

---

## Fully CpG Methylated Human Genomic DNA

(Catalog No. X10000)

---

### Description

Methylated human genomic DNA

### Source

Genomic DNA from female human cells was isolated by a standard genomic purification method. The DNA was enzymatically treated to methylate every CpG site.

### Concentration

10 µg in 0.1 ml

### Storage

-20°C, avoid multiple freeze/thaw cycles. Stable for 1 year from the date of shipment.

### Quality Assurance

Purified free of contaminating protein and RNA. A260/280 ratio: 1.8

### Application

Used as the positive control for methylation-specific PCR, bisulfite sequencing and various DNA methylation analyses.

### Research use

Research use only.

### Handling

For maximum recovery of the products, centrifuge the vial prior to opening the cap.

### Ordering Information:

Products	Size	Cat. No.
Fully CpG Methylated Human Genomic DNA	10 µg/100 µl	X10000 -1

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