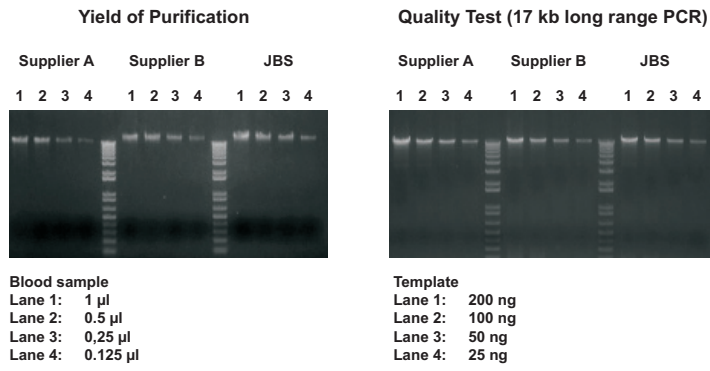
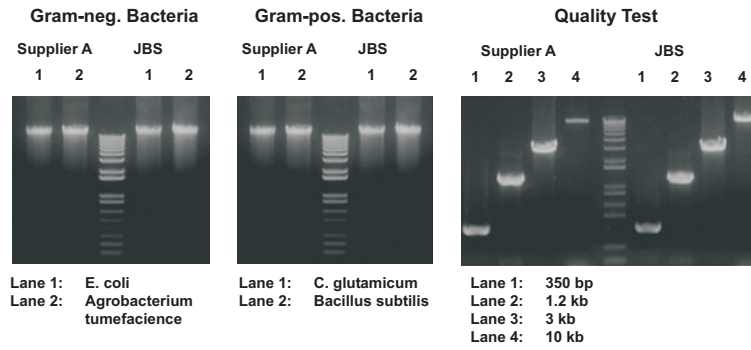


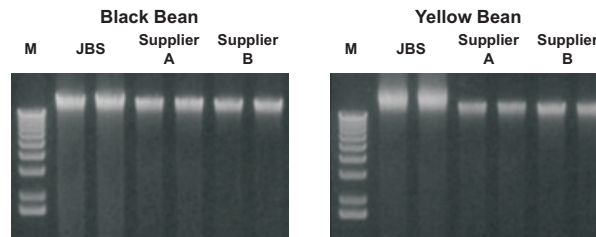
## Genomic DNA purification from whole blood



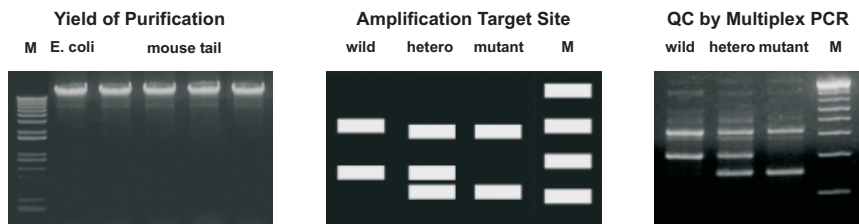
## Genomic DNA purification from bacteria



## Genomic DNA purification from plant tissue



## Genomic DNA purification from animal tissue and fungi



## Genomic DNA purification from yeast

