

---

## Human Genomic DNA

(Catalog No. X11000)

---

### Description

Human Genomic DNA is highly purified and isolated from human placenta. The average length of the DNA is greater than 50Kb and is suitable for Southern blotting, genomic library construction and as a control template.

### Concentration

The Human Genomic DNA is supplied at a concentration of 200ng/μl.

### Size

0.2 ml

### Storage Buffer

10 mM Tris-HCl pH 8.0, 1 mM EDTA

### Storage

Human Genomic DNA should be stored at -20°C for up to 12 months. Repeated freeze/thaw cycles should be avoided.

### Extract Origin

Human placenta tissue

### Applications

Southern blotting; Genomic library construction; Control template

### Ordering Information:

Products	Size	Cat. No.
Human Genomic DNA	0.2 ml	X11000 -1

*This product is for research purposes only. Not intended for use in diagnostic procedures.*