

Topoisomerase II α (p170) (human) (recombinant)

Contains purified DNA Topoisomerase II (p170) overexpressed and purified to homogeneity to a single band on SDS-PAGE. Included with the enzyme are two buffers (10X Incomplete Buffer A and 10X ATP Buffer B) ([TOP-4040](#)) that are used to make a final working stock of 10X topo II assay buffer, as described below. The dilution buffer ([TOP-4090](#)) is sold separately.

Ordering Information

[Order Online »](#)

TOP-TG2000H-1	250U
---------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	Stable for 1-2 days at +4°C. Buffers A and B should be kept at -20°C. The enzyme needs to be kept at -80°C.
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 TOP2A

Specific Activity One unit of topoisomerase II is defined as the amount of enzyme which can decatenate 0.2µg of KDNA in 30 min. at 37°C.

Technical Info / Product Notes

Two buffers that are used to prepare a final 10X working stock of assay buffer are provided:

Note that the final 10X Complete Buffer contains both A and B and should be made fresh daily!

1. 10X Incomplete Buffer A contains the following: 0.5M TRIS-HCl (pH 8.0), 1.2M KCl, 0.1M MgCl₂, 5mM dithiothreitol). [0.5mL provided]
2. 10X ATP Buffer B contains 50mM ATP in water [50µl provided].

These two buffers must be made fresh and mixed to make the final 10X Complete Topo II Assay Buffer (A+B) as follows. Simply add 0.1 volume of Buffer B to Buffer A. Prepare only as much of the Complete Topo II Assay Buffer (A+B) as needed for each day or set of experiments. For example: If you are performing 10 assays (20 µl/assay each) in a single day, you will need roughly 20 µl of the 10X Complete Buffer (20 µl/reaction x 10 reactions). You should prepare 25 µl of 10X Complete Buffer (A+B) by mixing 25 µl of Incomplete Buffer A (-ATP) and 2.5 µl of ATP Buffer B. Note that this buffer should be kept on ice and used the same day, then discarded.



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