

Etoposide

Topoisomerase inhibitor. p53 activator.

Antitumor reagent. Topoisomerase II inhibitor. Induces apoptosis by FasL. Activates p53 resulting in upregulated expression of TRAIL-R2 (DR5) and Bak to overcome TRAIL resistance in Bax-deficient human colon carcinoma cancer cells.

Citations: 32

[View Online »](#)

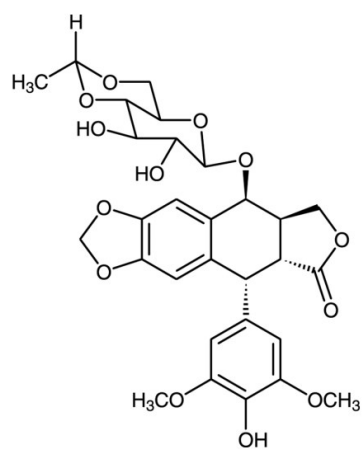
Ordering Information

[Order Online »](#)

BML-GR307-0100	100mg
BML-GR307-0500	500mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store solutions in DMSO at 4°C. For long term storage prepare aliquots and store at -20°C.
Handling	Protect from light and moisture.
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	4-Desmethylepipodophyllotoxin 9-(4,6-O-ethylidene)-β-D-glucopyranoside, VP-16-213
Appearance	White to off-white powder.
CAS	33419-42-0
Couple Target	p53, Topoisomerase
Couple Type	Activator, Inhibitor
Formula	$C_{29}H_{32}O_{13}$
MI	14: 3886
MW	588.6
Purity	≥95% (HPLC)
RTECS	KC0190000
Solubility	Soluble in DMSO, dimethyl formamide, ethyl ether or chloroform:methanol (1:1); slightly soluble in methanol or acetone; 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-
be@enzolifesciences.com](mailto:info-be@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)