

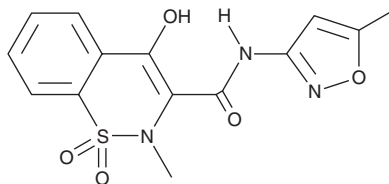
PRODUCT INFORMATION



Isoxicam

Item No. 41340

CAS Registry No.: 34552-84-6
Formal Name: 4-hydroxy-2-methyl-N-(5-methyl-3-isoxazolyl)-12H-1,2b-enzothiazine-3-carboxamide, 1-dioxide
Synonyms: W-8495
MF: C₁₄H₁₃N₃O₅S
FW: 335.3
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

Isoxicam is supplied as a solid. A stock solution may be made by dissolving the isoxicam in the solvent of choice, which should be purged with an inert gas. Isoxicam is slightly soluble (0.1-1 mg/ml) in acetonitrile and sparingly soluble (1-10 mg/ml) in DMSO.

Description

Isoxicam is a non-steroidal anti-inflammatory drug (NSAID).¹ It inhibits the production of prostaglandin E₂ (PGE₂; Item No. 14010) in HSDM1C1 mouse fibrosarcoma cells (IC₅₀ = 2 μM).² Isoxicam (12.5, 25, and 50 mg/kg) reduces carrageenan-induced paw edema in rats.¹ It also reduces hind paw weight in a rat model of adjuvant-induced arthritis when administered at doses of 1.56, 6.25, or 25 mg/kg per day.

References

1. DiPasquale, G., Rassaert, C., Richter, R., *et al.* The anti-inflammatory properties of isoxicam (4-hydroxy-2-methyl-N-[5-Methyl-3-isoxolyl-2H-1,2-benzothiazine-3-carboxamide 1,1-dioxide]). *Agents Actions* **5(3)**, 256-263 (1975).
2. Pugsley, T.A., Spencer, C., Bactor, A.M., *et al.* Selective inhibition of the cyclooxygenase pathway of the arachidonic acid cascade by the nonsteroidal antiinflammatory drug isoxicam *Drug Develop. Res.* **5**, 171-178 (1985).

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, 08/27/2024

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