

# PRODUCT INFORMATION



## Saikosaponin C

Item No. 34599

**CAS Registry No.:** 20736-08-7  
**Formal Name:** (3 $\beta$ ,16 $\beta$ )-13,28-epoxy-16-hydroxyolean-11-en-3-yl O-6-deoxy- $\alpha$ -L-mannopyranosyl-(1 $\rightarrow$ 4)-O- $[\beta$ -D-glucopyranosyl-(1 $\rightarrow$ 6)]- $\beta$ -D-glucopyranoside

**MF:** C<sub>48</sub>H<sub>78</sub>O<sub>17</sub>

**FW:** 927.1

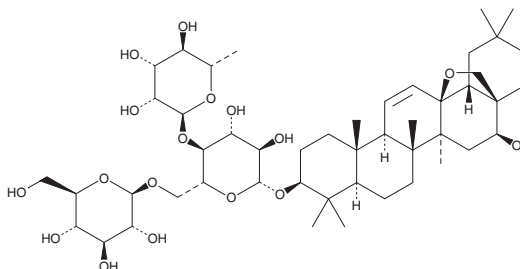
**Purity:**  $\geq$ 98%

**Supplied as:** A solid

**Storage:** -20°C

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

**Item Origin:** Plant/*Radix Bupleuri*



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

### Laboratory Procedures

Saikosaponin C is supplied as a solid. A stock solution may be made by dissolving the saikosaponin C in the solvent of choice, which should be purged with an inert gas. Saikosaponin C is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of saikosaponin C in ethanol is approximately 1 mg/ml and approximately 10 mg/ml in DMSO and DMF.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of saikosaponin C can be prepared by directly dissolving the solid in aqueous buffers. The solubility of saikosaponin C in PBS (pH 7.2) is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Saikosaponin C is a saikosaponin that has been found in *B. falactum* and has anti-inflammatory and analgesic activities.<sup>1,2</sup> It reduces basal production of prostaglandin E<sub>2</sub> (PGE<sub>2</sub>; Item No. 14010) in isolated rat peritoneal macrophages when used at a concentration of 100  $\mu$ g/ml.<sup>1</sup> Saikosaponin C (5 mg/kg per day) decreases acetic acid-induced writhing in repeated cold stressed (SART) mice.<sup>2</sup>

### References

- Ohuchi, K., Watanabe, M., Ozeki, T., *et al.* Pharmacological influence of saikosaponins on prostaglandin E<sub>2</sub> production by peritoneal macrophages. *Planta Med.* **51(3)**, 208-212 (1985).
- Kita, T., Hata, T., Itoh, E., *et al.* Analgesic and other pharmacologic actions of saiko-saponin in repeated cold stressed (SART stressed) animals. *J. Pharmacobiodyn.* **3(6)**, 269-280 (1980).

#### 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, 11/02/2022

#### 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