

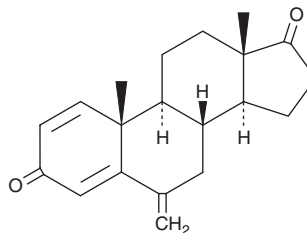
PRODUCT INFORMATION



Exemestane

Item No. 15008

CAS Registry No.: 107868-30-4
Formal Name: 6-methylene-androsta-1,4-diene-3,17-dione
Synonym: FCE 24304
MF: C₂₀H₂₄O₂
FW: 296.4
Purity: ≥95%
UV/Vis.: λ_{max}: 246 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Exemestane is supplied as a crystalline solid. A stock solution may be made by dissolving the exemestane in the solvent of choice, which should be purged with an inert gas. Exemestane is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of exemestane in ethanol is approximately 20 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Exemestane is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, exemestane should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Exemestane has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Exemestane is an irreversible inhibitor of aromatase (K_i = 26 nM).¹ It is selective for aromatase over 5α-reductase when used at a concentration of 30 μM. Subcutaneous administration of exemestane (3-100 mg/kg per day) decreases tumor proliferation in rats with mammary tumors induced by the carcinogenic polycyclic aromatic hydrocarbon 7,12-dimethylbenz(a)anthracene (DMBA; Item No. 30383).² Formulations containing exemestane have been used in the treatment of estrogen receptor-positive breast cancers in postmenopausal women.

References

1. Giudici, D., Ornati, G., Briatico, G., et al. 6-Methylenandrosta-1,4-diene-3,17-dione (FCE 24304): A new irreversible aromatase inhibitor. *J. Steroid Biochem.* **30(1-6)**, 391-394 (1988).
2. Zaccheo, T., Giudici, D., Lombardi, P., et al. A new irreversible aromatase inhibitor, 6-methylenandrosta-1,4-diene-3,17-dione (FCE 24304): Antitumor activity and endocrine effects in rats with DMBA-induced mammary tumors. *Cancer Chemother. Pharmacol.* **23(1)**, 47-50 (1989).

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, 03/06/2023

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