

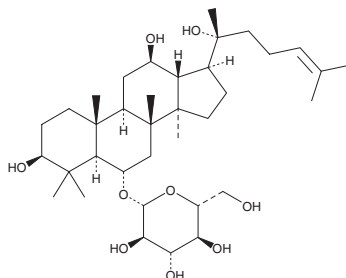
# PRODUCT INFORMATION



## 20(R)-Ginsenoside Rh<sub>1</sub>

Item No. 35493

CAS Registry No.: 80952-71-2  
Formal Name: 3 $\beta$ ,12 $\beta$ ,20R-trihydroxydammar-24-en-6 $\alpha$ -yl  $\beta$ -D-glucopyranoside  
MF: C<sub>36</sub>H<sub>62</sub>O<sub>9</sub>  
FW: 638.9  
Purity:  $\geq$ 98%  
Supplied as: A solid  
Storage: -20°C  
Stability:  $\geq$ 4 years  
Item Origin: Plant/*Ginseng Radix Et Rhizoma*



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

### Laboratory Procedures

20(R)-Ginsenoside Rh<sub>1</sub> is supplied as a solid. A stock solution may be made by dissolving the 20(R)-ginsenoside Rh<sub>1</sub> in the solvent of choice, which should be purged with an inert gas. 20(R)-Ginsenoside Rh<sub>1</sub> is soluble in the organic solvent DMSO.

### Description

20(R)-Ginsenoside Rh<sub>1</sub> is a ginsenoside that has been found in *P. ginseng* and has pro-inflammatory activity.<sup>1</sup> It induces the production of IL-12 and enhances the DNA binding activity of NF- $\kappa$ B in isolated LPS-stimulated mouse peritoneal macrophages.

### Reference

1. Lu, J.-L., Dou, D.-Q., Chen, X.-H., *et al.* Protopanaxatriol-type ginsenosides differentially modulate type 1 and type 2 cytokines production from murine splenocytes. *Planta Med.* **71**(3), 202-207 (2005).

#### 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, 11/23/2022

#### 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