

Paclitaxel

Ultra-pure; stabilizes microtubules

Antitumor agent. Stabilizes microtubules and lowers the critical concentration for tubulin polymerization. Reversibly binds to polymerized tubulin. Inhibits protein isoprenylation and induces apoptosis in PC-3 human prostate cancer cells. The checkpoint of mitotic spindle assembly, aberrant activation of cyclin-dependent kinases (CKDs) and the c-Jun N-terminal kinase/stress activated protein kinase (JNK/SAPK) are involved in Paclitaxel-induced apoptosis.

Citations: 9

[View Online »](#)

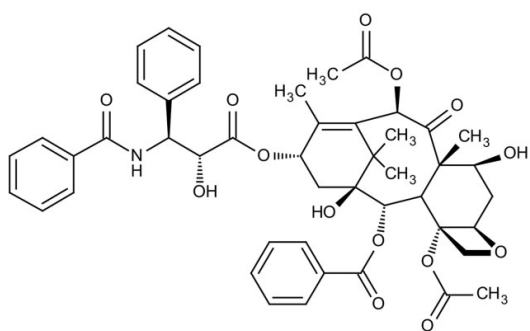
Ordering Information

[Order Online »](#)

BML-T104-0005	5mg
BML-T104-0025	25mg
BML-T104-0250	250mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store solutions at -20°C for up to 6 months.
Handling	Keep cool and dry.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	White solid.
CAS	33069-62-4
Formula	$C_{47}H_{51}NO_{14}$
MI	14: 6982
MW	853.9
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO or methanol at 10mg/ml.
Source	Isolated from the bark of the Pacific Yew tree (<i>Taxus brevifolia</i>).

Last modified: May 29, 2024



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-fr@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