

PRODUCT INFORMATION



Digitoxin

Item No. 27825

CAS Registry No.: 71-63-6
Formal Name: (3 β ,5 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-hydroxy-card-20(22)-enolide

Synonyms: Digitoxoside, Lanatoxin, NSC 7529

MF: C₄₁H₆₄O₁₃

FW: 765.0

Purity: \geq 98%

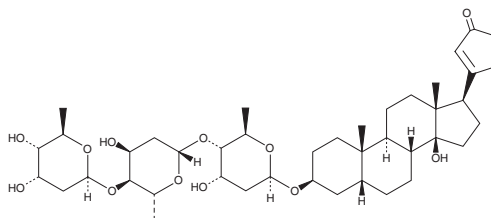
UV/Vis.: λ_{\max} : 217 nm

Supplied as: A solid

Storage: -20°C

Stability: \geq 4 years

Item Origin: Plant/*Digitalis lanata*



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

Laboratory Procedures

Digitoxin is supplied as a solid. A stock solution may be made by dissolving the digitoxin in the solvent of choice, which should be purged with an inert gas. Digitoxin is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of digitoxin in these solvents is approximately 5, 20, and 25 mg/ml, respectively.

Digitoxin is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, digitoxin should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Digitoxin has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Digitoxin is a cardiac glycoside that has been found in *Digitalis* and has diverse biological activities.¹⁻⁵ It inhibits human recombinant $\alpha_1\beta_1$, $\alpha_2\beta_2$, and $\alpha_3\beta_1$ subunit-containing Na⁺/K⁺-ATPases with K_i values of 250, 63, and 136 nM, respectively.¹ Digitoxin inhibits the human-ether-a-go-go (hERG) potassium channel, also known as K_v11.1, in HEK293 cells expressing hERG (IC₅₀ = 11.1 nM).² It enhances developed tension and contractile force in electrically stimulated isolated guinea pig left atrial muscle when used at concentrations of 0.2 and 0.4 μ M, respectively.³ Dietary administration of digitoxin (~1 mg/kg per day) attenuates congestive heart failure and reduces myocardial hypertrophy in a rat model of myocardial infarction induced by coronary artery ligation.⁴ Digitoxin is also cytotoxic to a panel of 10 human cancer cell lines, including myeloma, lymphoma, and leukemia cancer cells, with IC₅₀ values ranging from 12 to 76 nM.⁵ Formulations containing digitoxin have previously been used in the treatment of congestive heart failure and cardiac arrhythmias.

References

1. Katz, A., Lifshitz, Y., Bab-Dinitz, E., et al. *J. Biol. Chem.* **285**(25), 19582-19592 (2010).
2. Wang, L., Wible, B.A., Wan, X., et al. *J. Pharmacol. Exp. Ther.* **320**(2), 525-534 (2007).
3. Temma, K., Akeru, T., and Brody, T.M. *Eur. J. Pharm.* **76**(4), 361-370 (1981).
4. Picollo, C.T., dos Santos, A.A., Antonio, E.L., et al. *Pharmacol. Ther.* **25**(3), 265-272 (2019).
5. Johansson, S., Lindholm, P., Gullbo, J., et al. *Anticancer Drugs* **12**(5), 475-483 (2001).

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, 10/20/2022

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