

# PRODUCT INFORMATION



## Chonglou Saponin VII

Item No. 39251

CAS Registry No.: 68124-04-9

Formal Name: (3 $\beta$ ,25R)-17-hydroxyspirost-5-en-3-yl O-6-deoxy- $\alpha$ -L-mannopyranosyl-(1 $\rightarrow$ 2)-O-[O-6-deoxy- $\alpha$ -L-mannopyranosyl-(1 $\rightarrow$ 4)]- $\beta$ -D-glucopyranoside

Synonyms: Paris Saponin VII, PS VII

MF: C<sub>51</sub>H<sub>82</sub>O<sub>21</sub>

FW: 1,031.2

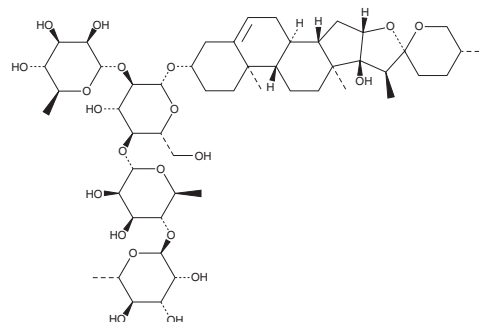
Purity:  $\geq$ 98%

Supplied as: A solid

Storage: -20°C

Stability:  $\geq$ 4 years

Item Origin: Plant/Paris polyphylla



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

### Laboratory Procedures

Chonglou saponin VII is supplied as a solid. A stock solution may be made by dissolving the chonglou saponin VII in the solvent of choice, which should be purged with an inert gas. Chonglou saponin VII is soluble in methanol.

### Description

Chonglou saponin VII is a steroid saponin that has been found in *P. polyphylla* and has anticancer activities.<sup>1</sup> It also inhibits the proliferation of MDA-MB-231, MDA-MB-436, and MCF-7 breast cancer cells (IC<sub>50</sub>s = 3.16, 3.45, and 2.86  $\mu$ M, respectively). It induces apoptosis and autophagosome formation in the same cells. Chonglou saponin VII (1.5 mg/kg) induces intratumoral autophagy and reduces tumor growth in an MDA-MB-231 mouse xenograft model.

### Reference

1. Xiang, Y.-C., Peng, P., Liu, X.-W., *et al.* Paris saponin VII, a Hippo pathway activator, induces autophagy and exhibits therapeutic potential against human breast cancer cells. *Acta Pharmacol. Sin.* **43**(6), 1568-1580 (2022).

#### 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, 06/18/2025

#### 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