

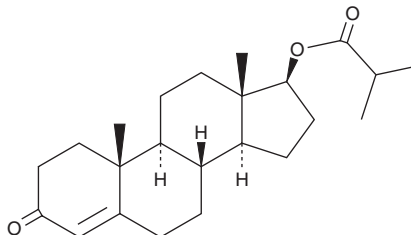
PRODUCT INFORMATION



Testosterone Isobutyrate

Item No. 41937

CAS Registry No.: 1169-49-9
Formal Name: 17 β -(2-methyl-1-oxopropoxy)-androst-4-en-3-one
Synonyms: Testosterone 2-methylpropanoate, Testosterone 17-isobutyrate
MF: C₂₃H₃₄O₃
FW: 358.5
Purity: \geq 98%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years



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

Description

Testosterone isobutyrate (Item No. 41937) is an analytical reference standard categorized as an anabolic androgenic steroid. Testosterone isobutyrate is a prodrug form of testosterone (Item Nos. ISO60154 | ISO00154 | 15645). Testosterone isobutyrate is regulated as a Schedule III compound in the United States. This product is intended for research and forensic applications.

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, 05/13/2026

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