

E. coli Bacterial DNA Gyrase (2'000 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-7 | 2000U |
|---------------|-------|

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

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