

Harringtonine

Cephalotaxine alkaloid originally found in *Cephalotaxus*; ribosomal protein synthesis inhibitor.

Alkaloid isolated from *Cephalotaxus*. Antitumor agent.

Ordering Information

[Order Online »](#)

LKT-H0169-M005	5mg
----------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	Light sensitive.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Cephalotaxine, 4-methyl (2R)-2-hydroxy-2-(3-hydroxy-3-methylbutyl)butanedioate (ester) (9CI)
CAS	26833-85-2
Formula	$C_{28}H_{37}NO_9$
MW	531.59
Purity	≥97%
Solubility	Does not mix well with water. Soluble in DMSO to 100 mM and in ethanol to 50 mM.



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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