

# Topoisomerase II Immunoprecipitation Kit

The immunoprecipitation (IP) kit contains the key reagents and protocol for immunoprecipitating p170 human topo II $\alpha$  ([Prod. No. TOP-2000H](#)). The method should work well with other cells besides human cells. For example, mouse, rat and monkey cells will cross react with topo II antibody ([Prod. No. TOP-2011-1](#)). The antibody is specific for p170 (topo II $\alpha$ ) in human (HeLa) cells and although it is presumed that it will behave similarly in other species with cross reactive topo II isozymes, this should be carefully evaluated.

## Ordering Information

[Order Online »](#)

TOP-TG1035	1Kit
------------	------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Short Term Storage -20°C

Long Term Storage -20°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

Quantity >50 analyses



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)