

# Agilent Femto Pulse System Kit Portfolio

## Quantitative DNA

Genomic DNA 165 kb  
1.3 kb to 165 kb

**FP-1002**

Smear  
5 pg/μL –  
500 pg/μL

Fragment  
0.3 pg/μL –  
30 pg/μL

Genomic Quality Number (GQN)

US NGS  
100 bp to 6,000 bp

**FP-1101**

Smear  
25 pg/μL –  
250 pg/μL

Fragment  
0.1 pg/μL –  
5 pg/μL

Fragment  
LOD  
50 fg/μL

## Qualitative DNA

55 kb BAC  
75 bp to 48.5 kb

**FP-1003**

Single Fragment  
1 pg/μL –  
25 pg/μL

Max input  
100 pg/μL

## Quantitative RNA

US RNA  
200 nt to 6,000 nt

**FP-1201**

15 pg/μL –  
250 pg/μL

LOD  
2.5 pg/μL

RNA Quality Number (RQN)  
DV<sub>200</sub>