

# PRODUCT INFORMATION



## Aszonapyrone A

Item No. 29940

**CAS Registry No.:** 83103-08-6  
**Formal Name:** 3-[[[(1R,4aS,4bR,7S,8aR,10aS)-7-(acetyloxy)tetradecahydro-4b,8,8,10a-tetramethyl-2-methylene-1-phenanthrenyl]methyl]-4-hydroxy-6-methyl-2H-pyran-2-one

**MF:** C<sub>28</sub>H<sub>40</sub>O<sub>5</sub>

**FW:** 456.6

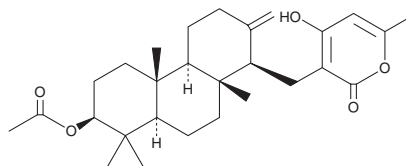
**Purity:** ≥95%

**Supplied as:** A solid

**Storage:** -20°C

**Stability:** ≥2 years

**Item Origin:** Fungi/*Aspergillus* sp.



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

### Laboratory Procedures

Aszonapyrone A is supplied as a solid. A stock solution may be made by dissolving the aszonapyrone A in the solvent of choice, which should be purged with an inert gas. Aszonapyrone A is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide.

### Description

Aszonapyrone A is a meroditerpene fungal metabolite that has been found in *Neosartorya* and has diverse biological activities.<sup>1-3</sup> It inhibits the growth of MCF-7, NCI H460, and A375-C5 cancer cells (GI<sub>50</sub>s = 13.6, 11.6, and 10.2 μM, respectively).<sup>1</sup> Aszonapyrone A is active against multidrug-resistant isolates of *S. aureus*, *E. faecalis*, and *E. faecium* (MICs = 8, 16, and 16 μg/ml, respectively) and inhibits *S. aureus* biofilm formation.<sup>2</sup> It is also active against *P. falciparum* *in vitro* (IC<sub>50</sub> = 1.34 μg/ml).<sup>3</sup>

### References

1. Eamvijarn, A., Gomes, N.M., Dethoup, T., *et al.* Bioactive meroditerpenes and indole alkaloids from the soil fungus *Neosartorya fischeri* (KUFC 6344), and the marine-derived fungi *Neosartorya lacinososa* (KUFC 7896) and *Neosartorya tsunodae* (KUFC 9213). *Tetrahedron* **69(40)**, 8583-8591 (2013).
2. Gomes, N.M., Bessa, L.J., Buttachon, S., *et al.* Antibacterial and antibiofilm activities of tryptoquivalines and meroditerpenes isolated from the marine-derived fungi *Neosartorya paulistensis*, *N. lacinososa*, *N. tsunodae*, and the soil fungi *N. fischeri* and *N. siamensis*. *Mar. Drugs* **12(2)**, 822-839 (2014).
3. Yim, T., Kanokmedhakul, K., Kanokmedhakul, S., *et al.* A new meroterpenoid tatenic acid from the fungus *Neosartorya tatenoi* KKKU-2NK23. *Nat. Prod. Res.* **28(21)**, 1847-1852 (2014).

#### 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, 03/05/2020

#### 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