

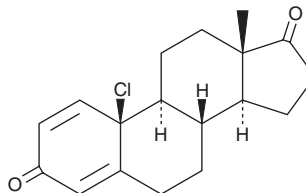
PRODUCT INFORMATION



10-Chloroestra-1,4-diene-3,17-dione

Item No. 39758

CAS Registry No.: 91413-55-7
Formal Name: 10-chloro-estra-1,4-diene-3,17-dione
Synonym: 10-chloro-1,4-Estradiene
MF: C₁₈H₂₁ClO₂
FW: 304.8
Purity: ≥95%
UV/Vis.: λ_{max}: 241 nm
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥1 year



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

Description

10-Chloroestra-1,4-diene-3,17-dione is a sterone agonist of estrogen receptor α (ER α).¹ It increases reporter gene expression in *S. cerevisiae* (EC₅₀ = 0.36 nM for the human receptor). It also is an aromatase inhibitor (IC₅₀ = 2.4 μ M).²

References

1. Nakamura, H., Shiozawa, T., Terao, Y., *et al.* By-products produced by the reaction of estrogens with hypochlorous acid and their estrogen activities. *J. Health Sci.* **52(2)**, 124-131 (2006).
2. Jójárt R., Traj P., Kovács É., *et al.* Synthesis, biological evaluation and docking studies of 13-epimeric 10-fluoro- and 10-chloroestra-1,4-dien-3-ones as potential aromatase inhibitors. *Molecules* **24(9)**, 1783 (2019).

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, 10/16/2023

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