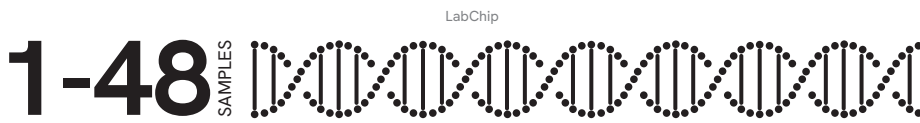
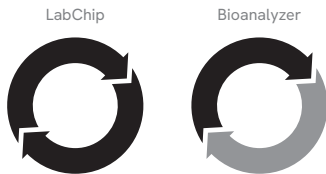


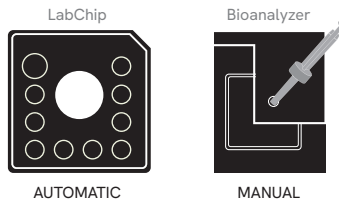
THROUGHPUT



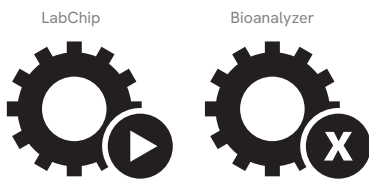
CHIP REUSABILITY



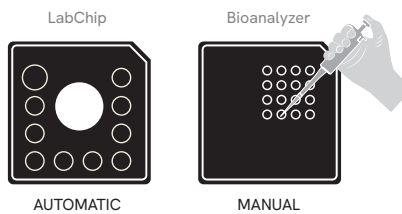
CHIP PRIMING



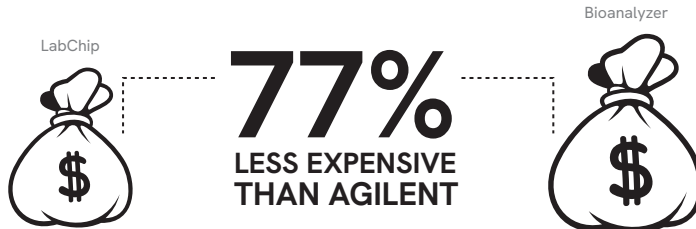
AUTOMATED SOLUTION



SAMPLE LOADING



COST PER SAMPLE
(12 SAMPLES)



Get more without spending more.

A comparison of the LabChip® GX Touch™ 24 Nucleic Acid Analyzer versus the Agilent Bioanalyzer instrument.

Learn more

