

# PRODUCT INFORMATION



## Limonin

Item No. 16932

**CAS Registry No.:** 1180-71-8  
**Formal Name:** (2aR,4aR,4bR,5aS,8S,8aS,10aR,10bR,14aS)-8-(3-furanyl)decahydro-2,2,4a,8a-tetramethyl-11H,13H-oxireno[d]pyrano[4',3':3,3a]isobenzofuro[5,4-f][2]benzopyran-4,6,13(2H,5aH)-trione di- $\delta$ -lactone Limonoic Acid

**Synonym:**

**MF:** C<sub>26</sub>H<sub>30</sub>O<sub>8</sub>

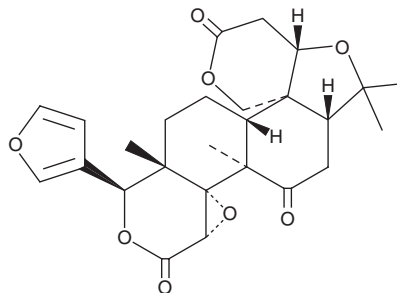
**FW:** 470.5

**Purity:**  $\geq$ 98%

**Supplied as:** A crystalline solid

**Storage:** -20°C

**Stability:**  $\geq$ 4 years



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

### Laboratory Procedures

Limonin is supplied as a crystalline solid. A stock solution may be made by dissolving the limonin in the solvent of choice. Limonin is soluble in organic solvents such as DMSO and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of limonin in these solvents is approximately 10 and 25 mg/ml, respectively.

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

### Description

Limonin is the first described limonoid, which are naturally occurring tetranortriterpenoids that impart bitter flavor in citrus fruits. Compounds belonging to this group have reported anti-proliferative and proapoptotic activity on several cancer cell lines as well as pesticidal, anti-malarial, anti-microbial, and anti-HIV activity.<sup>1</sup>

### Reference

1. Roy, A. and Saraf, S. Limonoids: Overview of significant bioactive triterpenes distributed in plants kingdom. *Biol. Pharm. Bull.* **29(2)**, 191-201 (2006).

#### 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, 11/30/2022

#### 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