

**Human Genomic DNA**

Positive control template for PCR

Cat. No.	Amount
PCR-261	20 µg

For general laboratory use.**Shipping:** shipped on gel packs**Storage Conditions:** store at -20 °C**Shelf Life:** 12 months**Form:** liquid**Description:**

Human genomic DNA is recommended for genomic analysis i.g. as template in positive control PCRs. The purified DNA comes from multiple donors. DNA fragments are longer than 50 kb.

Quality control:

Quality and concentration is confirmed by UV/Vis spectrometry.