



Optimized low input library prep workflow for improved complexity and target coverage

Sample pre-qualification is not sufficient to increase the probability of successful preparation of sequencing libraries. To maximize the information output from FFPE samples, the SureSelect^{XT} protocols have been optimized, providing specific recommendations on amplification of pre-capture libraries as well as amount of sequencing to allocate per library based on the sample quality. These modifications ensure that there is sufficient representation of the molecules present in the starting sample going into the hybridization step which is critical to efficient enrichment of the targets. In addition, once these targets are enriched, the recommendations on sequencing depth should enable enough reads to ensure deep target coverage whether the starting sample is of higher or lower quality, for sensitive and accurate variant detection (**Figure 2**).

Complete cancer research solutions from sample to data

Optimized workflows are critical to providing comprehensive variant detection and reduced turn-around time from sample to data. The Agilent NGS FFPE QC kit completes the sample to sequencing workflow for FFPE samples. Combined with the efficient SureSelect^{XT} workflows that generate high complexity libraries appropriate for highly heterogenous samples, capture libraries optimized for enrichment of targets relevant in cancer research such as the ClearSeq Comprehensive Cancer panel and the SureSelect Human All Exon V6 + COSMIC, deep target coverage is achieved. For data analysis, SureCall enables analysis of paired tumor and normal samples, typical in cancer research, allowing for sensitive variant profiling of FFPE samples. (**Figure 3**).

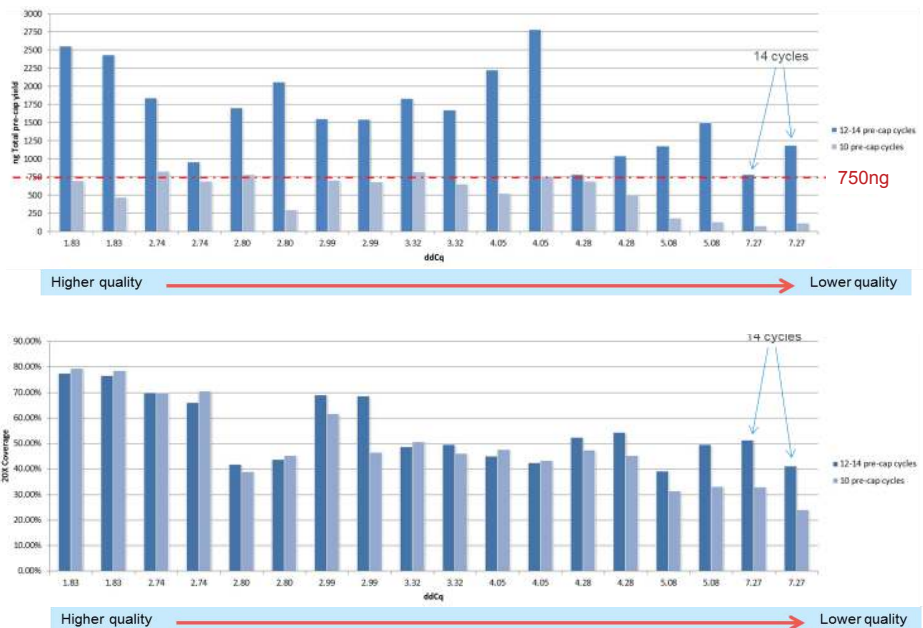


Figure 2. The Agilent NGS FFPE QC protocol provides recommendations to optimize preparation of enriched libraries for sequencing based on the quality score ($\Delta\Delta Cq$). Optimizations for lower quality FFPE samples include increasing amplification cycles for the pre-capture library to ensure sufficient template molecules are introduced into the hybridization, and increasing sequencing depth to enable better target coverage.

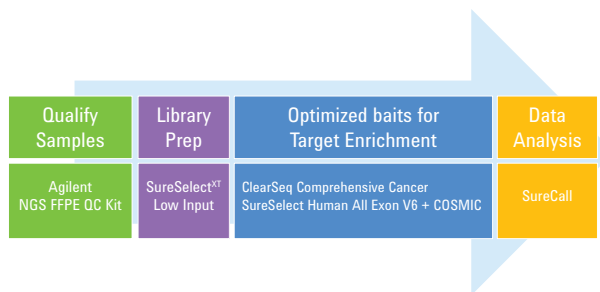


Figure 3. The Agilent NGS FFPE QC kit is part of the complete workflow solution that enables optimal data generation from even low quality FFPE samples. Complementary solutions for library prep, enrichment and data analysis are available providing support from sample pre-qualification all the way to variant calling.

Ordering Information:

	Part Number
Agilent NGS FFPE QC Kit, 16	G9700A
Agilent NGS FFPE QC Kit, 96	G9700B



Find an Agilent customer center in your country:

www.agilent.com/genomics/FFPEQC
U.S. and Canada: 1-800-227-9770
cag_sales-na@agilent.com

For Research Use Only. Not for Use in Diagnostic Procedures.

© Agilent Technologies, Inc. 2015
 Printed in USA, May 22, 2015
 5991-5941EN

