

PRODUCT INFORMATION

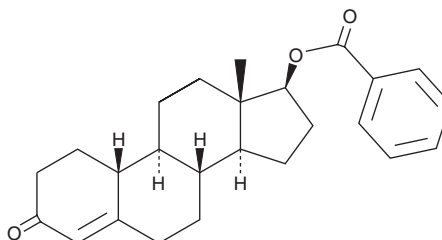


Nandrolone Benzoate

Item No. 31714

CAS Registry No.: 1474-55-1
Formal Name: (17 β)-17-(benzoyloxy)-estr-4-en-3-one
Synonyms: 4-Estren-17 β -ol-3-one benzoate,
Nandrolone 17-acetate,
19-Norandrostenolone benzoate,
19-Nortestosterone benzoate,
19-Nortestosterone-17-benzoate,
NSC 19858

MF: C₂₅H₃₀O₃
FW: 378.5
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 234 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥ 5 years



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

Description

Nandrolone benzoate (Item No. 31714) is an analytical reference standard categorized as an anabolic androgenic steroid.¹ Nandrolone benzoate is regulated as a Schedule III compound in the United States. This product is intended for research and forensic applications.

Reference

1. Lerner, L.J., Holthaus, F.J., Jr., and Thompson, C.R. A myotrophic agent and gonadotrophin inhibitor, 19-nortestosterone-17-benzoate. *Endocrinology* **64**(6), 1010-1016 (1959).

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, 06/30/2023

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