

Topoisomerase I (Y723F mutant) (human) (recombinant) (His)

Catalytically inactive mutant form of human topoisomerase I purified to homogeneity.

Ordering Information

[Order Online »](#)

TOP-TG2005H-Y723F	10µg
-------------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Short Term Storage -80°C

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Activity Enzyme cannot relax supercoiled DNA, but still binds to DNA.

Technical Info / Product Notes Topo I Reaction Buffer (10X) is supplied. Supercoiled plasmid DNA is included.



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