PRODUCT INFORMATION



TAN 420E

Item No. 28120

CAS Registry No.: Formal Name:	91700-93-5 (15R)-18,21-didehydro-17- demethoxy-18,21-dideoxo-18,21- dihydroxy-15-methoxy-11-O- methyl-geldanamycin	
Synonym:	Dihydroherbimycin A	
MF:	$C_{30}H_{44}N_{2}O_{9}$	
FW:	576.7	
Purity:	≥95%	
Supplied as:	A lyophilisate	н Г
Storage:	-20°C	
Stability:	≥4 years	
Item Origin:	Bacterium/Streptomyces sp.	
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.		

Laboratory Procedures

TAN 420E is supplied as a lyophilisate. A stock solution may be made by dissolving the TAN 420E in the solvent of choice, which should be purged with an inert gas. TAN 420E is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide.

Description

TAN 420E is a bacterial metabolite originally isolated from Streptomyces.¹⁻³ It scavenges 2,2-diphenyl-1picrylhydrazyl (DPPH; Item No. 14805) radicals in a cell-free assay (IC₅₀ = 1.3μ M) and reduces thiobarbituric acid reactive substances (TBARS) in rat liver microsomes by 72% when used at a concentration of 100 µg/ml.³ TAN 420E is active against B. brevis, B. cereus, M. flavus, and S. aureus (MICs = 50-100 µg/ml).¹ It is cytotoxic to P388 and KB cancer cells (EC₅₀s = 0.022 and 0.3 μ g/ml, respectively).²

References

- 1. Tanida, S., Muroi, M., and Hasegawa, T. Antibiotic TAN-420, its production and use. Takeda Pharmaceutical Co. Ltd. EP0110710 A2, (1990).
- 2. Lin, L.-Z., Blaskó, G., and Cordell, G.A. H-NMR analysis of herbimycins and dihydro-herbimycins. J. Nat. Prod. 51(6), 1161-1165 (1988).
- 3. Chang, H.B. and Kim, J.H. Antioxidant properties of dihydroherbimycin A from a newly isolated Streptomyces sp. Biotechnol. Lett. 29(4), 599-603 (2007).

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, 04/18/2023

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