

Paclitaxel (98%)

Diterpene found in *Taxus yunnanensis*;
microtubule depolymerization inhibitor.

Natural diterpenoid isolated from the stem bark of the Pacific yew tree (*Taxus brevifolia* Nutt.). It promotes the assembly of stable microtubules and inhibits the disassembly process of microtubules to tubulin by calcium ions and by low temperature. Taxol was shown to have activity against cancers, such as P388 and L1210 leukemia, B16 melanoma, WM-256 carcinosarcoma, Sarcoma 180, Lewis lung tumor, and several xenografts of human tumors including MX-1 mammary, LX-11.

Ordering Information

[Order Online »](#)

LKT-P0092-M025	25mg
----------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Taxol

CAS 33069-62-4

Couple Type Inhibitor

Formula $C_{47}H_{51}NO_{14}$

MI 13: 7052

MW 853.91

Purity ≥98%

Solubility Insoluble in water (4 mg/L). Soluble in ethanol (18 mg/mL), DMSO (50 mg/mL).



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