

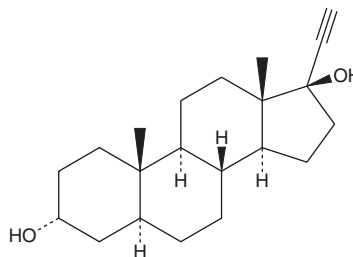
PRODUCT INFORMATION



HE-3235

Item No. 9002769

CAS Registry No.: 183387-50-0
Formal Name: (3 α ,5 α ,17 α)-pregn-20-yne-3,17-diol
Synonyms: Apoptone,
17 α -Ethynyl-5 α -androstane-3 α ,17 β -diol
MF: C₂₁H₃₂O₂
FW: 316.5
Purity: $\geq 95\%$
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥ 2 years



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

Laboratory Procedures

HE-3235 is supplied as a crystalline solid. A stock solution may be made by dissolving the HE-3235 in the solvent of choice, which should be purged with an inert gas. HE-3235 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of HE-3235 in these solvents is approximately 10 mg/ml.

HE-3235 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, HE-3235 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. HE-3235 has a solubility of approximately 0.14 mg/ml in a 1:6 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

HE-3235 is a synthetic adrenal hormone, an androstenediol analog, that acts as an androgen receptor antagonist.¹ In an androgen receptor-dependent manner, it downregulates anti-apoptotic genes, while increasing the expression of pro-apoptotic genes, thus stimulating apoptosis.¹ HE-3235 has been shown to inhibit androstenediol-dependent LNCaP cell tumor growth both *in vitro* (IC₅₀ = 6 nM) and in xenograft models *in vivo*.^{1,2}

References

1. Trauger, R., Corey, E., Bell, D., *et al.* Inhibition of androstenediol-dependent LNCaP tumour growth by 17 α -ethynyl-5 α -androstane-3 α ,17 β -diol (HE3235). *Br. J. Cancer* **100**(7), 1068-1072 (2009).
2. Koreckij, T. D., Trauger, R. J., Montgomery, R. B., *et al.* HE3235 inhibits growth of castration-resistant prostate cancer. *Neoplasia* **11**(11), 1216-1225 (2009).

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, 06/17/2024

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