

PRODUCT INFORMATION



Deslanoside Item No. 30872

CAS Registry No.: 17598-65-1
Formal Name: (3 β ,5 β ,12 β)-3-[(O- β -D-glucopyranosyl-(1 \rightarrow 4)-O-2,6-dideoxy- β -D-ribo-hexopyranosyl-(1 \rightarrow 4)-O-2,6-dideoxy- β -D-ribo-hexopyranosyl-(1 \rightarrow 4)-2,6-dideoxy- β -D-ribo-hexopyranosyl)oxy]-12,14-dihydroxy-card-20(22)-enolide

Synonym: Desacetyllanatoside C

MF: C₄₇H₇₄O₁₉

FW: 943.1

Purity: \geq 98%

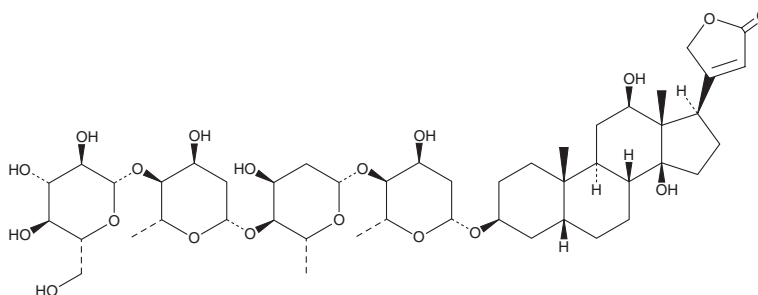
UV/Vis.: λ_{max} : 220 nm

Supplied as: A crystalline solid

Storage: -20°C

Stability: \geq 4 years

Item Origin: Plant/Unspecified sp.



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

Laboratory Procedures

Deslanoside is supplied as a crystalline solid. A stock solution may be made by dissolving the deslanoside in the solvent of choice, which should be purged with an inert gas. Deslanoside is soluble in the organic solvent DMSO at a concentration of approximately 1 mg/ml.

Deslanoside is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, deslanoside should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Deslanoside has a solubility of approximately 0.14 mg/ml in a 1:6 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Deslanoside is a cardiac glycoside that has been found in *D. lanata*.¹ It induces ventricular tachycardia in anesthetized dogs when administered intravenously at a dose of 0.18 mg/kg.² Formulations containing deslanoside have been used in the treatment of congestive heart failure and cardiac arrhythmias.

References

1. Koren, G. and Soldin, S.J. Cardiac glycosides. *Clin. Lab. Med.* **7**(3), 587-606 (1987).
2. Ghani, M.F. and Smith, J.R. The effectiveness of magnesium chloride in the treatment of ventricular tachyarrhythmias due to digitalis intoxication. *Am. Heart J.* **88**(5), 621-626 (1974).

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, 10/17/2022

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