

DNA Gyrase Assay Kit

Kit is designed to allow quick and specific detection of DNA gyrase. This kit facilitates the purification and characterization of type II enzymes (DNA gyrase) and contains all reagents necessary for routine assays of type II topoisomerases that either have or do not have the ability to supercoil. The gyrase assay is based upon a two step process: 1) decatenation of kinetoplast DNA (KDNA); 2) supercoiling of the resulting decatenated monomer KDNA species.

Ordering Information

[Order Online »](#)

TOP-TG1003	100 assays
------------	------------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling Avoid freeze/thaw cycles.

Short Term Storage -20°C

Long Term Storage -80°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