

PRODUCT INFORMATION



Gitoxin

Item No. 38888

CAS Registry No.: 4562-36-1

Formal Name: (3 β ,5 β ,16 β)-3-[[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]-14,16-dihydroxycard-20(22)-enolide

Synonyms: NSC 95099

MF: C₄₁H₆₄O₁₄

FW: 780.9

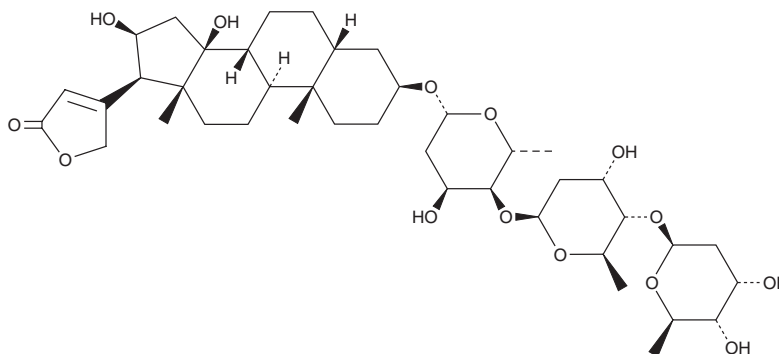
Purity: \geq 98%

Supplied as: A solid

Storage: -20°C

Stability: \geq 4 years

Item Origin: Synthetic



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

Laboratory Procedures

Gitoxin is supplied as a solid. A stock solution may be made by dissolving the gitoxin in the solvent of choice, which should be purged with an inert gas. Gitoxin is slightly soluble in chloroform and methanol.

Description

Gitoxin is a cardiac glycoside originally found in *Digitalis* that has cardiotonic and antiarrhythmic activities.¹⁻³ It inhibits Na⁺/K⁺-ATPase in an isoform-specific biphasic manner with IC₅₀ values of 0.298 and 410 μ M for the high- and low affinity human erythrocyte isoenzymes, respectively, and 1.18 and 28.5 μ M for the high- and low affinity porcine cerebral cortex isoenzymes, respectively.² Gitoxin increases contractility and rhythmicity in isolated guinea pig heart in a concentration-dependent manner.³ It is also a degradation product of the cardiac glycoside digoxin (Item No. 22266).⁴

References

1. Windaus, A. and Schwarte, G. A chloroform-insoluble glucoside from digitalis leaves, gitoxin. *Berichte der Deutschen Chemischen Gesellschaft [Abteilung] B: Abhandlungen* **58B**, 1515-1519 (1925).
2. Krstić, D., Krinulović, K., Spasojević-Tisma, V., et al. Effects of digoxin and gitoxin on the enzymatic activity and kinetic parameters of Na⁺/K⁺-ATPase. *J. Enzyme Inhib. Med. Chem.* **19(5)**, 409-415 (2004).
3. Haustein, K.O. and Hauptmann, J. Studies on cardioactive steroids. II. Structure-activity relationships in the isolated guinea-pig heart. *Pharmacology* **11(3)**, 129-138 (1974).
4. Desta, B., Kwong, E., and McErlane, K.M. Separation of digoxin, digitoxin and their potential metabolites, impurities or degradation products by high-performance liquid chromatography. *J. Chromatogr.* **240(1)**, 137-143 (1982).

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, 03/05/2024

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