

(S)-(+)-Camptothecin

Potent antitumor agent

Potent antitumor agent. Inhibitor of DNA-topoisomerase I. Activates p53 resulting in upregulated expression of TRAIL-R2 (DR5) and Bak to overcome TRAIL resistance in Bax-deficient human colon carcinoma cells. Inhibits Tat-mediated transactivation of HIV-1. Induces apoptosis in osteosarcoma and hepatoma cells. Suppresses nitric oxide (NO) biosynthesis.

Citations: 30

[View Online »](#)

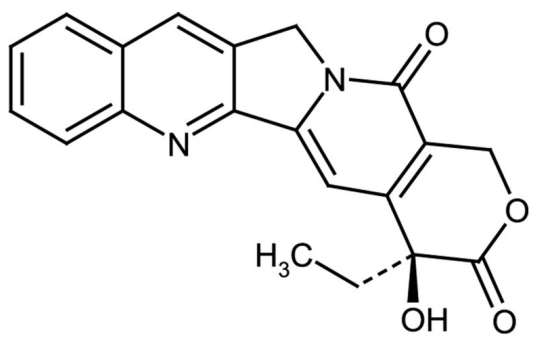
Ordering Information

[Order Online »](#)

ALX-350-015-M050	50mg
ALX-350-015-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 and moisture. After reconstitution, prepare aliquots and store at -20°C.
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	Light yellow powder or needle-shaped crystal.
CAS	7689-03-4
Couple Target	p53, Topoisomerase
Couple Type	Activator, Inhibitor
Formula	$C_{20}H_{16}N_2O_4$
Identity	Determined by IR.
MI	14: 1735
MW	348.4
Purity	≥98% (HPLC)
RTECS	UQ0492000
Solubility	Soluble in DMSO (10mg/ml), methanol (40mg/ml), 0.1N sodium hydroxide (50mg/ml) or acetic acid; insoluble in water.
Source	Isolated from <i>Camptotheca acuminata</i> seeds.



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

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info@enzolifesciences.com

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

Germany
Phone: +49 7621 5500 526
info@enzolifesciences.com

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

Last modified: May 29, 2024