



Get more without spending more.

LabChip® GX Touch™ 24 nucleic acid analyzer

A walk-away solution for your nucleic acid analysis

With a flexible throughput, low cost of analysis, and a completely automated protocol that allows for walk-away analysis, the LabChip® GX Touch™ 24 Nucleic Acid Analyzer is the best choice for automated, low throughput nucleic acid analysis.

The LabChip GX Touch 24 Nucleic Acid Analyzer offers high-resolution analysis providing exact sizing and quantitation of DNA fragments, and smears down to 5bp from samples as low as 2pg/ul.

By automating the tedious, labor intensive chip priming, sample loading, and analysis steps using microfluidic technology, the LabChip GX Touch 24 Nucleic Acid Analyzer is the only nucleic acid analysis solution with a completely walk-away workflow, greatly reducing required hands-on time.

The LabChip GX Touch 24 Nucleic Acid Analyzer also offers a flexible throughput which allows for the analysis of 1 – 48 samples per prime.

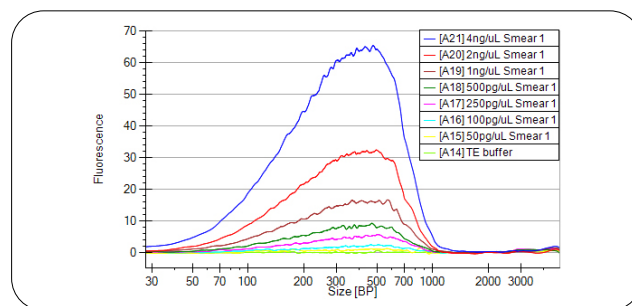
Because the chips are reusable and offer flexible throughput and affordable reagent pricing, the cost of analysis is lower using the LabChip GX Touch 24 Nucleic Acid Analyzer than other low throughput options.



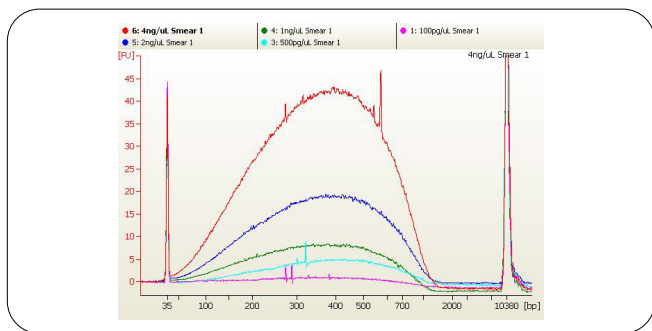
The LabChip GX Touch 24 Nucleic Acid Analyzer also offers the flexibility to grow with your lab. If your throughput increases as your lab grows, the LabChip GX Touch 24 Nucleic Acid Analyzer can easily be upgraded to analyze up to 384 samples at a time.

Contact your local Revvity sales representative to find out how you can get more, without spending more.

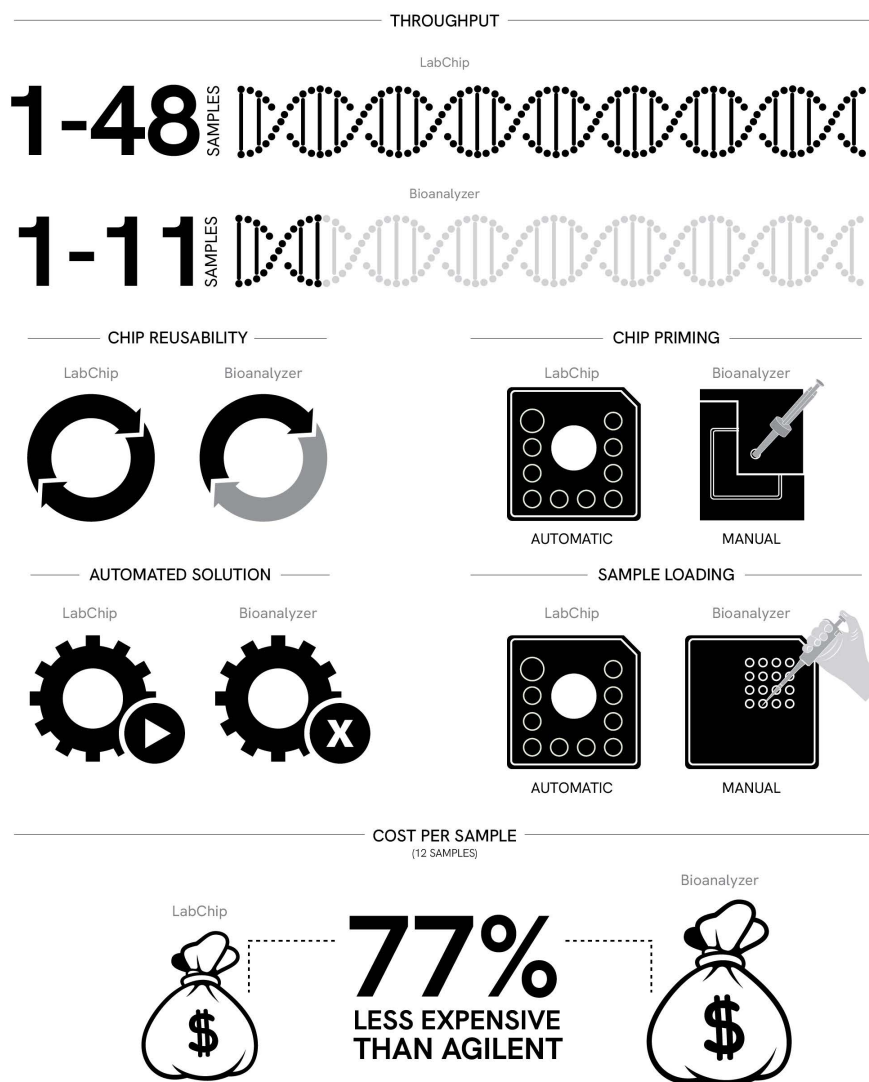
LabChip trace



Bioanalyzer trace



Nominal conc. (ng/μL)	GX conc. (ng/μL)	BA conc. (ng/μL)
0.05	0.033	-
0.1	0.096	0.158
0.5	0.472	0.461
1.0	1.0	0.958
2.0	2.1	1.8
4.0	4.3	3.8



revvity