

# E. coli Bacterial DNA Gyrase (100 U)

Highly purified *E. coli* gyrase offered as a holoenzyme (contains both A and B subunits). Excellent reagent for supercoiling plasmids *in vitro* or for novel drug screens.

A 5X assay buffer and dilution buffer is included with the enzyme.

## Ordering Information

[Order Online »](#)

TOP-TG2000G-1	100U
---------------	------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Handling	After opening, prepare aliquots and store at -80°C. Avoid freeze/thaw cycles.
Short Term Storage	-80°C
Long Term Storage	-80°C
Shipping	Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** gyr (Subunits A, B)



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)