

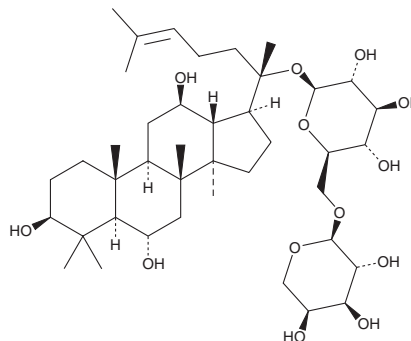
# PRODUCT INFORMATION



## Ginsenoside F<sub>3</sub>

Item No. 31409

**CAS Registry No.:** 62025-50-7  
**Formal Name:** (3 $\beta$ ,6 $\alpha$ ,12 $\beta$ )-3,6,12-trihydroxydammar-24-en-20-yl 6-O- $\alpha$ -L-arabinopyranosyl- $\beta$ -D-glucopyranoside  
**MF:** C<sub>41</sub>H<sub>70</sub>O<sub>13</sub>  
**FW:** 771.0  
**Purity:**  $\geq 95\%$   
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:**  $\geq 4$  years  
**Item Origin:** Plant/Ginseng root



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

### Laboratory Procedures

Ginsenoside F<sub>3</sub> is supplied as a crystalline solid. A stock solution may be made by dissolving the ginsenoside F<sub>3</sub> in the solvent of choice, which should be purged with an inert gas. Ginsenoside F<sub>3</sub> is soluble in methanol, ethanol, and DMSO.

### Description

Ginsenoside F<sub>3</sub> is a ginsenoside that has been found in *P. ginseng* and has immunomodulatory activity.<sup>1</sup> It increases proliferation of isolated mouse spleen cells and the production of IL-2 and IFN- $\gamma$ , as well as decreases the production of IL-4 and IL-10, when used at concentrations ranging from 0.1 to 100  $\mu$ M. Ginsenoside F<sub>3</sub> (10  $\mu$ M) also increases the DNA binding activity of NF- $\kappa$ B induced by concanavalin A (Item No. 14951) in isolated mouse spleen cells.

### Reference

1. Yu, J., Dou, D., Chen, X., *et al.* Immunoenhancing activity of protopanaxatriol-type ginsenoside-F<sub>3</sub> in murine spleen cells. *Acta Pharmacol. Sin.* **25(12)**, 1671-1676 (2004).

#### 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, 11/28/2022

#### 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](http://WWW.CAYMANCHEM.COM)