

Transcriptomics with Gene (DNA) Microarrays

- DNA oligonucleotides can be synthesized and immobilized on surfaces to permit the capture of mRNA from tissue and cell extracts
- Levels of gene expression (i.e. mRNA levels) reflect colour differences on chips

