

Topoisomerase IV (E. coli) (recombinant) (His)

Topo IV is a type II topoisomerase that is encoded by two genes *parC* and *parE* in *E. coli*. Topo IV is essential for decatenation and chromosomal segregation in *E. coli*; thus, topo IV tends to prefer intermolecular strand passing (catenating and decatenating reactions) over relaxation reactions. Topo IV is also the target for quinolone-based antibacterial agents which act by subverting the enzyme into a DNA damaging agent.

Ordering Information

[Order Online »](#)

TOP-TG2000EC-3	500U
----------------	------

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

Activity	Very active in decatenating KDNA (Prod. No. TOP-2013).
Alternative Name	parC



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