

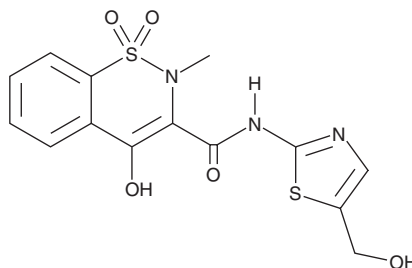
PRODUCT INFORMATION



5'-hydroxy Meloxicam

Item No. 22159

CAS Registry No.: 130262-92-9
Formal Name: 4-hydroxy-N-[5-(hydroxymethyl)-2-thiazolyl]-2-methyl-2H-1,2-benzothiazine-3-carboxamide, 1,1-dioxide
MF: C₁₄H₁₃N₃O₅S₂
FW: 367.4
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.

Laboratory Procedures

5'-hydroxy Meloxicam is supplied as a solid. A stock solution may be made by dissolving the 5'-hydroxy meloxicam in the solvent of choice, which should be purged with an inert gas. 5'-hydroxy Meloxicam is slightly soluble in DMSO and methanol.

Description

5'-hydroxy Meloxicam is a metabolite of the non-steroidal anti-inflammatory drug (NSAID) meloxicam (Item No. 14906).¹ It is formed *via* metabolism of meloxicam primarily by the cytochrome P450 (CYP) isoform CYP2C9 and to a lesser extent by CYP3A4.

References

1. Chesné, C., Guyomard, C., Guillouzo, A., *et al.* Metabolism of meloxicam in human liver involves cytochromes P450C9 and 3A4. *Xenobiotica* **28(1)**, (1998).

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, 12/22/2022

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