

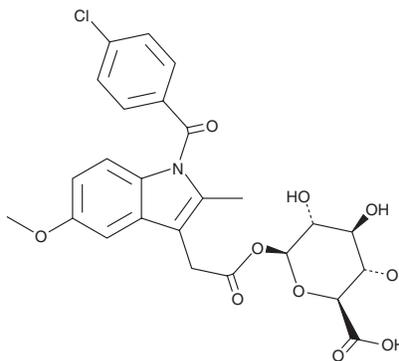
PRODUCT INFORMATION



Indomethacin Acyl Glucuronide

Item No. 34579

CAS Registry No.: 75523-11-4
Formal Name: β -D-glucopyranuronic acid, 1-[1-(4-chlorobenzoyl)-5-methoxy-2-methyl-1H-indole-3-acetate]
Synonym: Indomethacin Acyl- β -D-Glucuronide
MF: C₂₅H₂₄ClNO₁₀
FW: 533.9
Purity: \geq 90%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

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

Description

Indomethacin acyl glucuronide is a metabolite of the COX inhibitor indomethacin (Item No. 70270).¹

Reference

1. Moolenaar, F., Crancrinus, S., Visser, J., *et al.* Clearance of indomethacin occurs predominantly by renal glucuronidation. *Pharm. Weekbl. Sci.* **14(4)**, 191-195 (1992).

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, 10/10/2022

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