

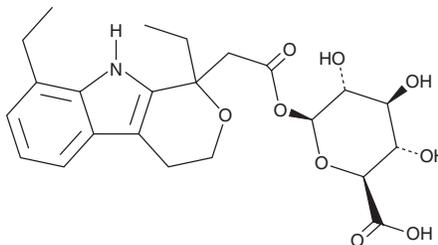
PRODUCT INFORMATION



Etodolac Acyl Glucuronide

Item No. 34072

CAS Registry No.: 79541-43-8
Formal Name: β -D-glucopyranuronic acid, 1-(1,8-diethyl-1,3,4,9-tetrahydropyrano[3,4-b]indole-1-acetate)
MF: $C_{23}H_{29}NO_9$
FW: 463.5
Purity: $\geq 90\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Etodolac acyl glucuronide is supplied as a solid. A stock solution may be made by dissolving the etodolac acyl glucuronide in the solvent of choice, which should be purged with an inert gas. Etodolac acyl glucuronide is soluble in methanol.

Description

Etodolac acyl glucuronide is a phase II metabolite of the non-steroidal anti-inflammatory drug (NSAID) and COX inhibitor etodolac (Item No. 20833).¹ It is formed *via* glucuronidation of etodolac by the UDP-glucuronosyltransferase (UGT) isoforms UGT1A9, UGT1A10, and UGT2B7.

Reference

1. Kutsuno, Y., Itoh, T., Tukey, R.H., *et al.* Glucuronidation of drugs and drug-induced toxicity in humanized UDP-glucuronosyltransferase 1 mice. *Drug Metab. Dispos.* **42(7)**, 1146-1152 (2014).

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, 12/05/2022

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