

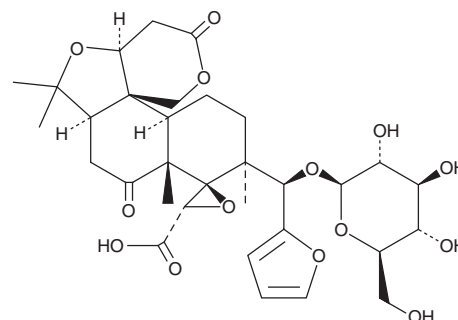
PRODUCT INFORMATION



Limonin 17-β-D-Glucopyranoside

Item No. 39147

CAS Registry No.: 123564-61-4
Formal Name: (2aR,3'S,4aS,5R,6S,8aR,8bR,12aS)-6-[(S)-3-furanyl(β-D-glucopyranosyloxy)methyl]decahydro-2,2,4a,6-tetramethyl-4,11-dioxo-spiro[9H,11H-naphtho[1',2':3,4]furo[3,2-c]pyran-5(2H),2'-oxirane]-3'-carboxylic acid
Synonym: Limonin Glucoside
MF: C₃₂H₄₂O₁₄
FW: 650.7
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Natural/Source unknown



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

Laboratory Procedures

Limonin 17-β-D-glucopyranoside is supplied as a solid. A stock solution may be made by dissolving the limonin 17-β-D-glucopyranoside in the solvent of choice, which should be purged with an inert gas. Limonin 17-β-D-glucopyranoside is soluble in water and methanol.

Description

Limonin 17-β-D-glucopyranoside is a furanolactone that has been found in *C. reticulata* and has anticancer activity.^{1,2} It induces apoptosis and disrupts the mitochondrial membrane potential in SW480 colon adenocarcinoma cells when used at a concentration of 100 μM.² Limonin 17-β-D-glucopyranoside inhibits the proliferation of MDA-Panc-28 pancreatic adenocarcinoma cells (IC₅₀ = 20.49 μM).¹

References

1. Patil, J.R., Jayaprakasha, G.K., Murthy, K.N.C., *et al.* Characterization of *Citrus aurantifolia* bioactive compounds and their inhibition of human pancreatic cancer cells through apoptosis. *Microchem. J.* **94**(2), 108-117 (2010).
2. Murthy, K.N.C., Jayaprakash, G.K., Kumar, V., *et al.* Citrus limonin and its glucoside inhibit colon adenocarcinoma cell proliferation through apoptosis. *J. Agric. Food Chem.* **59**(6), 2314-2323 (2011).

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, 07/14/2023

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