

# Docetaxel

Semi-synthetic analog of taxol; microtubule depolymerization inhibitor.

A derivative of paclitaxel. Microtubule inhibitor. Induces apoptosis and inhibits angiogenesis.

## Ordering Information

[Order Online »](#)

LKT-D5709-M010	10mg
LKT-D5709-M025	25mg
LKT-D5709-M100	100mg

## 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 N-debenzoyl-N-(tert-butoxycarbonyl)-10-deacetyltaxol

CAS 114977-28-5

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

MI **13:** 3431

MW 807.89

Purity ≥98%

Solubility Soluble in DMSO, ethanol, chloroform, dichloromethane, ethyl acetate, methanol.  
Insoluble in water.



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)