

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 ≥ 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	100 μ g
-----------------	-------------

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	<ul style="list-style-type: none">• 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 <i>AluI</i> and <i>RsaI</i> 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
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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