
Human Genomic DNA

(Catalog No. X11000)

Description

Human Genomic DNA is highly purified and isolated from human placenta. This DNA 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.