

PRODUCT INFORMATION



Tentoxin

Item No. 11442

CAS Registry No.: 28540-82-1
Formal Name: cyclo[N-methyl-L-alanyl-L-leucyl-(α Z)- α,β -didehydro-N-methylphenylalanylglycyl]

MF: C₂₂H₃₀N₄O₄

FW: 414.5

Purity: \geq 99%

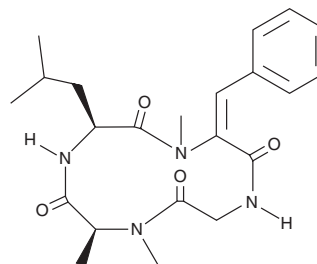
Supplied as: A powder

Storage: -20°C

Stability: \geq 4 years

Special Conditions: Protect from moisture; avoid strong, direct light

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



Laboratory Procedures

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

Description

Tentoxin is a cyclic tetrapeptide first isolated from the plant pathogen *A. alternata*. It is an herbicide that inhibits chloroplast development, resulting in chlorotic tissue.^{1,2} Tentoxin is selectively phytotoxic, as it kills a number of important weeds at concentrations that have no effects on several important crops.¹

References

1. Duke, S.O., Cantrell, C.L., Meepagala, K.M., *et al.* Natural toxins for use in pest management. *Toxins (Basel)* **2**(8), 1943-1962 (2010).
2. Duke, S.O. and Dayan, F.E. Modes of action of microbially-produced phytotoxins. *Toxins (Basel)* **3**(8), 1038-1064 (2011).

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, 12/20/2022

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