

Teniposide

Topoisomerase II inhibitor

Glycoside-derivative of podophyllotoxin. Has cytostatic effect. Inhibits topoisomerase II. Induces apoptosis.

Citations: 6

[View Online »](#)

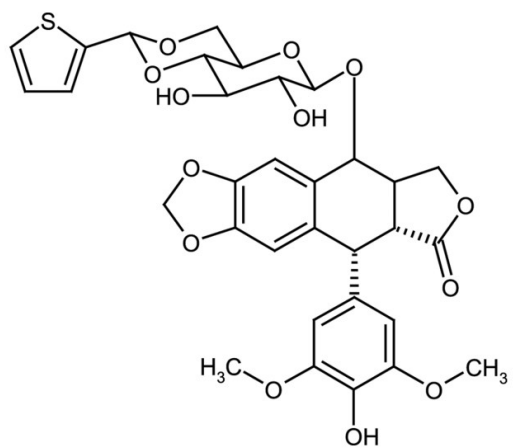
Ordering Information

[Order Online »](#)

ALX-400-053-M025	25mg
ALX-400-053-M100	100mg

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 (5R,5aR,8aR,9S)-5,8,8a,9-Tetrahydro-5-(4-hydroxy-3,5-dimethoxyphenyl)-9-[[4,6-O-[(R)-2-thienylmethylene]-β-D-glucopyranosyl]oxy]furo[3',4':6,7]naphtho[2,3-d]-1,3-dioxol-6(5aH)-one, 4'-Demethylepipodophyllotoxin-β-D-thenylidine glucoside

Appearance White to off-white powder.

CAS 29767-20-2

Couple Target Topoisomerase

Couple Type Inhibitor

Formula $C_{32}H_{32}O_{13}S$

Identity Determined by IR.

MI 14: 9145

MW 656.7

Purity ≥97% (HPLC)

Solubility Soluble in acetone or dimethyl formamide; slightly soluble in methanol; insoluble in water.

Source Semisynthetic derivative of podophyllotoxin.



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)

Last modified: May 29, 2024