

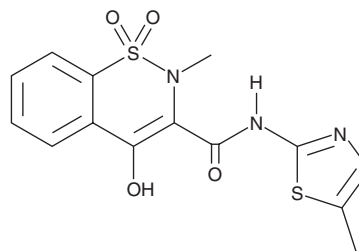
PRODUCT INFORMATION



Meloxicam

Item No. 40657

CAS Registry No.: 71125-38-7
Formal Name: 4-hydroxy-2-methyl-N-(5-methyl-2-thiazolyl)-2H-1,2-benzothiazine-3-carboxamide 1,1-dioxide
Synonym: UH-AC 62XX
MF: C₁₄H₁₃N₃O₄S₂
FW: 351.4
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.

Description

Meloxicam (Item No. 40657) is an analytical reference standard categorized as a non-steroidal anti-inflammatory drug (NSAID) and COX-2 inhibitor.¹ Meloxicam has been found as an adulterant in illicit samples distributed as food in capsule form.² Formulations containing meloxicam have been used in the treatment of osteoarthritis and rheumatoid arthritis. This product is intended for analytical forensic applications. This product is also available as a general research tool (Item No. 14906).

References

1. Ogino, K., Hatanaka, K., Kawamura, M., *et al.* Evaluation of pharmacological profile of meloxicam as an anti-inflammatory agent, with particular reference to its relative selectivity for cyclooxygenase-2 over cyclooxygenase-1. *Pharmacology* **55**, 44-53 (1997).
2. Park, H.-J., Ham, H.J., Yang, Y.J., *et al.* Liquid chromatography-quadrupole orbitrap and liquid chromatography-tandem mass spectrometry system for rapid identification and quantitation of thirty nonsteroidal anti-inflammatory drugs and acetaminophen in illegal products. *Rapid Commun. Mass Spectrom.* **37**(6), e9462 (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, 04/12/2024

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