

Gene (DNA) Microarrays

- Using thermal inkjet technology (10,000 fires/s at 10 m/s velocity) it is feasible to immobilize thousands of genes or oligonucleotides on a small surface such as a nylon membrane, glass slide or silicon chip
- Levels of gene expression (i.e. mRNA levels) reflect colour differences on chips

