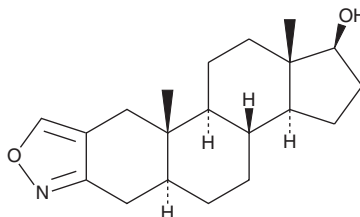


PRODUCT INFORMATION

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

Item No. 17929

CAS Registry No.: 884483-38-9
Formal Name: (5 α ,17 β)-androstano[3,2-c]isoxazol-17-ol
MF: C₂₀H₂₉NO₂
FW: 315.5
Purity: $\geq$ 98%
UV/Vis.: λ_{max} : 223 nm
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.

Description

17 β -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 17 β -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**(3), 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, 06/10/2024

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