

E. coli DNA Gyrase and Relaxed DNA Assay Kit (500 assays)

The E. coli DNA Gyrase and Relaxed DNA Assay Kit is designed to provide the investigator quick and specific assays for DNA gyrase. These kits facilitate the purification and characterization of gyrase by providing controlled reagents that we certify to work. These kits are also appropriate for the researcher who is interested in large scale screening for novel anti-gyrase active agents. We matched the amount of DNA gyrase enzyme with the amount of relaxed pHOT-1 DNA to get you going and keep you going on the hunt for new actives and leads.

Ordering Information

[Order Online »](#)

TOP-TG2000G-3KIT	500 assays
------------------	------------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	Note that ATP containing buffers can and will deactivate with repetitive freeze thawing cycles. Given the excellent stability and purity of our gyrase, we recommend that you first check buffer deterioration before assuming loss of enzyme activity.
Short Term Storage	-20°C
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Contents	Relaxed pHOT-1 DNA High purity E. coli DNA gyrase enzyme Assay buffer Dilution buffer for enzyme Detailed instruction manual
-----------------	--

Technical Info / Product Notes NOTE: Do not confuse this kit with the DNA Gyrase Assay kit (TOP-TG1003) which is designed for researchers who wish to assay for gyrase activity in extracts prepared by the investigator. The DNA Gyrase Assay Kit does not include enzyme.



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