

# PRODUCT INFORMATION



## Phomopsinamine

Item No. 33588

CAS Registry No.: 89085-54-1

Formal Name: (2E)-( $\beta$ S)-3-chloro- $\beta$ ,5-dihydroxy-N-methyl-L-tyrosyl-3,4-didehydro-L-valyl-3-hydroxy-L-isoleucyl-3,4-didehydro-L-prolyl-2,3-didehydro-isoleucinamide, cyclic (1<sup>5</sup>→3)-ether

Synonym: Phomopsinamine A

MF: C<sub>32</sub>H<sub>43</sub>ClN<sub>6</sub>O<sub>8</sub>

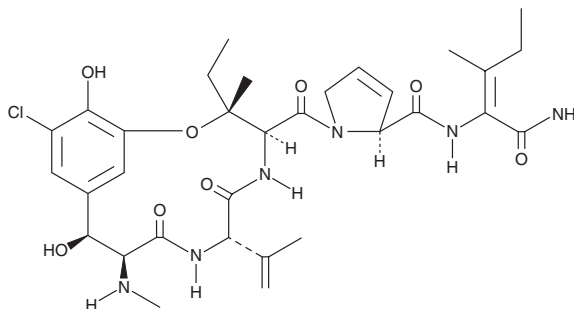
FW: 675.2

Purity:  $\geq$ 95%

Supplied as: A solid

Storage: -20°C

Stability:  $\geq$ 4 years



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

### Laboratory Procedures

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

### Description

Phomopsinamine is an inhibitor of microtubule polymerization and a derivative of phomopsin A (Item No. 19448).<sup>1</sup> It inhibits the rate and extent of polymerization of tubulin isolated from ovine brain (IC<sub>50</sub>s = 0.53 and 0.59  $\mu$ M, respectively).

### Reference

1. Lacey, E., Edgar, J.A., and Culvenor, C.C. Interaction of phomopsin A and related compounds with purified sheep brain tubulin. *Biochem. Pharmacol.* **36(13)**, 2133-2138 (1987).

#### 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, 11/15/2022

#### 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