

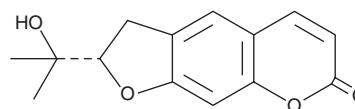
PRODUCT INFORMATION



Marmesin

Item No. 25049

CAS Registry No.:	13849-08-6
Formal Name:	(2S)-2,3-dihydro-2-(1-hydroxy-1-methylethyl)-7H-furo[3,2-g][1]benzopyran-7-one
Synonyms:	(+)-Marmesin, (S)-Marmesin, NSC 340840
MF:	C ₁₄ H ₁₄ O ₄
FW:	246.3
Purity:	≥98%
UV/Vis.:	λ _{max} : 224, 336 nm
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

Marmesin is supplied as a crystalline solid. A stock solution may be made by dissolving the marmesin in the solvent of choice. Marmesin is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of marmesin in these solvents is approximately 2, 50, and 30 mg/ml, respectively.

Marmesin is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, marmesin should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Marmesin has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Marmesin is a coumarin originally isolated from the mature bark of *A. marmelos*.¹ Marmesin decreases proliferation and cell invasion induced by fetal bovine serum (FBS) in H1299 non-small cell lung cancer cells (NSCLCs) when used at a concentration of 10 μM.² It decreases VEGF secretion in A549 and H1299 cells and inhibits capillary-like structure formation by human umbilical vein endothelial cells (HUVECs). Marmesin also decreases proliferation of U937 human leukemia cells (IC₅₀ = 40 μM).³ It halts the cell cycle at the G₂/M phase, increases the ratio of Bax to Bcl-2 protein, induces apoptosis, and inhibits cell migration. Marmesin reduces tumor growth in a U937 mouse xenograft model when administered at a dose of 30 mg/kg.

References

1. Chatterjee, A. and Mitra, S.S. On the constitution of the active principles isolated from the mature bark of *Aegle marmelos*, Correã. *J. Am. Chem. Soc.* **71**(2), 606-609 (1949).
2. Kim, J.H., Kim, M.S., Lee, B.H., et al. Marmesin-mediated suppression of VEGF/VEGFR and integrin β1 expression: Its implication in non-small cell lung cancer cell responses and tumor angiogenesis. *Oncol. Rep.* **37**(1), 91-97 (2017).
3. Dong, L., Xu, W.W., Li, H., et al. *In vitro* and *in vivo* anticancer effects of marmesin in U937 human leukemia cells are mediated via mitochondrial-mediated apoptosis, cell cycle arrest, and inhibition of cancer cell migration. *Oncol. Rep.* **39**(2), 597-602 (2018).

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, 12/19/2022

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