

PRODUCT INFORMATION



Territrem A

Item No. 29941

CAS Registry No.: 70407-19-1

Formal Name: (4aR,6aR,12aS,12bS)-
4a,6,6a,12,12a,12b-hexahydro-
4a,12a-dihydroxy-9-(7-methoxy-
1,3-benzodioxol-5-yl)-4,4,6a,12b-
tetramethyl-4H,11H-naphtho[2,1-b]
pyrano[3,4-e]pyran-1,11(5H)-dione

MF: $C_{28}H_{30}O_9$

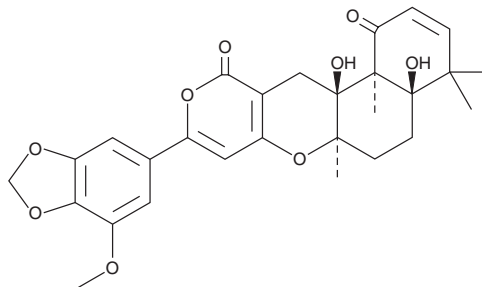
FW: 510.5

Purity: $\geq 95\%$

Supplied as: A solid

Storage: -20°C

Stability: ≥ 4 years



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

Laboratory Procedures

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

Description

Territrem A is a mycotoxin originally isolated from *A. terreus* that irreversibly inhibits acetylcholinesterase (AChE; $IC_{50} = 24$ nM for the electric eel enzyme).¹ It is toxic to mice, inducing tremors with a median tremulous dose of 0.31 mg/kg, and has an LD_{50} value of 17.6 mg/kg.²

References

1. Chen, J.-W. and Ling, K.-H. Territrems: Naturally occurring specific irreversible inhibitors of acetylcholinesterase. *J. Biomed. Sci.* **3**(1), 54-58 (1996).
2. Ling, K.H., Liou, H.-H., Yang, C.-M., et al. Isolation, chemical structure, acute toxicity, and some physicochemical properties of territrem C from *Aspergillus terreus*. *Appl. Environ. Microbiol.* **47**(1), 98-100 (1984).

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, 12/08/2022

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