

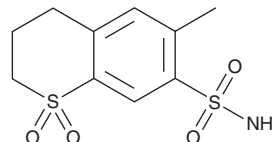
PRODUCT INFORMATION



Meticrane

Item No. 44226

CAS Registry No.: 1084-65-7
Formal Name: 3,4-dihydro-6-methyl-2H-1-benzothiopyran-7-sulfonamide, 1,1-dioxide
MF: C₁₀H₁₃NO₄S₂
FW: 275.3
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

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

Description

Meticrane is a diuretic.¹ It decreases the viability of Jurkat and K562 leukemia and SK-HEP-1 liver cancer cells when used at a concentration of 1 mM.² Meticrane (0.125 mM) enhances cell death induced by 5-azacytidine (Item No. 11164) or CUDC-101 (Item No. 16426) in the same cells. Formulations containing meticrane have been used in the treatment of hypertension.

References

1. Tanaka, R. and Hirayama, N. Crystal structure of meticrane. *Anal. Sci.* **23**, 119-120 (2007).
2. Wang, Y., Sharma, A., Ge, F., et al. Non-oncology drug (meticrane) shows anti-cancer ability in synergy with epigenetic inhibitors and appears to be involved passively in targeting cancer cells. *Front. Oncol.* **13**, 1157366 (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, 10/07/2025

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