

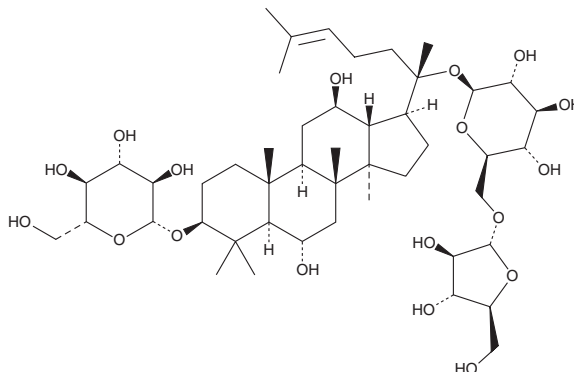
# PRODUCT INFORMATION



## Notoginsenoside Fe

Item No. 34297

**CAS Registry No.:** 88105-29-7  
**Formal Name:** (3 $\beta$ ,12 $\beta$ )-3-( $\beta$ -D-glucopyranosyloxy)-12-hydroxydammar-24-en-20-yl 6-O- $\alpha$ -L-arabinofuranosyl- $\beta$ -D-glucopyranoside  
**Synonyms:** Ginsenoside Fe, Ginsenoside Mb  
**MF:** C<sub>47</sub>H<sub>80</sub>O<sub>17</sub>  
**FW:** 917.1  
**Purity:**  $\geq 90\%$   
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:**  $\geq 4$  years  
**Item Origin:** Plant/*Panax notoginseng*



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

### Laboratory Procedures

Notoginsenoside Fe is supplied as a solid. A stock solution may be made by dissolving the notoginsenoside Fe in the solvent of choice, which should be purged with an inert gas. Notoginsenoside Fe is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of notoginsenoside Fe in these solvents is approximately 1, 5, and 10 mg/ml, respectively.

### Description

Notoginsenoside Fe is a saponin and metabolite of ginsenoside Rc (Item No. 29088) that has been found in *P. notoginseng*.<sup>1,2</sup> It is cytotoxic to L1210 murine skin lymphocytic leukemia cells (IC<sub>50</sub> = 80  $\mu$ M).<sup>1</sup>

### References

1. Bae, E.-A., Choo, M.-K., Park, E.-K., *et al.* Metabolism of ginsenoside R<sub>c</sub> by human intestinal bacteria and its related antiallergic activity. *Biol. Pharm. Bull.* **25(6)**, 743-747 (2002).
2. Liu, F., Ma, N., Xia, F.-B., *et al.* Preparative separation of minor saponins from *Panax notoginseng* leaves using biotransformation, macroporous resins, and preparative high-performance liquid chromatography. *J. Ginseng Res.* **43(1)**, 105-115 (2019).

#### 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, 11/18/2022

#### 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