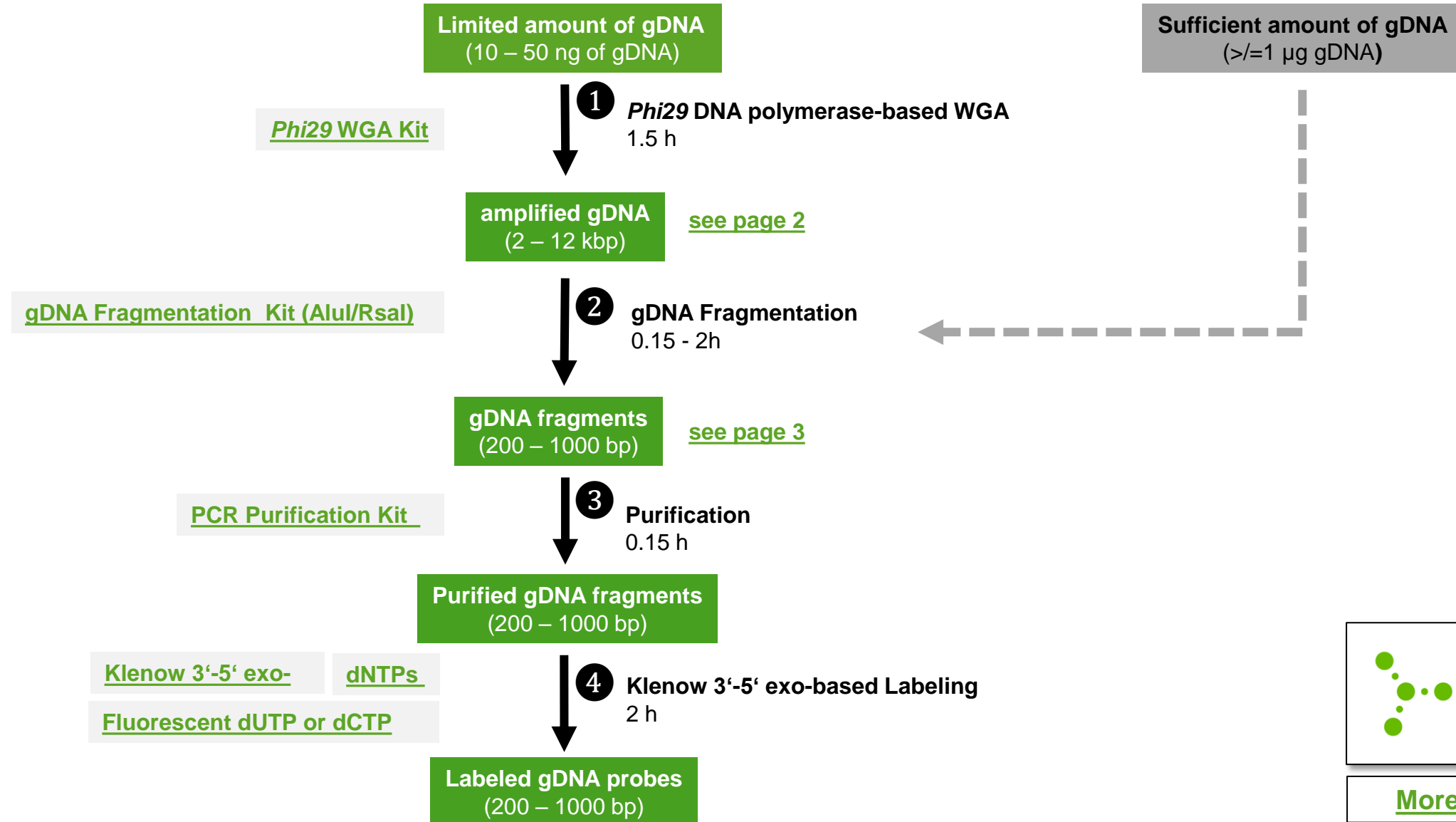
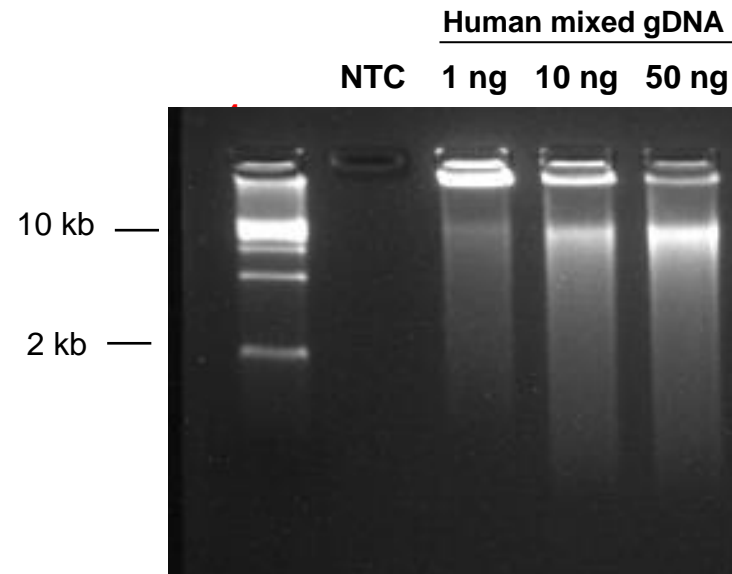


Labeled DNA probes from limited amount of genomic DNA (gDNA) are efficiently prepared via *Phi29*-based whole genome amplification (WGA)



Phi29 WGA Kit produces an average product length > 10 kb (major band at 10 – 12 kb, range between 2 kb and approximately 100 kb)

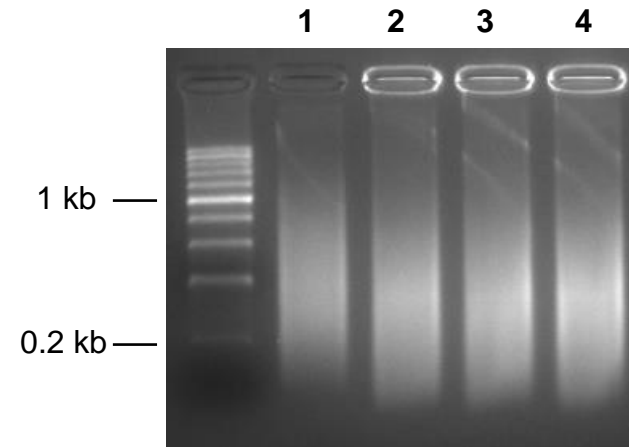


- WGA reaction: 1.5 h, 30°C
- 2 µl of 20 µl reaction analyzed via agarose gel electrophoresis (1% agarose gel, 1x TAE, EvaGreen staining)
- NTC: Non Template Control

gDNA is fragmented via AluI / RsaI digestion

Recommended starting point for individual AluI / RsaI fragmentation assay set-up:

50 U each AluI / RsaI in 10x Universal Buffer plus 0.5 µg BSA & 1 mM DTT, 2h at 37°C



1 µg of purified human mixed gDNA (lane 1) or 17 µl of WGA reaction mix (derived from Phi29 WGA Kit) (lane 2-4) have been incubated under different fragmentation conditions and analyzed via agarose gel electrophoresis (10 µl of 50 µl reaction mix, 2.4% agarose gel, 1x TAE, EvaGreen staining)

- 1: 50 U each AluI / RsaI, 2h at 37°C**
- 2: 50 U each AluI / RsaI, 2h at 37°C**
- 3: 50 U each AluI / RsaI, 0.5 µg BSA, 2h at 37°C**
- 4: 50 U each AluI / RsaI, 0.5 µg BSA, 1 mM DTT, 2h at 37°C**