## Human Genomic DNA, Female

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

Human Genomic DNA, Female, is comprised of a pool of  $\geq$  8 normal donors. It is intact DNA extracted from freshly harvested whole blood of healthy females. 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: 1

View Online »

**Ordering Information** 

**Order Online** »

ENZ-GEN118-0100

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 Alul and Rsal 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

European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com