

Docetaxel

Ultra-pure; stabilizes microtubules

Docetaxel is taxane-type antineoplastic agent structurally similar to taxol. It reversibly binds to polymerized tubulin inhibiting depolymerization. It induces apoptosis in a large number of tumor cells and is a clinically important anticancer agent.

Citations: 12

[View Online »](#)

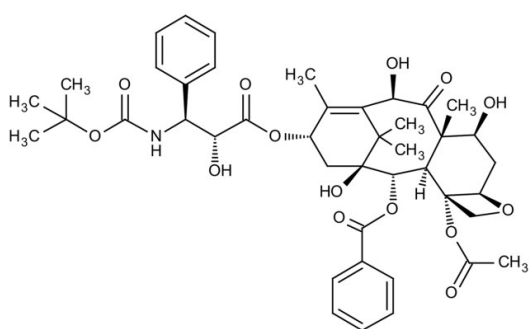
Ordering Information

[Order Online »](#)

BML-T129-0025	25mg
BML-T129-0250	250mg

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 Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Taxotere

Appearance White solid.

CAS 114977-28-5

Formula $C_{43}H_{53}NO_{14}$

MW 807.9

Purity ≥98% (HPLC)

Solubility Soluble in DMSO (25mg/ml) or 100% ethanol (25mg/ml).

Last modified: May 29, 2024



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)