

Doxorubicin . hydrochloride

**Anthracycline, DNA intercalator;
topoisomerase II inhibitor.**

An anthracycline antibiotic, clinically used for chemotherapy of malignant tumors, particularly solid tumors. Acts on DNA topoisomerase II.

Ordering Information

[Order Online »](#)

LKT-D5794-M005	5mg
LKT-D5794-M010	10mg
LKT-D5794-M050	50mg

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name DXR . HCl, 14-Hydroxydaunomycin . HCl, Adriamycin . HCl

CAS 25316-40-9

Couple Target Doxorubicin

Couple Type Inhibitor

Formula $C_{27}H_{29}NO_{11} \cdot HCl$

MI 13: 3473

MW 579.99

Purity ≥98%

Solubility Soluble in water, methanol or aqueous alcohols. Soluble in DMSO to 10 mg/mL.
Insoluble in acetone or chloroform.



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