

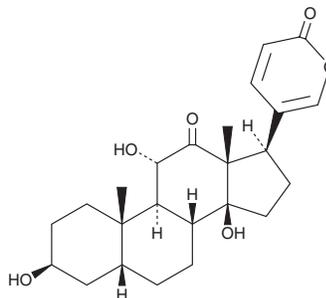
# PRODUCT INFORMATION



## Arenobufagin

Item No. 30695

**CAS Registry No.:** 464-74-4  
**Formal Name:** (3 $\beta$ ,5 $\beta$ ,11 $\alpha$ )-3,11,14-trihydroxy-12-oxo-bufa-20,22-dienolide  
**MF:** C<sub>24</sub>H<sub>32</sub>O<sub>6</sub>  
**FW:** 416.5  
**Purity:**  $\geq$ 95%  
**UV/Vis.:**  $\lambda_{\text{max}}$ : 299 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:**  $\geq$ 4 years  
**Item Origin:** Plant/*Bufo venenum*



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

### Laboratory Procedures

Arenobufagin is supplied as a crystalline solid. A stock solution may be made by dissolving the arenobufagin in the solvent of choice, which should be purged with an inert gas. Arenobufagin is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of arenobufagin in these solvents is approximately 20 and 10 mg/ml, respectively.

Arenobufagin is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, arenobufagin should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Arenobufagin has a solubility of approximately 0.11 mg/ml in a 1:8 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Arenobufagin is a cardiotoxic steroid that has been found in the skin secretions of the toad *B. gargarizans* and has anticancer activity.<sup>1-4</sup> It inhibits Na<sup>+</sup>/K<sup>+</sup>-ATPase activity in isolated guinea pig ventricular myocytes with an IC<sub>50</sub> value of 290 nM.<sup>1</sup> Arenobufagin (20-500 nM) induces cleavage of poly(ADP-ribose) polymerase (PARP), caspase-3, and caspase-9, as well as induces apoptosis in HepG2 hepatocellular carcinoma cells.<sup>2</sup> It inhibits proliferation of HepG2, Hep3B, BEL-7402, and MCF-7 cancer cells (IC<sub>50</sub>s = 101.9, 39.8, 416.3, and 143.7 nM, respectively), as well as multidrug-resistant HepG2/adm and MCF-7/adr cancer cells (IC<sub>50</sub>s = 730 and 67.5 nM, respectively). Arenobufagin (1 mg/kg once per day, i.v.) reduces tumor growth, the number and size of lung metastases, and tumor levels of the epithelial-to-mesenchymal transition (EMT) markers vimentin and  $\beta$ -catenin in a PC3 mouse xenograft model.<sup>3</sup> Arenobufagin (5 and 10  $\mu$ M/plug) inhibits VEGF-induced angiogenesis in a Matrigel™ plug assay in mice.<sup>4</sup>

### References

1. dos Santos Cruz, J. and Matsuda, H. Arenobufagin, a compound in toad venom, blocks Na<sup>+</sup>-K<sup>+</sup> pump current in cardiac myocytes *Eur. J. Pharmacol.* **239**(1-3), 223-226 (1993).
2. Zhang, D.-M., Liu, J.-S., Deng, L.-J., *et al.* Arenobufagin, a natural bufadienolide from toad venom, induces apoptosis and autophagy in human hepatocellular carcinoma cells through inhibition of PI3K/Akt/ mTOR pathway. *Carcinogenesis* **34**(6), 1331-1342 (2013).
3. Chen, L., Mai, W., Chen, M., *et al.* Arenobufagin inhibits prostate cancer epithelial-mesenchymal transition and metastasis by down-regulating  $\beta$ -catenin. *Pharmacol. Res.* **123**, 130-142 (2017).
4. Li, M.W., S., Liu, Z., Zhang, W., *et al.* Arenobufagin, a bufadienolide compound from toad venom, inhibits VEGF-mediated angiogenesis through suppression of VEGFR-2 signaling pathway. *Biochem. Pharmacol.* **83**(9), 1251-1260 (2012).

#### 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, 11/02/2022

#### 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