# **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 »

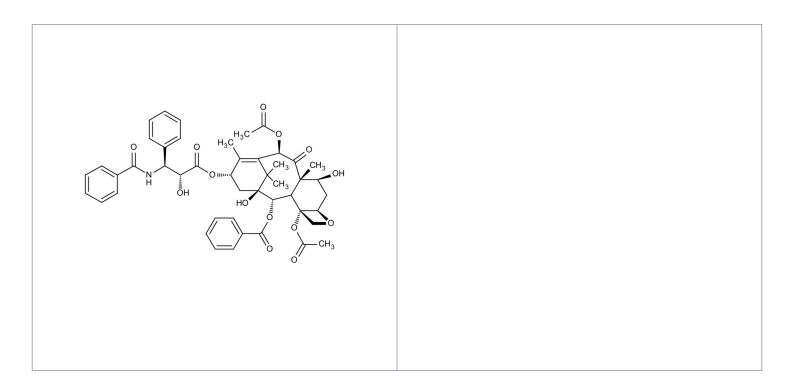
## **Ordering Information**

**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 (*Taxus* 

brevifolia).

Last modified: May 29, 2024

