

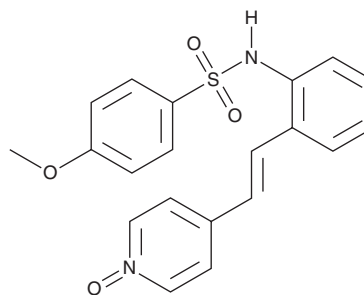
PRODUCT INFORMATION



HMN-176

Item No. 37424

CAS Registry No.: 173529-10-7
Formal Name: 4-methoxy-N-[2-[(1E)-2-(1-oxido-4-pyridinyl)ethenyl]phenyl]-benzenesulfonamide
MF: C₂₀H₁₈N₂O₄S
FW: 382.4
Purity: ≥98%
UV/Vis.: λ_{max}: 239, 343 nm
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

HMN-176 is supplied as a solid. A stock solution may be made by dissolving the HMN-176 in the solvent of choice, which should be purged with an inert gas. HMN-176 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of HMN-176 in these solvents is approximately 10 mg/ml.

Description

HMN-176 is an anticancer agent and active metabolite of the prodrug HMN-214 (Item No. 26210).¹ It is cytotoxic against a panel of cancer cell lines with a mean IC₅₀ value of 112 nM, as well as P388 leukemia cells that are resistant to the DNA-crosslinking agent cisplatin (Item No. 13119), anthracycline antitumor antibiotic doxorubicin (Item No. 15007), or antimetabolic vincristine (Item No. 11764) (IC₅₀s = 143, 557, and 265 nM, respectively). HMN-176 (3 μM) induces cell cycle arrest at the G₂/M phase in HeLa cervical cancer cells.

References

1. Takagi, M., Honmura, T., Watanabe, S., *et al.* *In vivo* antitumor activity of a novel sulfonamide, HMN-214, against human tumor xenografts in mice and the spectrum of cytotoxicity of its active metabolite, HMN-176. *Invest. New Drugs* **21**(4), 387-399 (2003).

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, 09/27/2023

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