

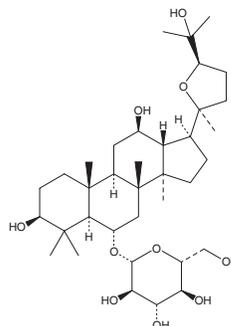
PRODUCT INFORMATION



Pseudoginsenoside RT₅

Item No. 42388

CAS Registry No.: 98474-78-3
Formal Name: (3β,6α,12β,24R)-20,24-epoxy-3,12,25-trihydroxydammaran-6-yl, β-D-glucopyranoside
Synonyms: Ginsenoside RT-5, 24(R)-Pseudoginsenoside RT₅
MF: C₃₆H₆₂O₁₀
FW: 654.9
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Pseudoginsenoside RT₅ is supplied as a solid. A stock solution may be made by dissolving the pseudoginsenoside RT₅ in the solvent of choice, which should be purged with an inert gas. Pseudoginsenoside RT₅ is soluble (≥10 mg/ml) in DMSO.

Description

Pseudoginsenoside RT₅ is an ocotillol-type ginsenoside that has been found in *P. quinquefolium*.¹

Reference

1. Li, Y., Hu, H., Li, Z., *et al.* Pharmacokinetic characterizations of ginsenoside ocotillol, RT₅ and F₁₁, the promising agents for Alzheimer's disease from American ginseng, in rats and beagle dogs. *Pharmacology* 104(1-2), 7-20 (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, 07/09/2025

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