

# A Large Genome Centre Core Pipeline Refresh

Scott Thurston\*, Peter Ellis, Lesley Shirley, Carol Scott, James Glover, Benjamin Farr, Jamieson Lovell, Richard Rance, Tony Cox

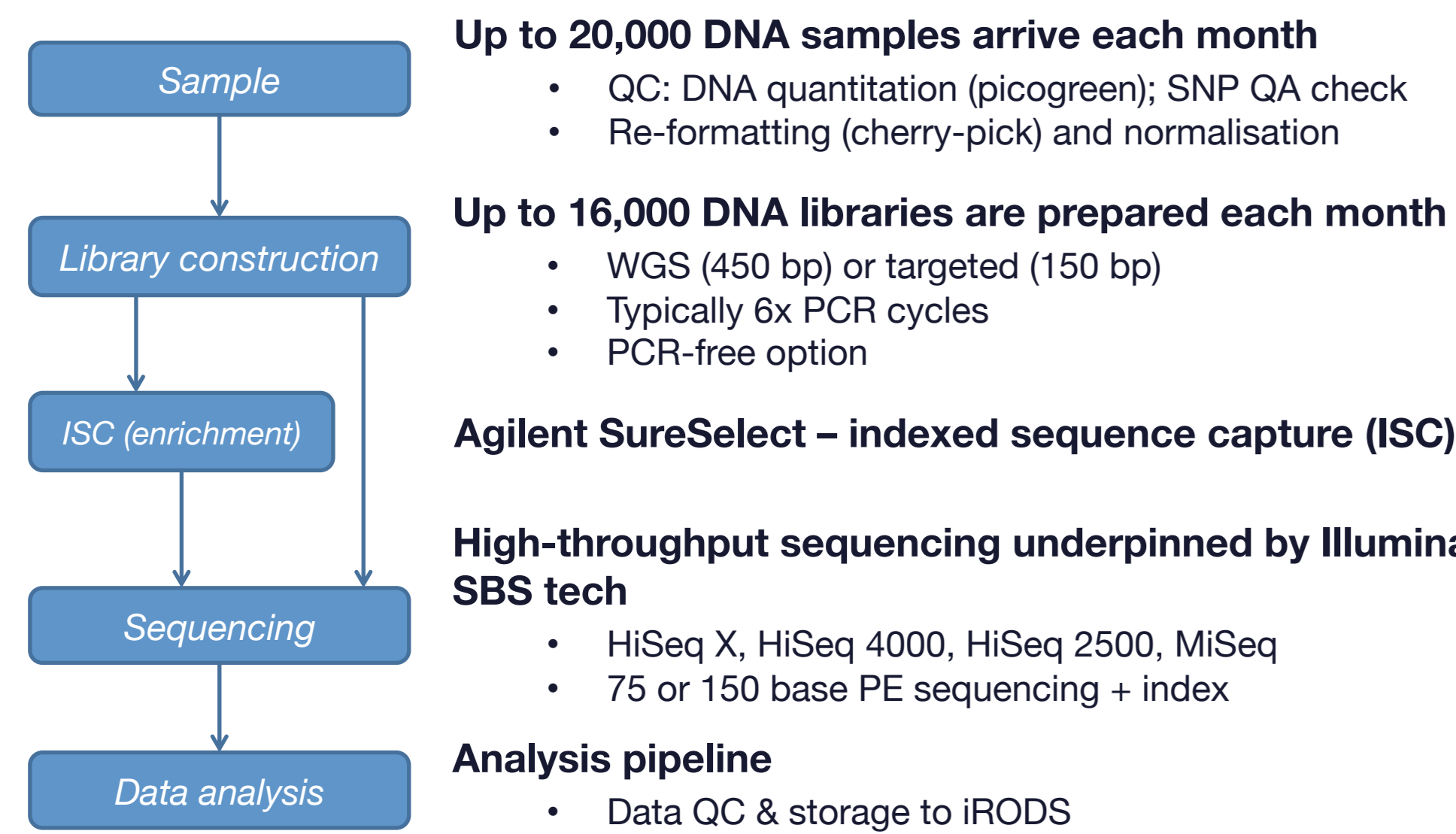
Wellcome Trust Sanger Institute, Hinxton, Cambridge, U.K.  
\*Presenting author



## Background

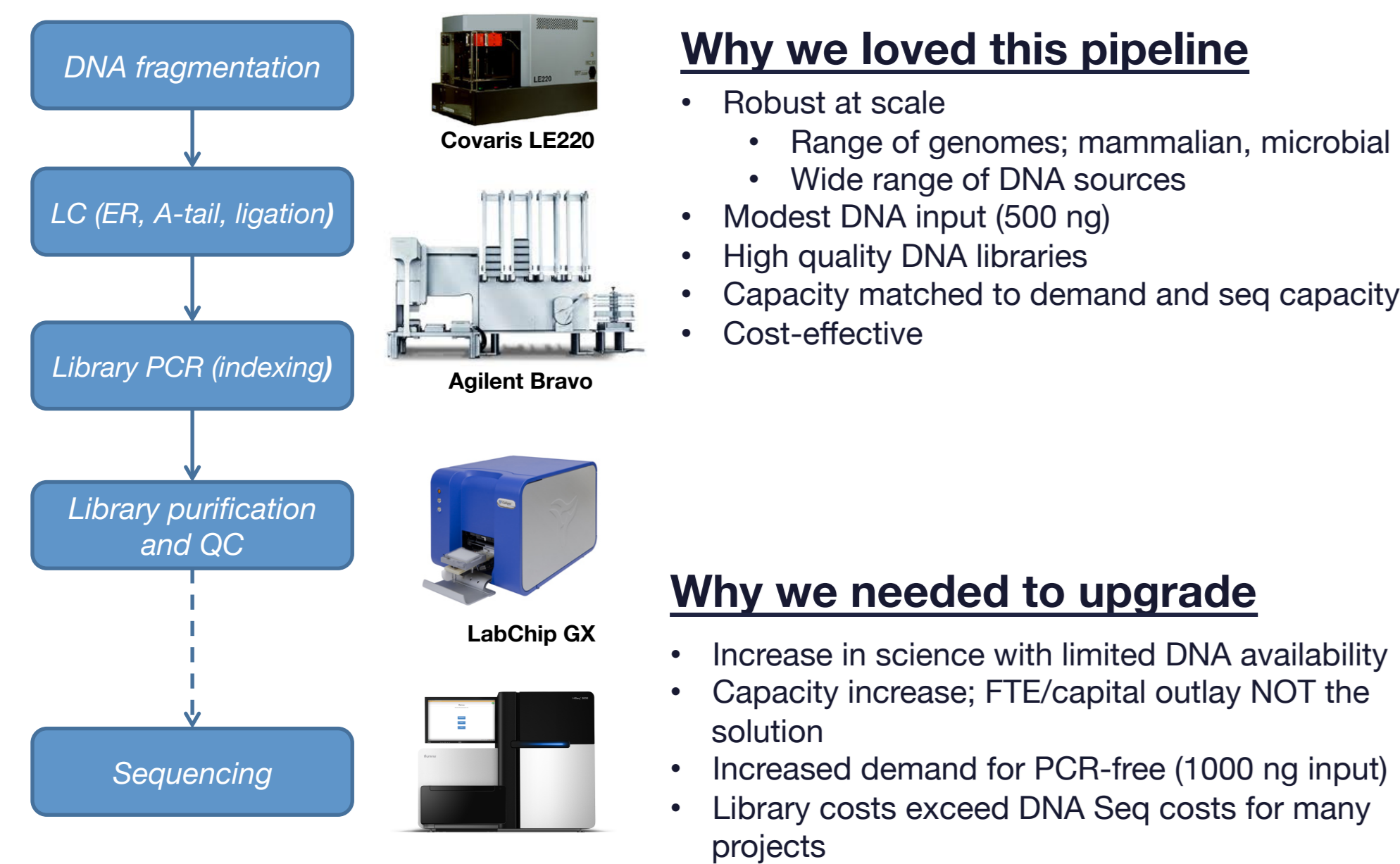
### Overview of High-throughput DNA Sequencing

The current operation employs 15 laboratory staff running multiple high-throughput Illumina sequencing pipelines to produce in excess of 1 petabase of data per year.



### DNA library construction (2010-2017)

Our high-throughput pipelines have been continually improved over the years but a step change in workflow design was necessary to prepare us for future projects.

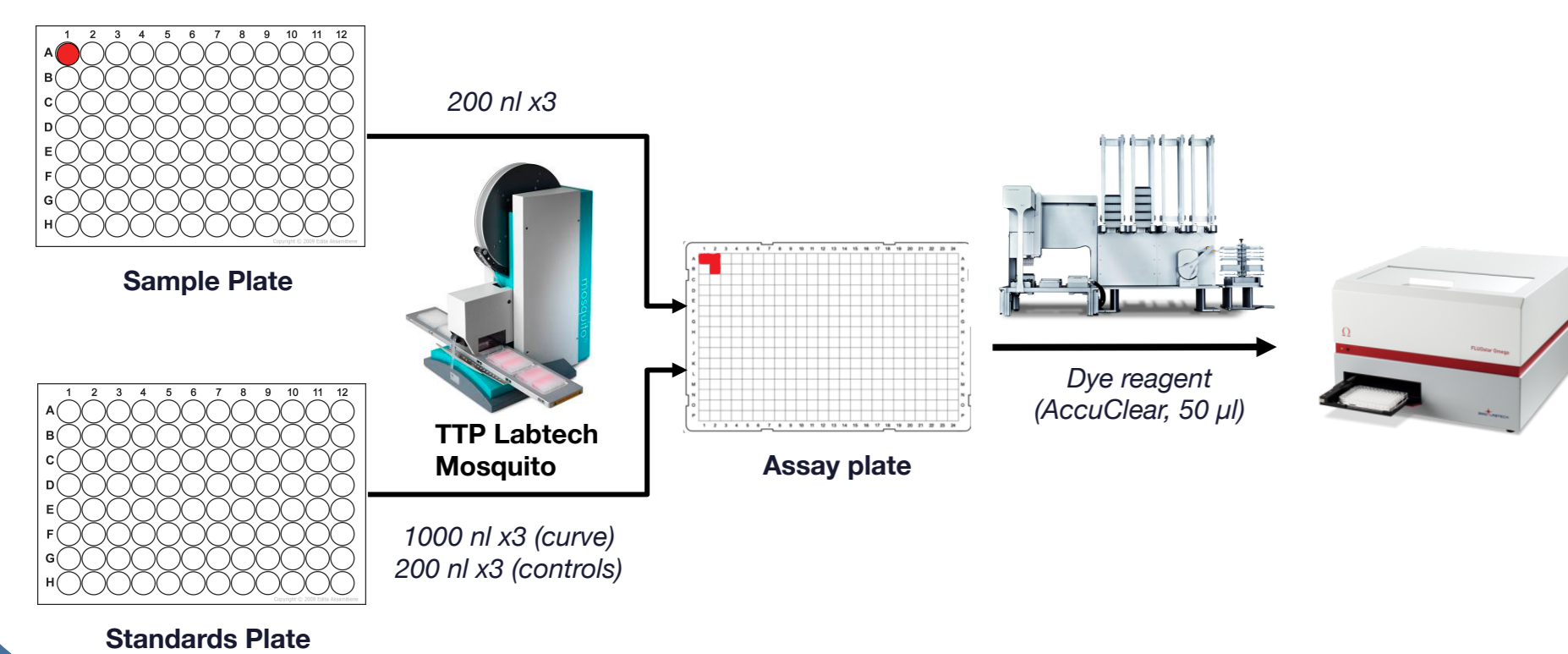


### Improving DNA quantitation: A key enabler

A key step towards overhauling and streamlining our workflows was to improve the quality of plates entering our high-throughput pipelines.

#### What are the benefits of this new QC?

- Less than 1 µl of sample is required
- Wider linear dynamic range than previous QC methods (0.03-200 ng/µl of stock DNA)
- Reduced turnaround time and costs compared to previous QC methods
- Positive displacement pipetting mitigates the viscoelastic properties of high molecular-weight DNA
- Improved cherry-picking and increased first-time pass rate



## PCR-Based LC

### The challenge

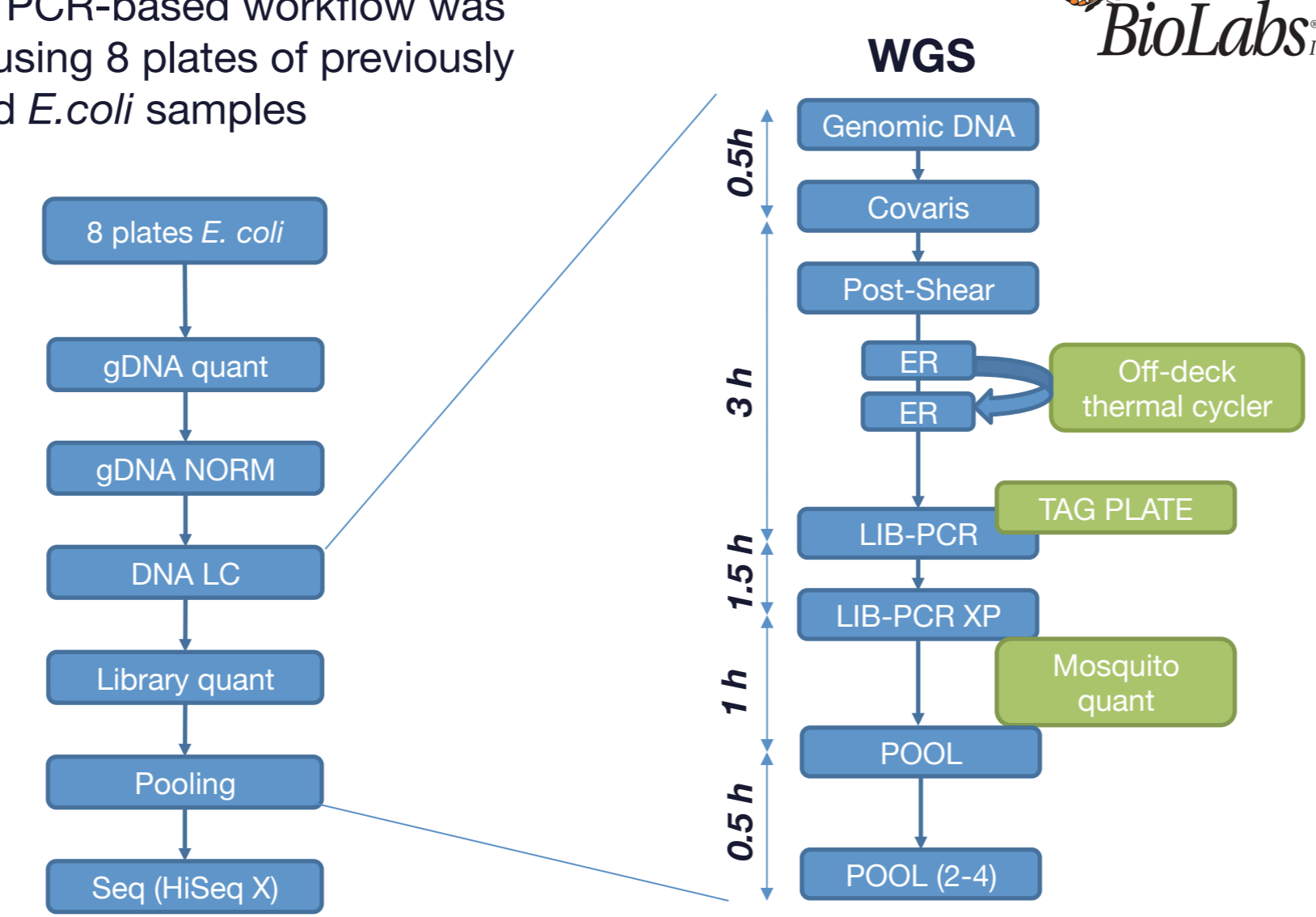
- Develop and implement a streamlined PCR-based LC process
- Reduce DNA input requirements by at least 50%
- Reduce per sample costs by ~50%
- No reduction in capacity nor increase in FTE requirement

### Development of a streamlined, automated PCR-based workflow

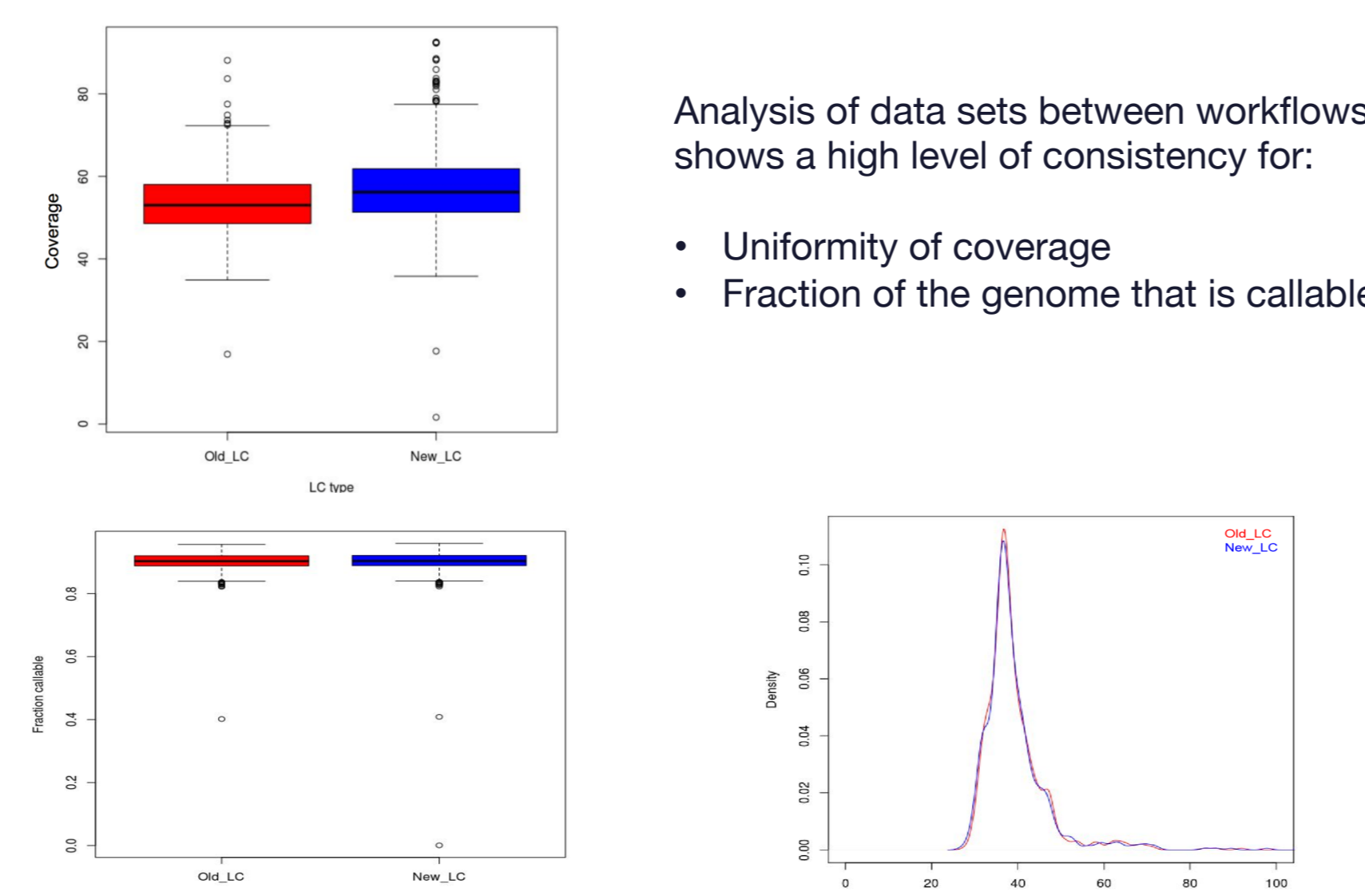
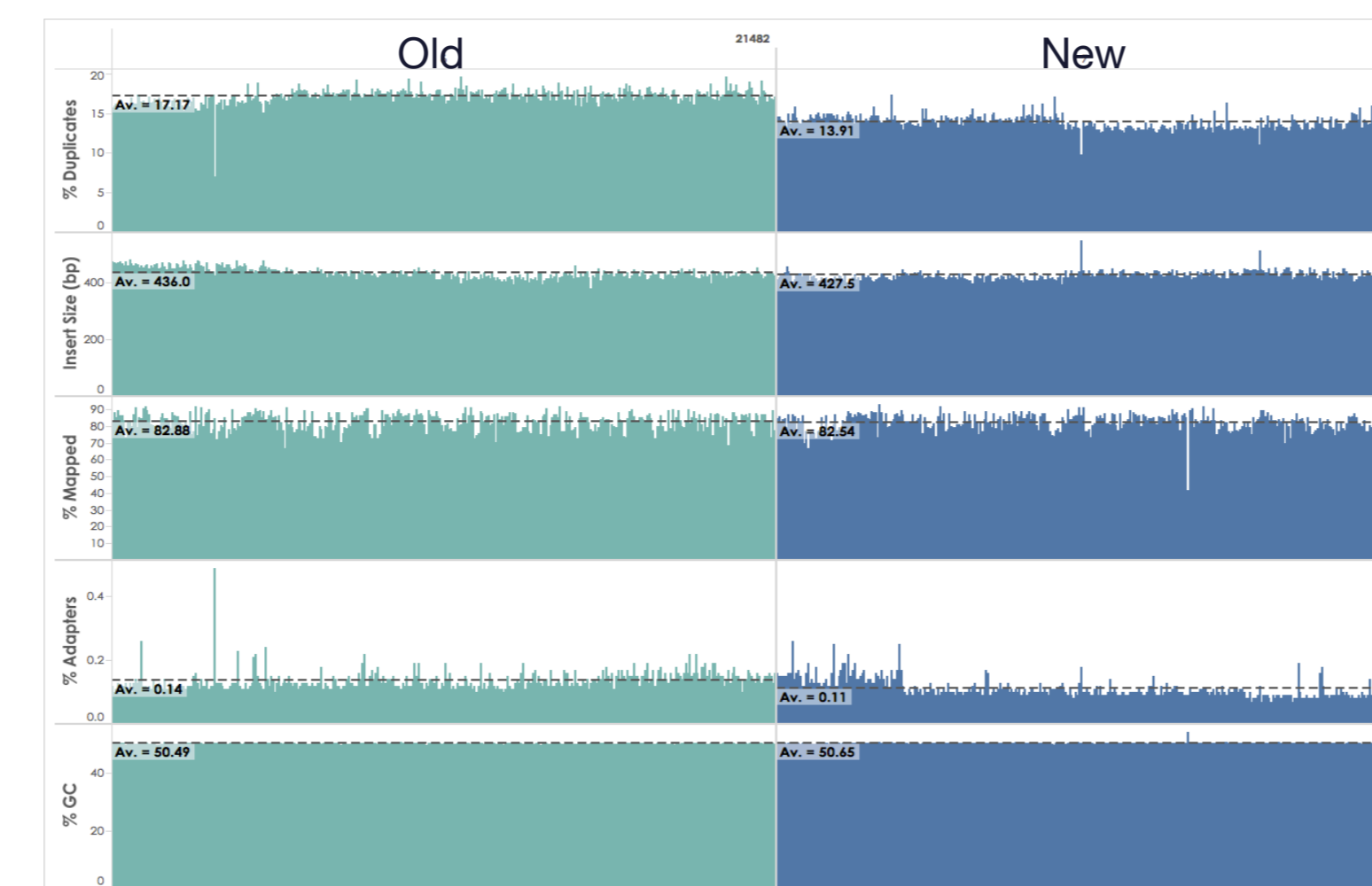
- New NEBNext UltraII workflow is ~30% faster enabling 2x PCR-based LC runs per Bravo per day
- Less demanding of FTE resource than the previous Sanger workflow
- Standard DNA input requirement reduced 2.5x to 200 ng
- Workflow is compatible with whole genome and targeted sequencing
- Streamlined DNA library quantitation utilising same method as novel gDNA assay

### LC Workflow Validation

The entire PCR-based workflow was validated using 8 plates of previously sequenced *E.coli* samples



- The new workflow met all pipeline metric performance targets
- Data comparison indicated no loss of quality with new, streamlined workflow



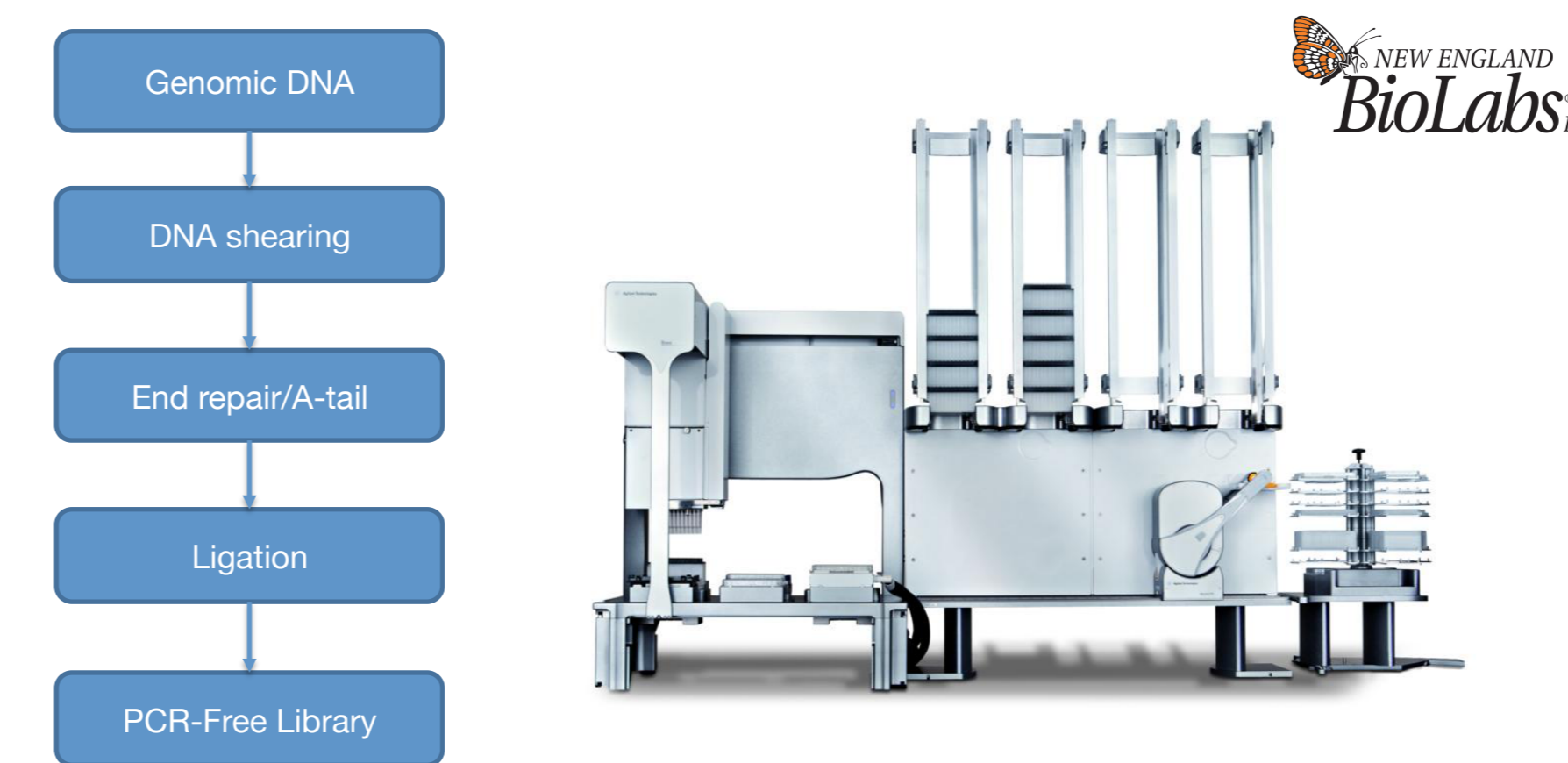
## PCR-Free LC

### The challenge

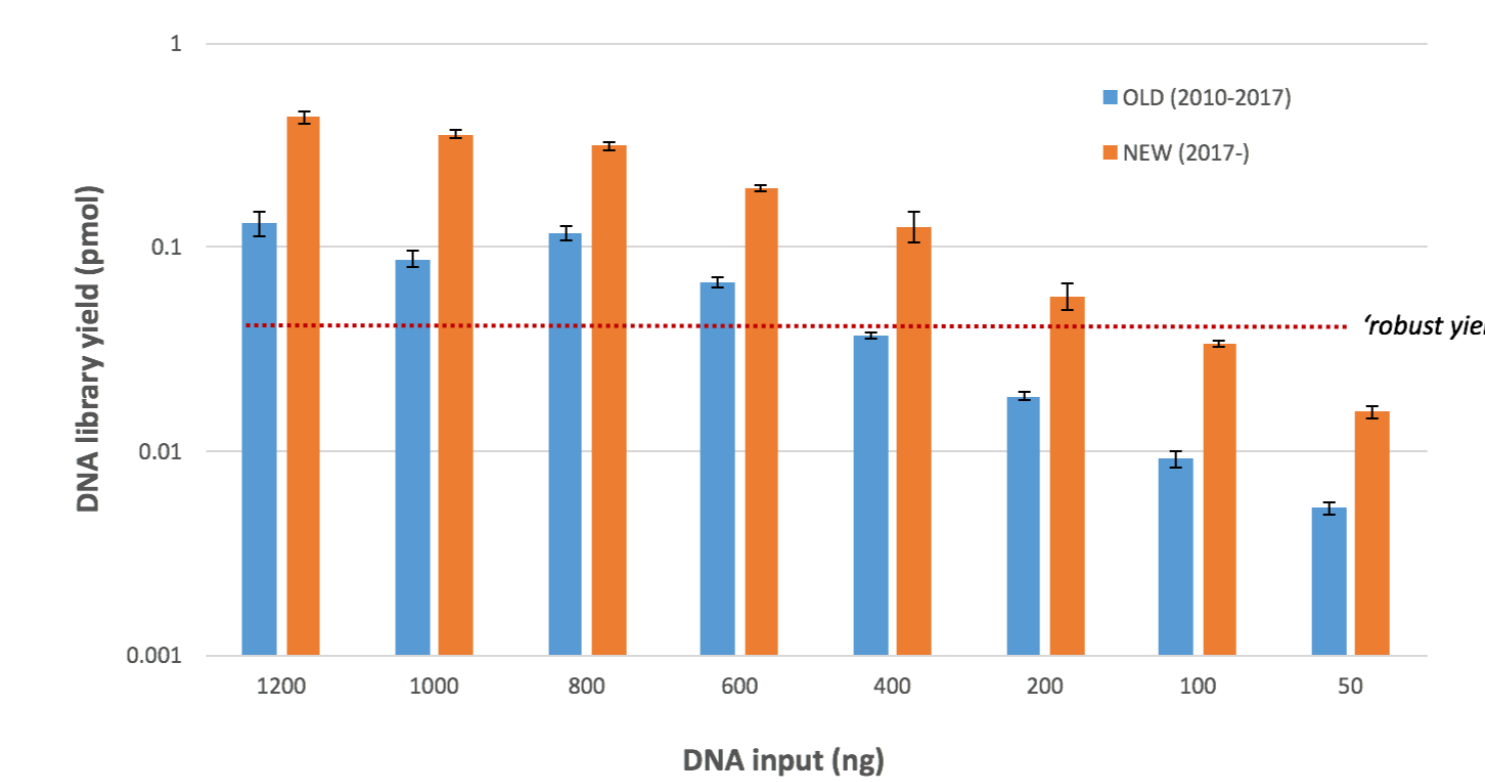
- Develop and implement a streamlined PCR-free LC process
- Reduce DNA input requirements by 50%
- No reduction in capacity nor increase in FTE requirement
- Align process with workflows implemented for PCR-based LC
- Implement Unique Dual Indexing (UDI)

### Development of a streamlined, automated PCR-Free workflow

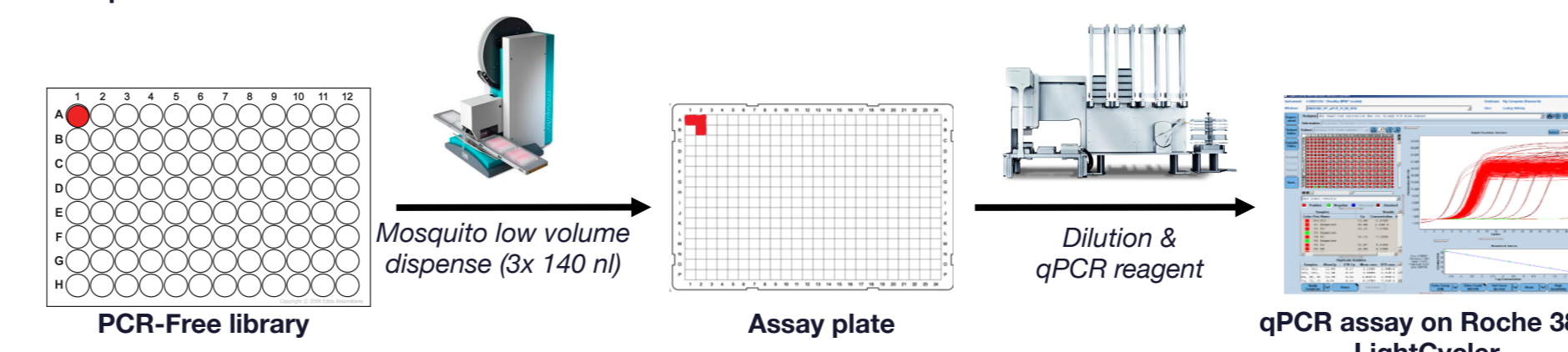
- New NEBNext UltraII PCR-Free workflow is 2x faster end-to-end and allows 2x PCR-Free LC runs per Bravo per day
- PCR-Free workflow adopts new gDNA quant and normalisation steps



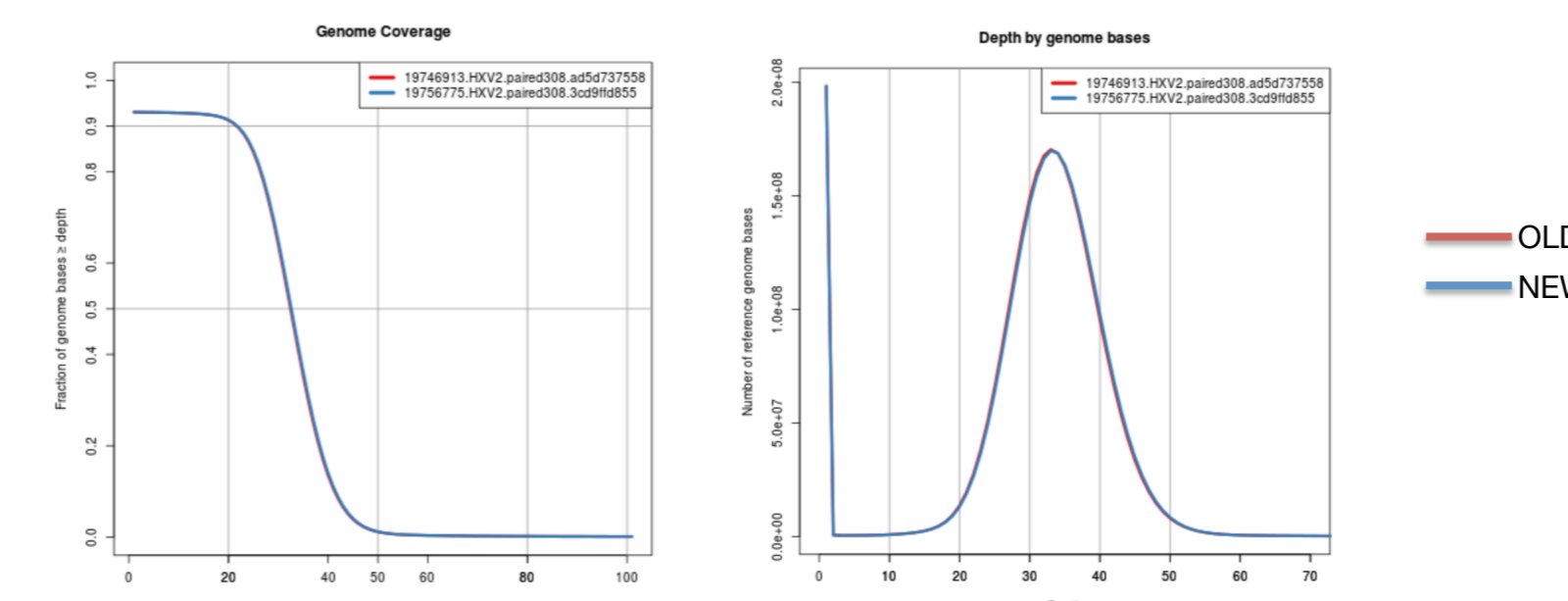
- DNA input requirements are reduced from 1000 ng to 500 ng



- Library QC is dramatically streamlined in new LC workflow (~5x faster)
- Low volume dispensing minimises loss of precious library to measurement process



- New PCR-Free workflow matches data quality of previous workflow
- Genotype concordance between methods >99%



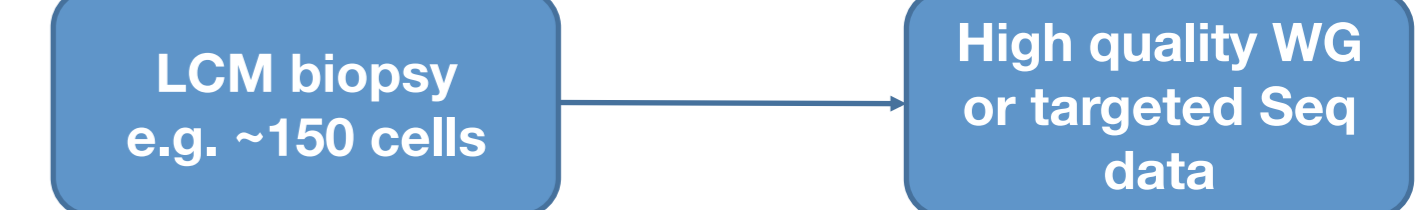
- Unique dual indexes (UDI) are incorporated in to each PCR-FREE library
- UDIs allow us to filter 'contaminating' reads generated by index hopping
- First release UDI set comprises 96 x 96 8-base barcoded adapters



## LCMB LC

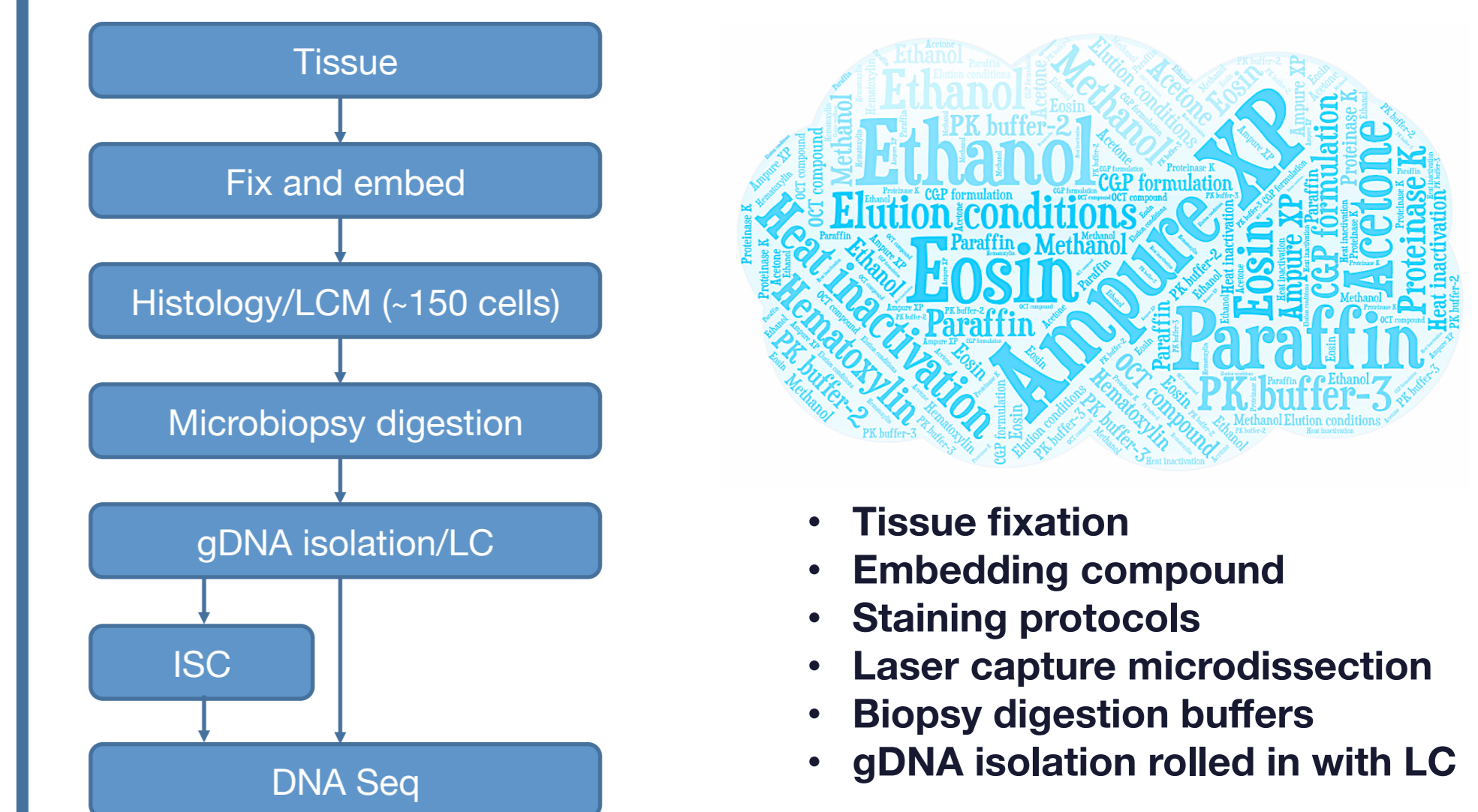
### The challenge

- Sanger's cancer and somatic mutation group aim to investigate clonal dynamics and mutational signatures in all tissues
- Requires laser capture microdissection (LCM) of just a few hundred cells
- DNA input is approximately 1 ng
- A high-throughput pipeline required for many thousands of samples
- Whole genome amplification methods NOT allowed



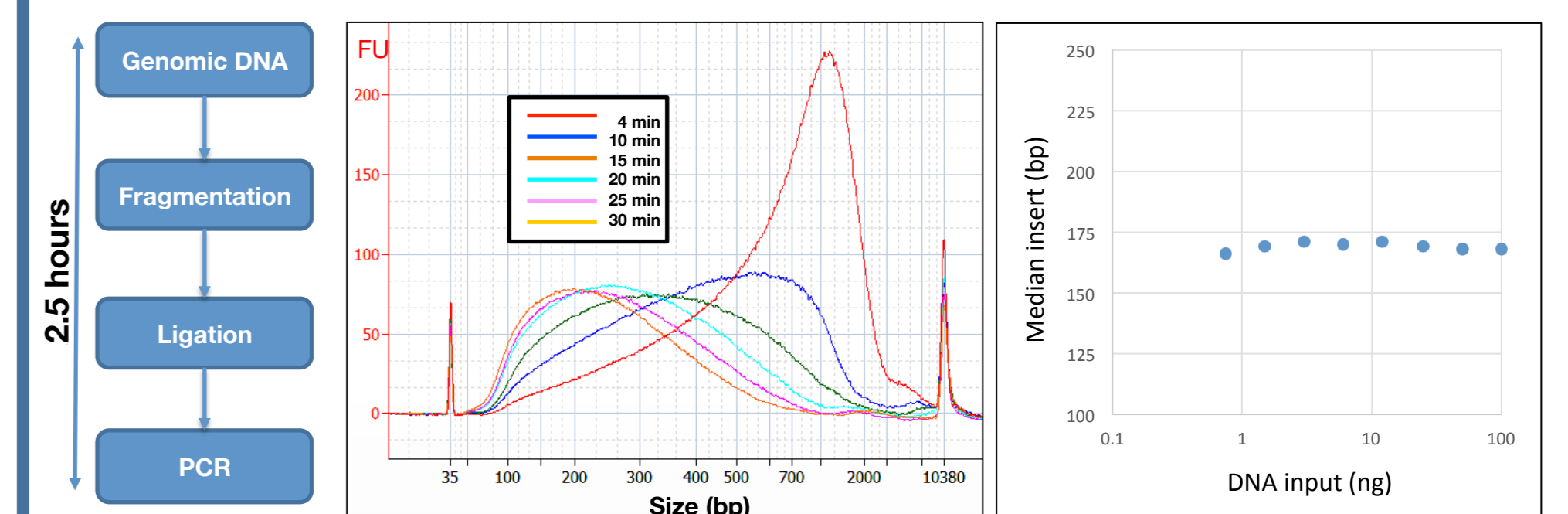
### Optimising LCM sample prep for low input LC workflow

- Sample prep and LC workflow were co-developed

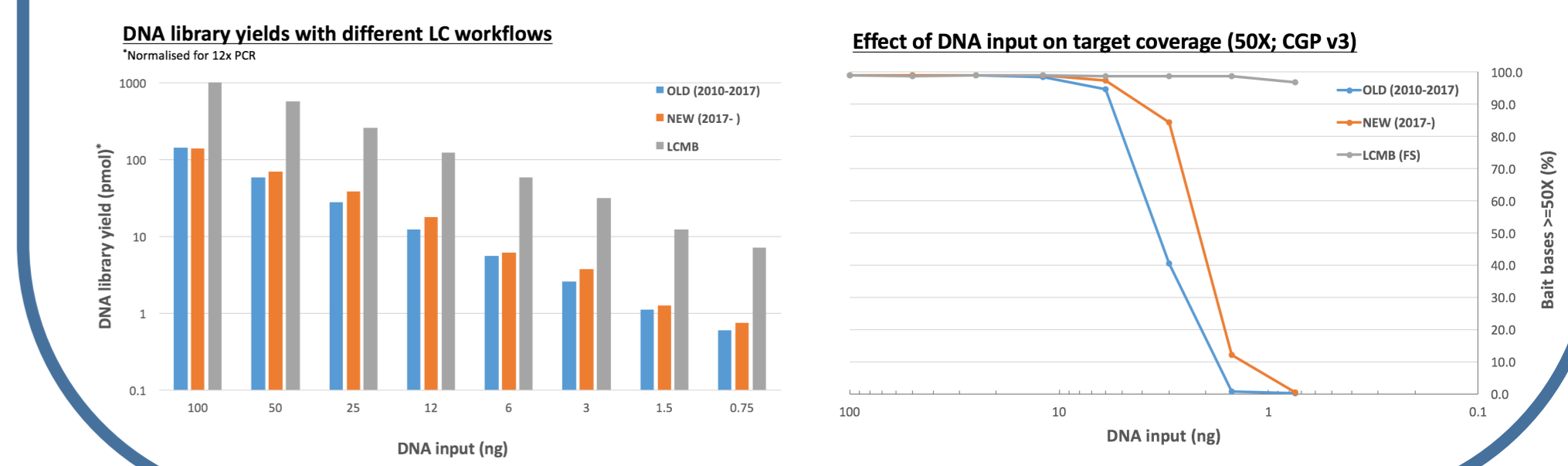


### Ultra low DNA input LC is enabled by NEBNext Ultra II FS

- Ultra II FS is a novel LC reagent from New England BioLabs (NEB)
- No requirement for Covaris shearing (enzymatic)
- Fragment profile is dependent on reaction time (e.g. 10 min for WGS profile)
- Fragment profile is independent of DNA input



- The LCMB pipeline produces 10-15x more DNA library compared with other methods
- DNA input requirements for high quality data are dramatically reduced



## Conclusions

2017 has seen a major overhaul of our high-throughput LC pipelines

- Improved quality of gDNA plates entering DNA library construction
- Faster, less expensive processes with no loss of data quality
- Reduction in DNA input requirements for PCR-based & PCR-Free workflows
- First steps to implementation of unique dual indexes for majority of DNA libraries
- Implementation of a new workflow (LCMB) capable of producing high quality whole genome and targeted human seq data from a few hundred cells