

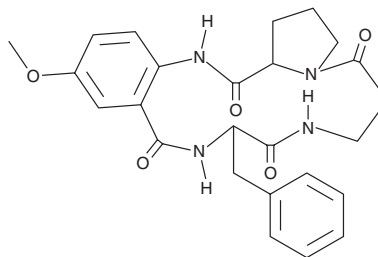
# PRODUCT INFORMATION



## Aspercolorin

Item No. 23144

**CAS Registry No.:** 29123-52-2  
**Formal Name:** cyclo( $\beta$ -alanylproyl-2-amino-5-methoxybenzoylphenylalanyl)  
**MF:** C<sub>25</sub>H<sub>28</sub>N<sub>4</sub>O<sub>5</sub>  
**FW:** 464.5  
**Purity:**  $\geq$ 95%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:**  $\geq$ 4 years  
**Item Origin:** Fungus/*Aspergillus versicolor*



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

### Laboratory Procedures

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

### Description

Aspercolorin is a fungal metabolite produced by *Aspergillus*.<sup>1</sup> It has been found in contaminated Ethiopian sorghum and finger millet.

### Reference

1. Chala, A., Taye, W., Ayalew, A., et al. Multimycotoxin analysis of sorghum (*Sorghum bicolor* L. Moench) and finger millet (*Eleusine coracana* L. Gärten) from Ethiopia. *Food Control* **45**, 29-35 (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, 02/23/2024

#### 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