

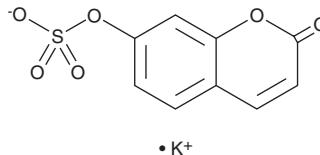
PRODUCT INFORMATION



7-hydroxy Coumarin sulfate (potassium salt)

Item No. 20892

CAS Registry No.: 1135316-80-1
Formal Name: 7-(sulfooxy)-2H-1-benzopyran-2-one, monopotassium salt
Synonyms: 7-HC, 7HCS
MF: C₉H₅O₆S • K
FW: 280.3
Purity: ≥98%
Supplied as: A crystalline 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

7-hydroxy Coumarin sulfate (potassium salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the 7-hydroxy coumarin sulfate (potassium salt) in water. 7-hydroxy coumarin sulfate (potassium salt) is slightly soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

7-hydroxy Coumarin sulfate is a phase II metabolite of coumarin that can be used as an internal standard for the analysis of 7-hydroxy coumarin metabolism using GC- or LC-MS.¹⁻³

References

1. Ekins, S., Williams, J.A., Murray, G.I., *et al.* Xenobiotic metabolism in rat, dog, and human precision-cut liver slices, freshly isolated hepatocytes, and vitrified precision-cut liver slices. *Drug Metab. Dispos.* **24**(9), 990-995 (1996).
2. Wang, Q., Jia, R., Ye, C., *et al.* Glucuronidation and sulfation of 7-hydroxycoumarin in liver matrices from human, dog, monkey, rat, and mouse. *In Vitro Cell. Dev. Biol. Anim.* **41**(3), 97-103 (2005).
3. Ho, M.D., Ring, N., Amaral, K., *et al.* Human enterocytes as an *in vitro* model for the evaluation of intestinal drug metabolism: characterization of drug-metabolizing enzyme activities of cryopreserved human enterocytes from twenty-four donors. *Drug Metab. Dispos.* **45**(6), 686-691 (2017).

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, 04/10/2025

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