

Topoisomerase II Assay Kit

Assay kit is designed to allow quick and specific assays for eukaryotic type II DNA topoisomerases. Facilitates the purification and characterization of type II enzymes and contains all reagents necessary for routine assays of type II topoisomerases.

Ordering Information

[Order Online »](#)

TOP-TG1001-1	100 assays
--------------	------------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Short Term Storage +4°C

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Contents

Kit Contents: (representative for 100 assay kit size)

-Kinetoplast DNA (KDNA) 20ug

-Marker linear KDNA in gel loading buffer

-Marker decatenated KDNA (topo II treated)

-10X Topoisomerase II assay buffer*

*NEW: To improve stability, the 10X Topo II Assay buffer is provided as two components: Buffer A

(no ATP) and Buffer B (ATP). The complete 10X Assay buffer is made fresh each time by combining

Buffer A and B.

-5X Stop buffer/ gel loading dye

-Detailed instruction manual



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)