

Teleocidin A1

Non-phorbol ester PKC activator

Teleocidin A1 is a selective protein kinase C activator representing a new class of non-phorbol ester tumor promoters. Teleocidin A1 mimics the effect of diacylglycerol (DAG), activating PKC, but unlike DAG causes persistent activation. It enhances actin redistribution, vacuole formation and c-fos expression in PLC/PRF/5 hepatoma cells. Occurs in toxic cyanobacterial blooms.

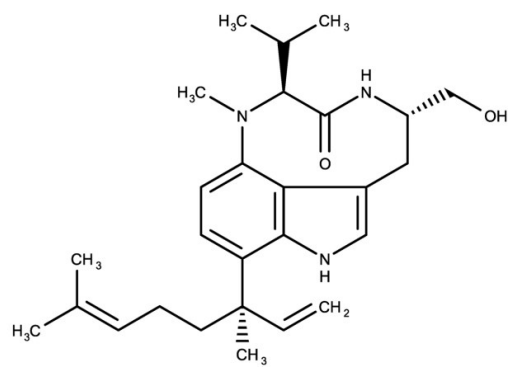
Ordering Information

[Order Online »](#)

ENZ-CHM443-0001	1mg
-----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name (10S,13S)-5-[(3R)-3,7-dimethylocta-1,6-dien-3-yl]-13-(hydroxymethyl)-9-methyl-10-propan-2-yl-3,9,12-triazatricyclo[6.6.1.0⁴,15]pentadeca-1,4,6,8(15)-tetraen-11-one

Appearance White solid/wax.

CAS 70497-14-2

Couple Type Activator

Formula C₂₇H₃₉N₃O₂

Identity Determined by NMR.

MW 437.63

Purity ≥95% (HPLC)

Solubility Soluble in DMSO (15 mg/ml).

Last modified: January 24, 2025



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)