

Topotecan . hydrochloride

Topoisomerase I inhibitor

Potent antitumor agent. Inhibitor of DNA-topoisomerase. Analog of (S)-(+)-camptothecin (Prod. No. ALX-350-015). Used in chemotherapy of several different cancers.

Citations: 9

[View Online »](#)

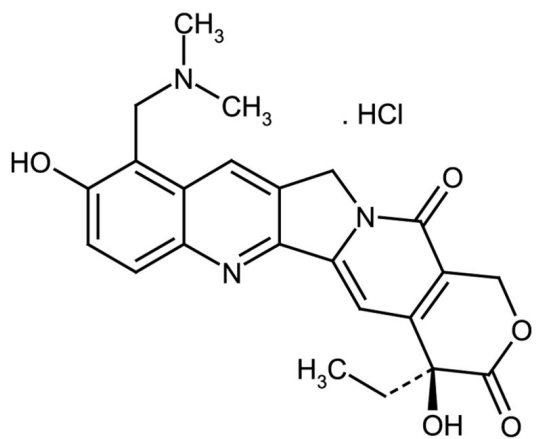
Ordering Information

[Order Online »](#)

ALX-350-133-M005	5mg
ALX-350-133-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

Alternative Name 9-[(Dimethylamino)methyl]-10-hydroxy-(20S)-camptothecin . HCl

Appearance Light yellow to greenish powder.

CAS 119413-54-6

Couple Target Topoisomerase

Couple Type Inhibitor

Formula $C_{23}H_{23}N_3O_5 \cdot HCl$

MI **14: 9548**

MW 458

Purity ≥98% (HPLC)

Solubility Soluble in water (1mg/ml) or methanol.

Source Semisynthetic.

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-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)