

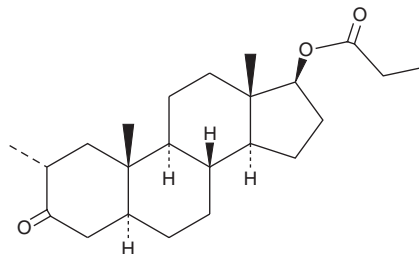
PRODUCT INFORMATION



Drostanolone Propionate

Item No. 32777

CAS Registry No.: 521-12-0
Formal Name: (2 α ,5 α ,17 β)-2-methyl-17-(1-oxopropoxy)-androstan-3-one
Synonyms: Dromostanolone Propionate, Masteron, Prometholon
MF: C₂₃H₃₆O₃
FW: 360.5
Purity: \geq 98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 5 years



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

Description

Drostanolone propionate (Item No. 32777) is an analytical reference standard categorized as an anabolic androgenic steroid.¹ Formulations containing drostanolone propionate have been used to enhance physical performance in racehorses. Drostanolone propionate is regulated as a Schedule III compound in the United States. This product is intended for research and forensic applications.

Reference

1. Kwok, K.Y., Choi, T.L., Kwok, W.H., *et al.* Detection of anabolic and androgenic steroids and/or their esters in horse hair using ultra-high performance liquid chromatography-high resolution mass spectrometry. *J. Chromatogr. A.* **1493**, 76-86 (2017).

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, 07/05/2023

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