

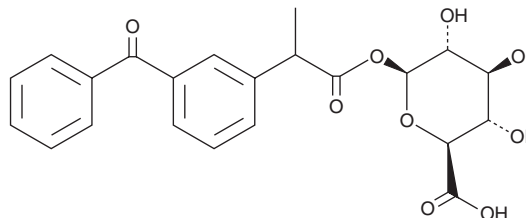
PRODUCT INFORMATION



(±)-Ketoprofen Glucuronide

Item No. 27180

CAS Registry No.: 76690-94-3
Formal Name: 1-(3-benzoyl- α -methylbenzeneacetate)
 β -D-glucopyranuronic acid
Synonyms: *rac*-Ketoprofen Acyl- β -D-glucuronide,
(R,S)-Ketoprofen Glucuronide
MF: C₂₂H₂₂O₉
FW: 430.4
Purity: \geq 95%
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

(±)-Ketoprofen glucuronide is supplied as a solid. A stock solution may be made by dissolving the (±)-ketoprofen glucuronide in the solvent of choice. (±)-Ketoprofen glucuronide is soluble in organic solvents such as methanol and DMSO, which should be purged with an inert gas.

Description

(±)-Ketoprofen glucuronide is a phase II metabolite of the non-steroidal anti-inflammatory drug (NSAID) ketoprofen (Item No. 10006661).¹ It is formed *via* glucuronidation by the UDP-glucuronosyltransferase (UGT) isoforms UGT1A3, UGT1A9, UGT1A10, and UGT2B7.

Reference

1. Sabolovic, N., Heydel, J.-M., Li, X., *et al.* Carboxyl nonsteroidal anti-inflammatory drugs are efficiently glucuronidated by microsomes of the human gastrointestinal tract. *Biochim. Biophys. Acta.* **1675(1-3)**, 120-129 (2004).

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, 12/19/2022

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