

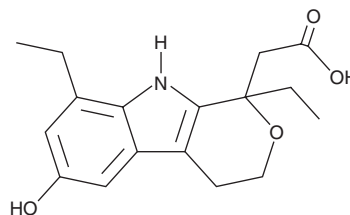
PRODUCT INFORMATION



6-hydroxy Etodolac

Item No. 34211

CAS Registry No.: 101901-06-8
Formal Name: 1,8-diethyl-1,3,4,9-tetrahydro-6-hydroxy-pyrano[3,4-b]indole-1-acetic acid
MF: C₁₇H₂₁NO₄
FW: 303.4
Purity: ≥90%
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

6-hydroxy Etodolac is supplied as a solid. A stock solution may be made by dissolving the 6-hydroxy etodolac in the solvent of choice, which should be purged with an inert gas. 6-hydroxy Etodolac is soluble in organic solvents such as DMSO and methanol.

Description

6-hydroxy Etodolac is a metabolite of the non-steroidal anti-inflammatory drug (NSAID) and COX inhibitor etodolac (Item No. 20833).¹ 6-hydroxy Etodolac induces false-positive reactions in diazo diagnostic tests for bilirubin in urine.

References

1. Ferdinandi, E.S., Sehgal, S.N., Demerson, C.A., *et al.* Disposition and biotransformation of ¹⁴C-etodolac in man. *Xenobiotica* **16**(2), 153-166 (1986).
2. Sho, Y., Ishiodori, T., Oketani, M., *et al.* Effects of urinary metabolites of etodolac on diagnostic tests of bilirubin in urine. *Arzneimittelforschung* **49**(7), 572-576 (1999).

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, 11/18/2022

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