

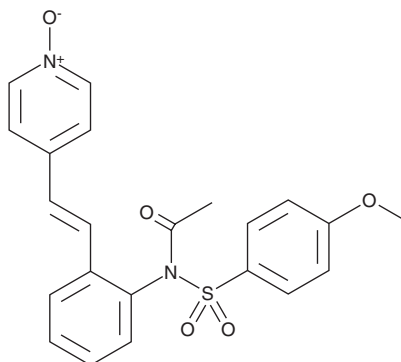
PRODUCT INFORMATION



HMN-214

Item No. 26210

CAS Registry No.: 173529-46-9
Formal Name: N-[(4-methoxyphenyl)sulfonyl]-
N-[2-[(1E)-2-(1-oxido-4-pyridinyl)
ethenyl]phenyl]-acetamide
Synonym: IVX-214
MF: C₂₂H₂₀N₂O₅S
FW: 424.5
Purity: ≥98%
UV/Vis.: λ_{max}: 244, 340 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

HMN-214 is supplied as a crystalline solid. A stock solution may be made by dissolving the HMN-214 in the solvent of choice, which should be purged with an inert gas. HMN-214 is soluble in the organic solvent DMSO at a concentration of approximately 12 mg/ml.

Description

HMN-214 is an orally bioavailable prodrug form of HMN-176, an indirect inhibitor of polo-like kinase (PLK) activity that inhibits proliferation of a variety of cancer cells.^{1,2} HMN-214 decreases the expression of multidrug resistance gene 1 (MDR1) in AB-A.1 cells and in tumors isolated from mice bearing multidrug-resistant KB-A1 xenografts.² HMN-214 (20 mg/kg per day) reduces tumor volume in PC3, WiDr, and A549 mouse xenograft models.¹ It does not decrease nerve conduction velocity or compound action potential amplitude in rabbit sciatic nerves *in vivo* when administered at a concentration of 30 mg/kg per day.

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).
2. Tanaka, H., Ohshima, N., Ikenoya, M., *et al.* HMN-176, an active metabolite of the synthetic antitumor agent HMN-214, restores chemosensitivity to multidrug-resistant cells by targeting the transcription factor NF- κ B. *Cancer Res.* **63**(20), 6942-6947 (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, 11/04/2022

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