

Oxaliplatin

Antitumor agent

Platinum analog which has wide spectrum anticancer activity. It is found to be more active against human melanoma cell lines, and has better biochemical, pharmacological and cytotoxic properties than cisplatin and carboplatin.

Citations: 3

[View Online »](#)

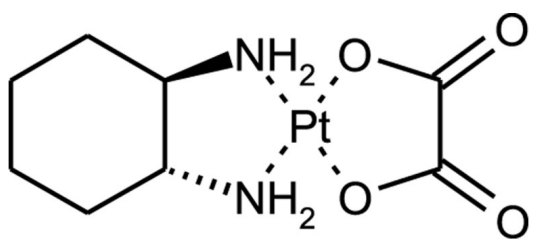
Ordering Information

[Order Online »](#)

| | |
|------------------|------|
| ALX-400-042-M005 | 5mg |
| ALX-400-042-M025 | 25mg |

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance White to off-white crystalline powder.

CAS 61825-94-3

Formula $C_8H_{14}N_2O_4Pt$

MI **14:** 6912

MW 397.3

Purity ≥98% (HPLC)

Solubility Slightly soluble in water (7.9mg/ml); very slightly soluble in methanol; almost insoluble in 100% ethanol.

Last modified: May 29, 2024



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)