

# PRODUCT INFORMATION



## Protogracillin

Item No. 11893

CAS Registry No.: 54848-30-5

Formal Name: (3 $\beta$ ,22 $\alpha$ ,25R)-26-( $\beta$ -D-glucopyranosyloxy)-22-hydroxyfurost-5-en-3-yl O-6-deoxy- $\alpha$ -L-mannopyranosyl-(1 $\rightarrow$ 2)-O-[ $\beta$ -D-glucopyranosyl-(1 $\rightarrow$ 3)]- $\beta$ -D-glucopyranoside

MF: C<sub>51</sub>H<sub>84</sub>O<sub>23</sub>

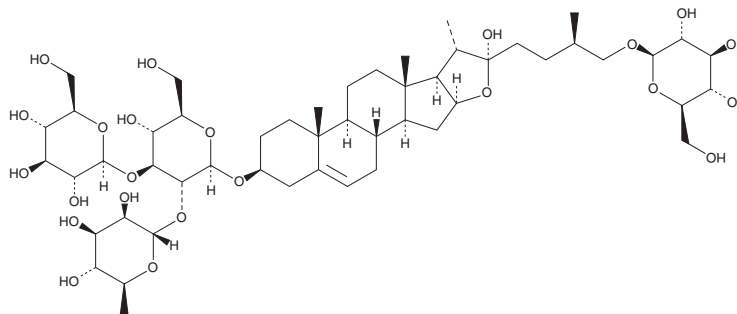
FW: 1,065.2

Purity:  $\geq$ 95%

Supplied as: A solid

Storage: -20°C

Stability:  $\geq$ 4 years



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

### Laboratory Procedures

Protogracillin is supplied as a solid. A stock solution may be made by dissolving the protogracillin in the solvent of choice, which should be purged with an inert gas. Protogracillin is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of protogracillin in these solvents is approximately 30 mg/ml.

### Description

Protogracillin is a steroid saponin originally isolated from *T. terrestris*.<sup>1</sup> It is cytotoxic to K562 cells (IC<sub>50</sub> = 3.3  $\mu$ M).<sup>2</sup>

### References

1. Tomova, M., Gyulemetova, R., Zarkova, S., et al. Steroidal saponins from *Tribulus terrestris* L. with a stimulating action on the sexual functions. *1st Int. Conf. Chem. Biotechnol. Biol. Act. Nat. Prod.* **3(1)**, 298-302 (1982).
2. Hu, K., Dong, A.-J., Yao, X.-S., et al. Antineoplastic steroidal saponins from rhizomes of *Dioscorea colleti* var. *Hypoglauca*. *Stud. Plant Sci.* **6**, 220-229 (1999).

#### 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, 10/06/2022

#### 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