

Topoisomerase I (human) (recombinant) (His)

Catalytically active wild type human topoisomerase I purified to homogeneity.

Ordering Information

[Order Online »](#)

TOP-TG2005H-RC1	250U
-----------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	Stable for 6 months when stored at -20°C.
Short Term Storage	-20°C
Long Term Storage	-20°C
Shipping	Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name TOP1

Specific Activity One unit of topoisomerase I is defined as the amount of enzyme which can relax 0.25µg of supercoiled plasmid DNA in 30 min. at 37°C.

Technical Info / Product Notes 0.25ml Topo I dilution buffer is provided as well as the assay buffer ([TOP-4010](#)).



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