

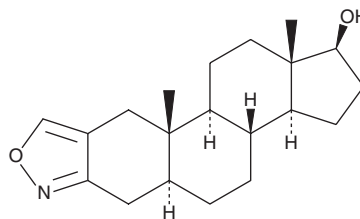
PRODUCT INFORMATION



17β-Hydroxy-androstano[3,2-c]isoxazole

Item No. 17929

CAS Registry No.: 884483-38-9
Formal Name: (5α,17β)-androstano[3,2-c]isoxazol-17-ol
MF: C₂₀H₂₉NO₂
FW: 315.5
Purity: ≥98%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid
UV/Vis.: λ_{max}: 223 nm



Laboratory Procedures

For long term storage, we suggest that 17β-hydroxy-androstano[3,2-c]isoxazole be stored as supplied at -20°C. It should be stable for at least two years.

17β-Hydroxy-androstano[3,2-c]isoxazole is supplied as a crystalline solid. A stock solution may be made by dissolving the 17β-hydroxy-androstano[3,2-c]isoxazole in the solvent of choice. 17β-Hydroxy-androstano[3,2-c]isoxazole is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of 17β-hydroxy-androstano[3,2-c]isoxazole in these solvents is approximately 15, 20, and 25 mg/ml, respectively.

17β-Hydroxy-androstano[3,2-c]isoxazole is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 17β-hydroxy-androstano[3,2-c]isoxazole should first be dissolved in DMF and then diluted with the aqueous buffer of choice. 17β-Hydroxy-androstano[3,2-c]isoxazole has a solubility of approximately 0.1 mg/ml in a 1:4 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

17β-Hydroxy-androstano[3,2-c]isoxazole is an analog of furazadrol and danazol, which are isoxazoles of testosterone (Item Nos. ISO60154 | 15645) that possess weak androgenic activity.¹ 17β-Hydroxy-androstano[3,2-c]isoxazole is a designer anabolic androgenic steroid that is classified as a Schedule III controlled substance in the United States. This compound is intended for forensic and research purposes only.

Reference

1. Parr, M.K., Gütschow, M., Daniels, J., *et al.* Identification of steroid isoxazole isomers marketed as designer supplement. *Steroids* **74**, 322-328 (2009).

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, 11/02/2015

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