

# Human Genomic DNA, Male

**A normal pool of male genomic DNA that can serve as a standard in molecular biology applications**

Human Genomic DNA, Male, is comprised of a pool of  $\geq 8$  normal donors. It is intact DNA extracted from freshly harvested whole blood of healthy males. DNA is treated with DNase free-RNase to remove residual contaminant RNA.

Human genomic DNA can be used for a variety of molecular applications including PCR, southern blotting, and DNA library construction. It can also be used as normal reference DNA for comparative genomic hybridizations (CGH).

Citations: 4

[View Online »](#)

## Ordering Information

[Order Online »](#)

ENZ-GEN117-0100	100µg
-----------------	-------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

**Handling** Avoid freeze/thaw cycles.

**Short Term Storage** +4°C

**Long Term Storage** -20°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** gDNA

**Application Notes** Genomic DNA is suitable for PCR, Southern blot analysis, DNA library construction, and used as reference DNA in CGH and other hybridizations.

**Concentration** 150 – 250 µg/mL in 10 mM Tris-HCl, pH 8.0, containing 1 mM EDTA

**Formulation** Liquid.

**Quality Control**

- A260/280 is between 1.8 and 2.0
- Sex specificity determined by PCR
- The integrity of each DNA sample is verified by agarose gel electrophoresis before and after incubation at 37°C for 16 hours
- Genomic DNA is successfully digested by *AluI* and *RsaI* and migration on agarose gel was as expected
- Molecular weight is greater than 20 kb

**Last modified: May 29, 2024**



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-be@enzolifesciences.com)

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