

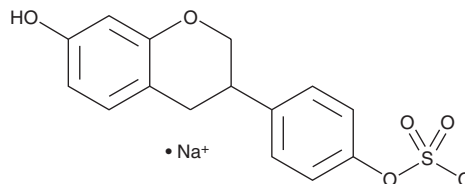
PRODUCT INFORMATION



(±)-Equol 4'-sulfate (sodium salt)

Item No. 21440

CAS Registry No.:	1189685-28-6
Formal Name:	3,4-dihydro-3-[4-(sulfooxy)phenyl]-2H-1-benzopyran-7-ol, monosodium salt
Synonym:	R,S-Equol 4'-Sulfate
MF:	C ₁₅ H ₁₃ O ₆ S • Na
FW:	344.3
Purity:	≥90%
Supplied as:	A solid
Storage:	-20°C
Stability:	≥4 years
Item Origin:	Synthetic



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

Laboratory Procedures

(±)-Equol 4'-sulfate (sodium salt) is supplied as a solid. A stock solution may be made by dissolving the (±)-equol 4'-sulfate (sodium salt) in the solvent of choice, which should be purged with an inert gas. (±)-Equol 4'-sulfate (sodium salt) is slightly soluble in methanol.

(±)-Equol 4'-sulfate (sodium salt) is slightly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

(±)-Equol 4'-sulfate is a metabolite of the isoflavonoid phytoestrogen (±)-equol (Item No. 13184).¹ It is produced from equol in the gut. (±)-Equol 4'-sulfate has been used as a standard for the quantification of unconjugated equol in rat plasma.²

References

1. Lee, D.-H., Kim, M.J., Lee, S.H., *et al.* Nutrikinetics of isoflavone metabolites after fermented soybean product (Cheonggukjang) ingestion in ovariectomized mice. *Mol. Nutr. Food Res.* **61(12)**, 1700322 (2017).
2. Gavina, J.M., Priem, J., Wood, C.M., *et al.* Determination of isoflavones in rat serum using liquid chromatography-tandem mass spectrometry with a highly efficient core-shell column. *Anal. Bioanal. Chem.* **405(8)**, 2643-2651 (2013).

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, 12/22/2022

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