
Human Brain Genomic DNA

(Catalog No. X11001-1)

Description

This human brain genomic DNA is highly purified and carefully isolated from human adult normal brain tissues. Evidence shows that DNA isolated from human brain tissues contains high amounts of 5-hydroxymethylcytosine (5-hmC), a modified cytosine which plays an important role in epigenetic regulation of gene function. This DNA is suitable for genomic library construction and as a control DNA for 5-hmC detection. A260/280 is between 1.8 and 2.0, and A260/230 is about 2.0 as tested by spectrophotometer.

Concentration

250 ng/μl

Size

10 μg in 40 μl

Storage Buffer

10 mM Tris-HCl pH 8.0, 1 mM EDTA

Quality Assurance

Purified free of contaminating protein and RNA. A260/280 ratio: 1.8, and A260/230 ratio: ~2.0

Storage

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

Extract Origin

Human adult normal brain tissue

Research Use

For research use only.

Applications

Genomic library construction; Control for 5-hmC detection

Products

Human Brain Genomic DNA

Size

10 μg

Cat. No.

X11001 -1

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