

# PRODUCT INFORMATION



## HE-3235

Item No. 9002769

**CAS Registry No.:** 183387-50-0  
**Formal Name:** (3 $\alpha$ ,5 $\alpha$ ,17 $\alpha$ )-pregn-20-yne-3,17-diol  
**Synonyms:** Apoptone, 17 $\alpha$ -Ethyne-5 $\alpha$ -androstane-3 $\alpha$ ,17 $\beta$ -diol

**MF:** C<sub>21</sub>H<sub>32</sub>O<sub>2</sub>

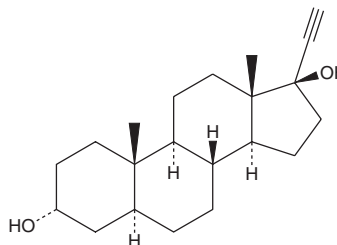
**FW:** 316.5

**Purity:**  $\geq$ 95%

**Supplied as:** A crystalline solid

**Storage:** -20°C

**Stability:** As supplied, 2 years from the QC date provided on the Certificate of Analysis, when stored properly



### 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. HE-3235 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. 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.<sup>1</sup> In an androgen receptor-dependent manner, it downregulates anti-apoptotic genes, while increasing the expression of pro-apoptotic genes, thus stimulating apoptosis.<sup>1</sup> HE-3235 has been shown to inhibit androstenediol-dependent LNCaP cell tumor growth both *in vitro* (IC<sub>50</sub> = 6 nM) and in xenograft models *in vivo*.<sup>1,2</sup>

### References

1. Trauger, R., Corey, E., Bell, D., *et al.* Inhibition of androstenediol-dependent LNCaP tumour growth by 17 $\alpha$ -ethynyl-5 $\alpha$ -androstane-3 $\alpha$ ,17 $\beta$ -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, 09/26/2016

#### 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