

Camptothecin, -9- Amino (10mM)

9-Amino-20-(R,S)-camptothecin (also known as 9-amino campthecin) is an analogue of camptothecin. Like other CPT congeners, it shows some promising activity against a range of tumour types. This compound causes cell toxicity by stabilizing a ternary complex between the nuclear enzyme topoisomerase I (topo I) and double-stranded DNA, leading to single and double strand breaks. This reagent is provided as a control poison/inhibitor of topoisomerase I.

Ordering Information [Order Online »](#)

TOP-TG4104	0.25ml
------------	--------

Manuals, SDS & CofA [View Online »](#)

Handling & Storage

Short Term Storage -20°C

Long Term Storage -20°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 9-Amino-CPT (10mM)



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
info-eu@enzolifesciences.com

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

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

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

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