

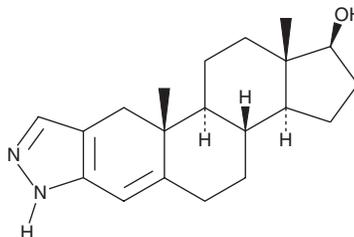
PRODUCT INFORMATION



[3,2-c]Pyrazole-androst-4-en-17 β -ol

Item No. 17930

CAS Registry No.: 929042-39-7
Formal Name: 17 β -2'H-androsta-2,4-dieno[3,2-c]pyrazol-17-ol
MF: C₂₀H₂₈N₂O
FW: 312.5
Purity: \geq 95%
UV/Vis.: λ_{max} : 212, 260 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.

Laboratory Procedures

[3,2-c]Pyrazole-androst-4-en-17 β -ol is supplied as a crystalline solid. A stock solution may be made by dissolving the [3,2-c]pyrazole-androst-4-en-17 β -ol in the solvent of choice, which should be purged with an inert gas. [3,2-c]Pyrazole-androst-4-en-17 β -ol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of [3,2-c]pyrazole-androst-4-en-17 β -ol in ethanol is approximately 10 mg/ml and approximately 30 mg/ml in DMSO and DMF.

[3,2-c]Pyrazole-androst-4-en-17 β -ol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, [3,2-c]pyrazole-androst-4-en-17 β -ol should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. [3,2-c]Pyrazole-androst-4-en-17 β -ol has a solubility of approximately 0.1 mg/ml in a 1:6 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

[3,2-c]Pyrazole-androst-4-en-17 β -ol is a steroidal heterocyclic compound that is classified as a controlled substance in the United States through the Designer Anabolic Steroid Control Act of 2014.^{1,2} This compound is intended for forensic and research purposes only.

References

1. Singh, C.B. and Singh, L.W. A facile synthesis of steroidal pyrazoles, isoxazoles and pyrimidines from testosterone. *Indian J. Chem.* **45B**, 959-966 (2006).
2. Designer Anabolic Steroid Control Act of 2014, Washington, One Hundred Thirteenth Congress of the United States of America, (2014).

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, 02/27/2023

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