# **PRODUCT** INFORMATION



**Zingibroside R1** 

Item No. 34130

CAS Registry No.:	80930-74-1		$\setminus$ /
Formal Name:	(3β)-17-carboxy-28-norolean-12-en-	QH	X
	3-yl 2-O-β-D-glucopyranosyl-β-D-	ОН	
	glucopyranosiduronic acid	HO	
Synonyms:	Ginsenoside Z-R <sub>1</sub> , Polysciasaponin P <sub>5</sub>		
MF:	C <sub>42</sub> H <sub>66</sub> O <sub>14</sub>	ОН ОН	
FW:	795.0		
Purity:	≥95%		H Y I V
Supplied as:	A solid		
Storage:	-20°C		
Stability:	≥4 years	O H	
Item Origin:	Plant/Panax pseudoginseng	/ \	

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

# Laboratory Procedures

Zingibroside R1 is supplied as a solid. A stock solution may be made by dissolving the zingibroside R1 in the solvent of choice, which should be purged with an inert gas. Zingibroside R1 is soluble in ethanol and DMSO.

# Description

Zingibroside R1 is a triterpene saponin and an active metabolite of ginsenoside Ro (Item No. 25137) that has been found in *P. japonicus*.<sup>1,2</sup> It is cytotoxic to B16/F10 murine melanoma cells ( $IC_{50}$  = 24.52 µg/ml) and inhibits tube formation by human umbilical vein endothelial cells (HUVECs) when used at a concentration of 40 μg/ml. Zingibroside R1 (25 mg/kg) reduces tumor growth in a B16/F10 murine melanoma model.

# References

- 1. Yoshizaki, K., Devkota, H.P., Fujino, H., et al. Saponins composition of rhizomes, taproots, and lateral roots of Satsuma-ninjin (Panax japonicus). Chem. Pharm. Bull. (Tokyo) 61(3), 344-350 (2013).
- 2. Zheng, S.-W., Xiao, S.-Y., Wang, J., et al. Inhibitory effects of ginsenoside ro on the growth of B16F10 melanoma via its metabolites. Molecules 24(16), 2985 (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.

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, 11/17/2022

# 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