

# Topoisomerase IV Assay Buffer (10X)

Topo IV Assay Buffer is supplied as a 10X Stock solution in two parts that must be mixed prior to use:

1. Incomplete Buffer A contains the following: 200mM HEPES-KOH [pH7.6], 500mM potassium glutamate, 50mM magnesium acetate, 50mM dithiothreitol, 250µg BSA/ml.
2. ATP Buffer B contains 100mM ATP in water.

These two buffers must be mixed to prepare a final 10X complete Topo IV Assay Buffer (A+B) as follows: add 0.1 volume of Buffer B to 1.0 volume Buffer A. Prepare only as much of the complete Topo IV Assay Buffer (A+B) as needed for each day or set of experiments. Note that this buffer should be kept on ice and used the same day, then discarded.

## Ordering Information

[Order Online »](#)

TOP-TG4032	1ml
------------	-----

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

**Use/Stability** The two-part buffer is shipped on blue ice (or on dry ice if included with enzyme) and should be stored at -20°C. The complete Topo IV Assay Buffer (A+B) is stable for 24 hours at 4°C. The complete buffer should be made fresh for each experiment.

**Short Term Storage** -20°C

**Long Term Storage** -20°C

**Shipping** Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details



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)