

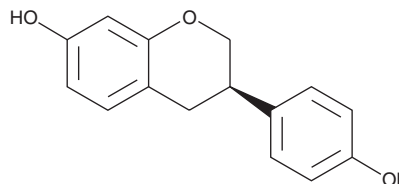
PRODUCT INFORMATION



(R)-Equol

Item No. 10010172

CAS Registry No.:	221054-79-1
Formal Name:	3,4-dihydro-3R-(4-hydroxyphenyl)-2H-1-benzopyran-7-ol
Synonyms:	(+)-Equol, Isoequol
MF:	C ₁₅ H ₁₄ O ₃
FW:	242.3
Purity:	≥98%
UV/Vis.:	λ _{max} : 225, 282 nm
Supplied as:	A crystalline 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

(R)-Equol is supplied as a crystalline solid. A stock solution may be made by dissolving the (R)-equol in the solvent of choice, which should be purged with an inert gas. (R)-Equol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of (R)-equol in ethanol and DMSO is approximately 20 mg/ml and approximately 10 mg/ml in DMF.

(R)-Equol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, (R)-equol should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. (R)-Equol has a solubility of approximately 0.1 mg/ml in a 1:10 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

(R)-Equol is an agonist of estrogen receptor α (ER α) and active metabolite of the polyketide isoflavone daidzein (Item No. 10005166).^{1,2} It is formed from daidzein by gut microbiota.¹ (R)-Equol selectively induces ER α -mediated gene transcription over ER β -mediated transcription in HEC-1 cells expressing the human receptors in a reporter assay (EC₅₀s = 66 and 330 nM, respectively).²

References

1. Setchell, K.D.R., Clerici, C., Lephart, E.D., *et al.* S-equol, a potent ligand for estrogen receptor β , is the exclusive enantiomeric form of the soy isoflavone metabolite produced by human intestinal bacterial flora. *Am. J. Clin. Nutr.* **81**(5), 1072-1079 (2005).
2. Muthyala, R.S., Ju, Y.H., Sheng, S., *et al.* Equol, a natural estrogenic metabolite from soy isoflavones: Convenient preparation and resolution of R- and S-equols and their differing binding and biological activity through estrogen receptors alpha and beta. *Bioorg. Med. Chem.* **12**(6), 1559-1567 (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, 01/26/2024

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