

PRODUCT INFORMATION



Digitonin

Item No. 14952

CAS Registry No.: 11024-24-1

Formal Name: (2 α ,3 β ,5 α ,15 β ,25R)-2,15-dihydroxyspirostan-3-yl O- β -D-glucopyranosyl-(1 \rightarrow 3)-O- β -D-galactopyranosyl-(1 \rightarrow 2)-O-[β -D-xylopyranosyl-(1 \rightarrow 4)- β -D-galactopyranoside]

Synonyms:

MF: C₅₆H₉₂O₂₉

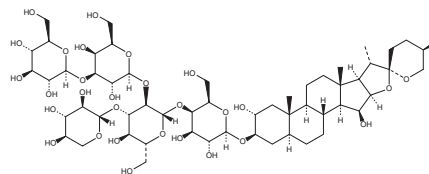
FW: 1,229.3

Supplied as: A crystalline solid

Storage: -20°C

Stability: \geq 4 years

Item Origin: Plant/*Digitalis purpurea*



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

Laboratory Procedures

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

Description

Digitonin is a glycoside obtained from *D. purpurea* that is used as a non-ionic detergent to solubilize membrane-bound proteins, to precipitate cholesterol, and to permeabilize cell membranes.¹⁻³

References

1. Nazari, M., Kurdi, M., and Heerklotz, H. Classifying surfactants with respect to their effect on lipid membrane order. *Biophys. J.* **102**(3), 498-506 (2012).
2. Gardner, J.A., Gainsborough, H., and Murray, R. Studies in the cholesterol content of normal human plasma: An improved macromethod for the estimation of cholesterol by digitonin. *Biochem. J.* **32**(1), 15-18 (1938).
3. Knight, D.E. and Scrutton, M.C. Gaining access to the cytosol: The technique and some applications of electroporeabilization. *Biochem. J.* **234**(3), 497-506 (1986).

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