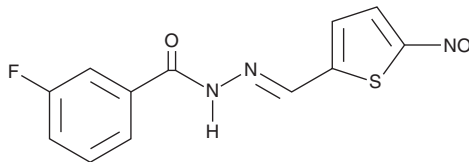
PRODUCT INFORMATION



MitoBloCK-10

Item No. 42054

CAS Registry No.: 394694-98-5
Formal Name: 3-fluoro-benzoic acid 2-[(5-nitro-2-thienyl)methylene]hydrazide
Synonym: MB-10
MF: C₁₂H₈FN₃O₃S
FW: 293.3
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

MitoBloCK-10 is supplied as a solid. A stock solution may be made by dissolving the MitoBloCK-10 in the solvent of choice, which should be purged with an inert gas. MitoBloCK-10 is soluble (≥10 mg/ml) in the organic solvent DMSO.

Description

MitoBloCK-10 is an inhibitor of mitochondrial protein import.¹ It inhibits the import of cytochrome c1 into the mitochondrial intermembrane space in yeast cells when used at a concentration of 100 μM. MitoBloCK-10 (25 μM) inhibits migration and invasion of, and induces apoptosis and cell cycle arrest at the G₁ phase in, primary human bladder cancer cells.² It induces mitochondrial depolarization, decreases ATP levels, and increases reactive oxygen species (ROS) in primary human bladder cancer cells. MitoBloCK-10 (20 mg/kg every other day) decreases tumor volume in a patient-derived xenograft (PDX) mouse model of bladder cancer. Intravitreal administration of MitoBloCK-10 (0.25 nM/eye) reduces superoxide dismutase (SOD) activity and increases TBARS levels and acellular capillary numbers in mouse retina.³ It induces dorsal curvature and cardiac apoptosis in zebrafish embryos when used at a concentration of 10 μM.¹

References

1. Miyata, N., Tang, Z., Conti, M.A., *et al.* Adaptation of a genetic screen reveals an inhibitor for mitochondrial protein import component Tim44. *J. Biol. Chem.* **292**(13), 5429-5442 (2017).
2. Zhang, L., Shi, X., Zhang, L., *et al.* A first-in-class TIMM44 blocker inhibits bladder cancer cell growth. *Cell Death Dis.* **15**(3), 204 (2024).
3. Ma, Z.-R., Li, h.-P., Cai, S.-Z., *et al.* The mitochondrial protein TIMM44 is required for angiogenesis in vitro and in vivo. *Cell Death Dis.* **14**(5), 307 (2023).

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, 12/09/2024

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