

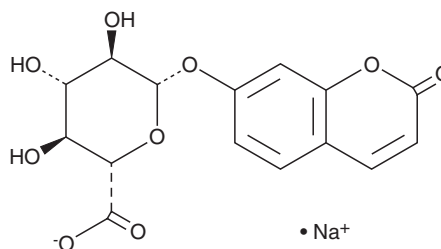
PRODUCT INFORMATION



7-hydroxy Coumarin Glucuronide (sodium salt)

Item No. 19882

CAS Registry No.: 168286-98-4
Formal Name: 2-oxo-2H-1-benzopyran-7-yl- β -D-glucopyranosiduronic acid, monosodium salt
MF: $C_{15}H_{13}O_9 \cdot Na$
FW: 360.2
Purity: $\geq 95\%$
UV/Vis.: λ_{max} : 316 nm
Supplied as: A crystalline solid
Storage: $-20^{\circ}C$
Stability: ≥ 2 years
Item Origin: Synthetic



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

Laboratory Procedures

7-hydroxy Coumarin glucuronide (sodium salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the 7-hydroxy coumarin glucuronide (sodium salt) in the solvent of choice, which should be purged with an inert gas. 7-hydroxy Coumarin glucuronide (sodium salt) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 7-hydroxy coumarin glucuronide (sodium salt) in these solvents is approximately 0.2, 16, and 14 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of 7-hydroxy coumarin glucuronide (sodium salt) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of 7-hydroxy coumarin glucuronide (sodium salt) in PBS (pH 7.2) is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

7-hydroxy Coumarin glucuronide is a benzopyrone and a metabolite of 7-hydroxy coumarin.¹ It is formed from 7-hydroxy coumarin by the UDP-glucuronosyltransferase (UGT) isoform UGT1A1.

Reference

1. Negoro, R., Tasaka, M., Deguchi, S., *et al.* Generation of HepG2 cells with high expression of multiple drug-metabolizing enzymes for drug discovery research using a PiTCh system. *Cells* **11(10)**, 1677 (2022).

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, 04/17/2023

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