

Doxorubicin HCl

Doxorubicin HCl is an anthracycline antibiotic cancer chemotherapy drug that inhibits topoisomerase II by induction of double-stranded DNA breaking. Inhibits nucleic acid synthesis by intercalating with DNA, and induces apoptosis via p53 tumor suppressor protein accumulation. Commonly used to treat leukemias and Hodgkin's lymphoma, as well as various other cancers.

Materials Provided

Catalog Number:	92-1101
Quantity Supplied:	1 vial(s), 10 mg per vial

Description

DiscoverX control ligands have been functionally tested and validated for optimal use with all cell line targets and Assay Ready kits.

Product Information

Molecular Weight:	579.99 g/mol
Source:	Synthetic
Purity:	≥ 99%
Endotoxin Level:	N/A
Formulation:	N/A
Storage:	≤ -20°C Please avoid multiple freeze/thaw cycles.

Reagent Preparation

To avoid condensation, equilibrate the vial to ambient temperature before opening.

Stock Concentration:	25 mM
Reconstitution Volume:	0.69 mL
Reconstitution Solvent:	H ₂ O

For Research Use Only

Ordering: +1.510.979.1415 option 4 or e-mail CustomerServiceDRX@eurofins.com
Technical support: +1.510.979.1415 option 5 or e-mail DRX_SupportUS@eurofinsUS.com
General product information: www.discoverx.com