

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-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

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