

PRODUCT INFORMATION



Cipecofol Glucuronide

Item No. 42168

Formal Name: (2S,3S,4S,5R,6S)-6-(2-((R)-1-cyclopropylethyl)-6-isopropylphenoxy)-3,4,5-trihydroxytetrahydro-2H-pyran-2-carboxylic acid

Synonym: Ciprofol Glucuronide

MF: C₂₀H₂₈O₇

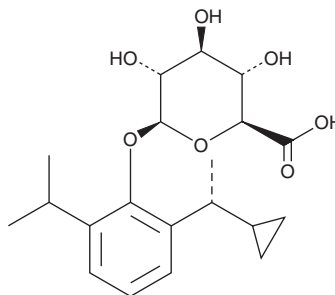
FW: 380.4

Purity: ≥95%

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

Cipecofol glucuronide is supplied as a solid. A stock solution may be made by dissolving the cipecofol glucuronide in the solvent of choice, which should be purged with an inert gas. Cipecofol glucuronide is slightly soluble (0.1-1 mg/ml) in methanol.

Description

Cipecofol glucuronide is a metabolite of cipecofol (Item No. 40951).¹ It is formed from cipecofol by the UDP-glucuronosyltransferase (UGT) isoform UGT1A9.

Reference

1. Hou, L., Zhao, Y., Zhao, S., *et al.* Ciprofol is primarily glucuronidated by UGT1A9 and predicted not to cause drug-drug interactions with typical substrates of CYP1A2, CYP2B6, and CYP2C19. *Chem. Biol. Interact.* **387**, 110811 (2024).

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, 01/23/2025

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