

Irinotecan . hydrochloride . trihydrate

Topoisomerase I inhibitor

A member of the camptothecin drug family, and an inhibitor of the nuclear enzyme topoisomerase I, which is involved in cellular DNA replication and transcription. During replication topoisomerase I mediates the relaxation of super coiled DNA, and its inhibition results in breakage of the DNA chain and likely induce apoptosis. Irinotecan is, therefore, an attractive target for anticancer drug development. Currently it is used for the treatment of small cell lung cancer and advanced colorectal cancer.

Citations: 6

[View Online »](#)

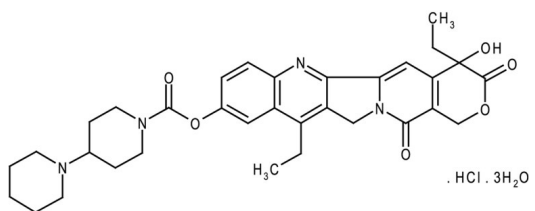
Ordering Information

[Order Online »](#)

ALX-430-139-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 +4°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Camptosar, 7-ethyl-10-hydroxycamptothecin, CPT-11

Appearance White solid.

CAS 136572-09-3

Couple Target Topoisomerase

Couple Type Inhibitor

Formula $C_{33}H_{38}N_4O_6 \cdot HCl \cdot 3H_2O$

Identity Determined by NMR.

MI 14: 5091

MW 586.7 . 36.5 . 54.0

Purity ≥97% (HPLC)

RTECS DW1061000

Solubility Slightly soluble in methanol or water.



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)