

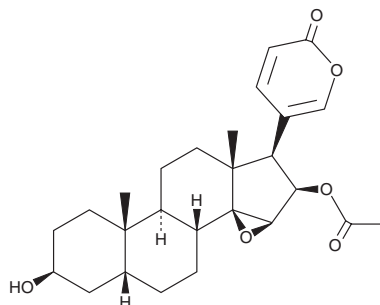
# PRODUCT INFORMATION



## Cinobufagin

Item No. 19844

**CAS Registry No.:** 470-37-1  
**Formal Name:** (5 $\beta$ )-16 $\beta$ -(acetyloxy)-14,15 $\beta$ -epoxy-3 $\beta$ -hydroxy-bufa-20,22-dienolide  
**Synonym:** NSC 90325  
**MF:** C<sub>26</sub>H<sub>34</sub>O<sub>6</sub>  
**FW:** 442.5  
**Purity:**  $\geq$ 98%  
**UV/Vis.:**  $\lambda_{\text{max}}$ : 295 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:**  $\geq$ 4 years  
**Item Origin:** Animal/*Bufois venenum*



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

### Laboratory Procedures

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

Cinobufagin is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, cinobufagin should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Cinobufagin has a solubility of approximately 0.33 mg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Cinobufagin is a cardiotonic steroid that has been found in the skin of toads of genus *Bufo* and has diverse biological activities.<sup>1-4</sup> It inhibits Na<sup>+</sup>/K<sup>+</sup>-ATPase activity in guinea pig heart ventricular muscle homogenates by 45% when used at a concentration of 0.3  $\mu$ M.<sup>1</sup> Cinobufagin is cytotoxic to HCT116 colorectal cancer cells *in vitro* with IC<sub>50</sub> values of less than 50 ng/ml at 48- and 72-hour time points and induces apoptosis in a concentration-dependent manner.<sup>2</sup> *In vivo*, cinobufagin (10 mg/kg) reduces tumor growth in an HCT116 mouse xenograft model. Cinobufagin (1  $\mu$ g/ml) inhibits LPS-induced expression of MHC class II, CD80, and CD86 and release of IL-6, IL-8, TNF- $\alpha$ , and IL-10 in human monocyte-derived dendritic cells.<sup>3</sup> It also increases expression of the antimicrobial peptides hBD2 and hBD3 in dendritic cells. Cinobufagin exhibits dose-dependent antinociceptive effects in the hot-plate, acetic acid writhing, and formalin tests in mice.<sup>4</sup>

### References

1. Tōma, S., Morishita, S., Kuronuma, K., *et al.* Metabolism and pharmacokinetics of cinobufagin. *Xenobiotica* **17(10)**, 1195-1202 (1987).
2. Lu, X.-s., Qiao, Y.-b., Li, Y., *et al.* Preclinical study of cinobufagin as a promising anti-colorectal cancer agent. *Oncotarget* **8(1)**, 988-998 (2017).
3. Xie, S., Spelmink, L., Codemo, M., *et al.* Cinobufagin modulates human innate immune responses and triggers antibacterial activity. *PLoS One* **11(8)**, e0160734 (2016).
4. Xu, L., Zhang, X., Feng, Q., *et al.* Alpha-7 nicotinic receptor-targeted cinobufagin induces antinociception and inhibits NF- $\kappa$ B signaling pathway in DRG neurons. *ACS Chem. Neurosci.* **10(1)**, 497-506 (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, 12/19/2022

#### 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