

# pRYG DNA (supercoiled)

pUC19 derivative in which the 206 bp SphI-NarI fragment of pUC19 was replaced with a 245 bp SphI-AccI fragment upstream of the human adult  $\beta$ -globin gene promoter. The 245 bp fragment contains a 54 bp alternating repeat of purine/pyrimidine sequence that constitutes a strong eukaryotic topoisomerase II cleavage site.

## Ordering Information

[Order Online »](#)

TOP-TG2023-2	20 $\mu$ g
--------------	------------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Short Term Storage +4°C

Long Term Storage +4°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)