

Topoisomerase I Assay Kit

Assay kit contains reagents for routine detection of type I topoisomerase activity. DNA markers (included) allow unequivocal detection of the enzyme. Assay is specific for topoisomerase I; (eukaryotic type II topoisomerases require $MgCl_2$ and ATP). The standard kit is supplied with EDTA containing reaction buffer (minus Mg^{++}) in order to prevent nucleolytic nicking of the supercoiled substrate (although we note that some eukaryotic type I activities are stimulated by Mg^{++}). Assay will work even with crude extracts.

Ordering Information

[Order Online »](#)

TOP-TG1015-2	250 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



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