

# Topoisomerase II Immunofluorescence Kit

Assay kit contains all the reagents for cytolocalization of human, p170 topo II $\alpha$  ([Prod. No. TOP-2000H](#)) using immunofluorescence. The system is optimized for use with FITC or Rhodamine labeled secondary antibody with adherent cells. The method may work with other cells besides human. 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 $\alpha$ ).

## Ordering Information

[Order Online »](#)

TOP-TG1030	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)