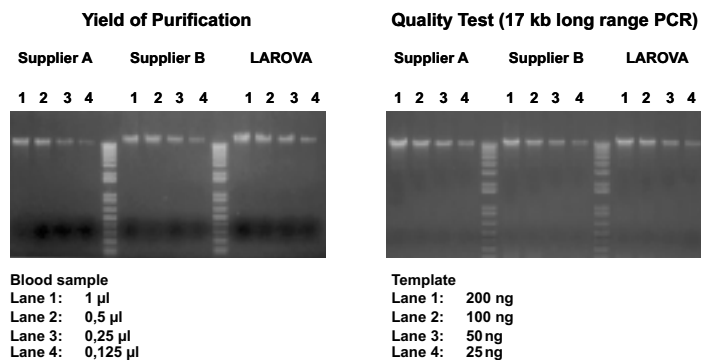
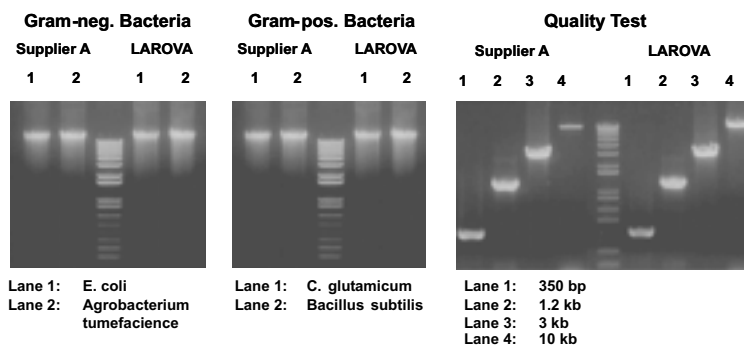


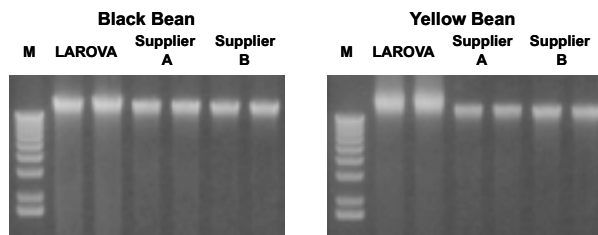
Genomic DNA purification from whole blood



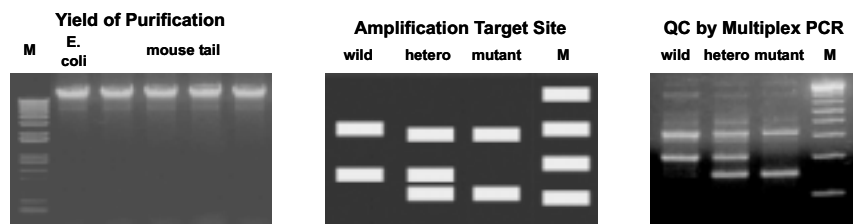
Genomic DNA purification from bacteria



Genomic DNA purification from plant tissue



Genomic DNA purification from animal tissue and fungi



Genomic DNA purification from yeast

