

Cisplatin

Antitumor agent

A potent platinum-based antineoplastic drug. Its mode of action involves formation of inter- and intrastrand DNA adducts which activate signaling pathways culminating in apoptosis. Platinated DNA adducts also enhance poisoning of DNA topoisomerase I. Cisplatin synergizes with a variety of anticancer agents. Tumor cells rapidly develop resistance to cisplatin and numerous agents are available to block or reverse resistance.

Citations: 51

[View Online »](#)

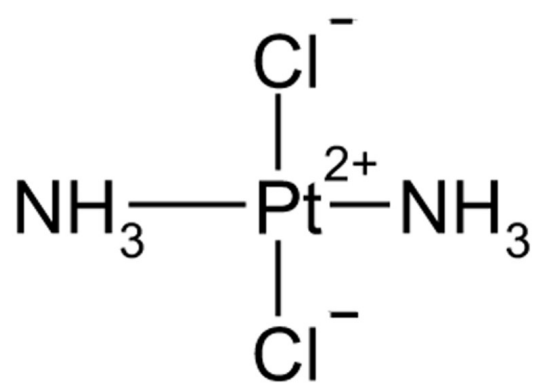
Ordering Information

[Order Online »](#)

ALX-400-040-M050	50mg
ALX-400-040-M250	250mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light.
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	cis-Platinum(II)diamine dichloride
Appearance	Off-white to orange powder.
CAS	15663-27-1
Formula	$\text{Cl}_2\text{H}_6\text{N}_2\text{Pt}$
MI	14: 2317
MW	300.0
Purity	≥98% (HPLC)
RTECS	TP2450000
Solubility	Soluble in dimethyl formamide (16mg/ml) or DMSO (10mg/ml); sparingly soluble in water; insoluble in 100% ethanol.
Technical Info / Product Notes	<p>In aqueous solution cis-trans isomerization of cisplatin occurs. Isomerization is increased at elevated temperatures. Stability of cisplatin in aqueous solutions was enhanced by increasing the sodium chloride (NaCl) concentration to 0.9% and was adversely affected in alkaline solutions such as sodium bicarbonate solutions. Further, even though soluble in DMSO we do not recommend to dissolve cisplatin in DMSO. The DMSO inserts itself into the ligand. It is recommended to prepare all solutions fresh and protect from light.</p>



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)