

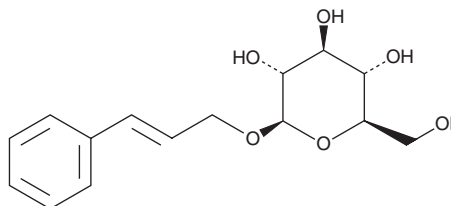
# PRODUCT INFORMATION



## Rosin

Item No. 34383

CAS Registry No.: 85026-55-7  
Formal Name: 3-phenyl-2E-propen-1-yl β-D-glucopyranoside  
MF: C<sub>15</sub>H<sub>20</sub>O<sub>6</sub>  
FW: 296.3  
Purity: ≥98%  
UV/Vis.: λ<sub>max</sub>: 252 nm  
Supplied as: A solid  
Storage: -20°C  
Stability: ≥4 years  
Item Origin: Plant/Rhodiola rosea



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

### Laboratory Procedures

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

### Description

Rosin is a cinnamic glycoside that has been found in *R. rosea* and has antidepressant and neurotropic activities.<sup>1,2</sup> *In vivo*, rosin (0.26 mg/kg) decreases immobility time in the forced swim test in rats.<sup>1</sup> It also reduces chloral hydrate-induced sleep in mice.<sup>2</sup>

### References

1. Panossian, A., Nikoyan, N., Ohanyan, N., *et al.* Comparative study of Rhodiola preparations on behavioral despair of rats. *Phytomedicine* **15(1-2)**, 84-91 (2008).
2. Sokolov, S.Y., Ivashin, V.M., Zapesochnaya, G.G., *et al.* Studies of neurotropic activity of new compounds isolated from Rhodiola rosea. *Khimiko-Farmatsevticheskii Zhurnal* **19(11)**, 1367-1371 (1985).

#### 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, 12/08/2022

#### 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