

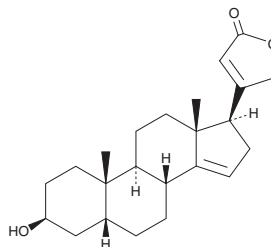
PRODUCT INFORMATION



14-Anhydrodigitoxigenin

Item No. 30439

CAS Registry No.: 4321-20-4
Formal Name: (5 β)-3 β -hydroxy-carda-
14,20(22)-dienolide
MF: C₂₃H₃₂O₃
FW: 356.5
Purity: \geq 98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 2 years



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

Laboratory Procedures

14-Anhydrodigitoxigenin is supplied as a crystalline solid. A stock solution may be made by dissolving the 14-anhydrodigitoxigenin in the solvent of choice, which should be purged with an inert gas. 14-Anhydrodigitoxigenin is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 14-anhydrodigitoxigenin in these solvents is approximately 5, 20, and 25 mg/ml, respectively.

14-Anhydrodigitoxigenin is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 14-anhydrodigitoxigenin should first be dissolved in DMF and then diluted with the aqueous buffer of choice. 14-Anhydrodigitoxigenin 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

14-Anhydrodigitoxigenin is a cardenolide and a derivative of digitoxin.¹ It reduces the activity of guinea pig heart Na⁺/K⁺-ATPase by 15% when used at a concentration of 10 μ M.²

References

1. Kamano, Y., Tozawa, M., and Pettit, G.R. Steroids and related natural products. 90. 15 β -Hydroxydigitoxigenin. *J. Org. Chem.* **40**(6), 793-796 (1975).
2. Megges, R., Weiland, J., Schön, R., *et al.* Cardenamides from cardenolides: Cardiac and anticancer activities. *Collect. Czech. Commun.* **67**(3), 336-352 (2002).

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, 03/13/2020

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