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

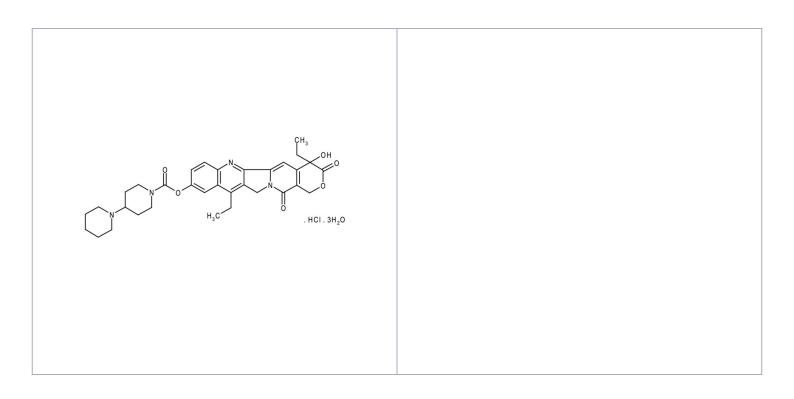
Order Online »

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 ${\rm C_{33}H_{38}N_4O_6}$. HCl . ${\rm 3H_2O}$

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.

Last modified: May 29, 2024



ENZO LIFE SCIENCES, INC.
Phone: 800.942.0430 info-usa@enzolifesciences.com

European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany Phone: +49 7621 5500 526 infode@enzolifesciences.com UK & Ireland Phone (UK customers): 0845 601 1488 Phone: +44 1392 825900 info-

uk@enzolifesciences.com