

## Jurkat genomic DNA

**Catalog No:** 55007

**Lot No:** 05421005

**Format:** 10 µg

**Quality Control:** Contents: 10 µg Jurkat genomic DNA (50 ng/µl)

A260/280 ratio: 1.83

Jurkat cells were grown to confluency in RPMI media containing 10% FBS. Genomic DNA was isolated and purified. 300 ng of Jurkat genomic DNA is quality control tested by restriction digest with Mse I and run on a 2% agarose gel in TAE buffer. Undigested Jurkat DNA should produce a large molecular weight band, while digested Jurkat genomic DNA should produce a smear of DNA fragments, as shown in Figure 1.

Jurkat genomic DNA is provided undigested. To use Jurkat genomic DNA as a control in the MethylCollector™ Ultra Kit (Cat. No. 55005), it must first be digested with a restriction enzyme of choice like Mse I, Bfa I, Tas I or Csp6 I.

**Storage and Guarantee:** Store at -20°C. This product is guaranteed stable for 6 months from date of receipt when stored properly.

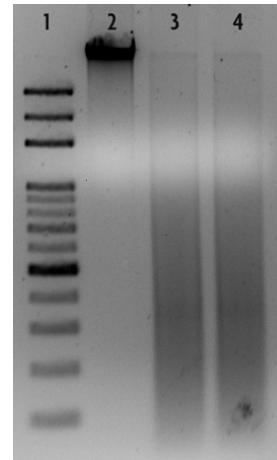


Figure 1.

1. 100 bp DNA ladder
2. 300 ng undigested Jurkat genomic DNA
3. 300 ng Mse I digested Jurkat genomic DNA
4. 300 ng Mse I digested Fully Methylated Jurkat DNA (Catalog No. 55003)