

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 formatting, 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 5 bp from samples as small as 2 pg/μL.

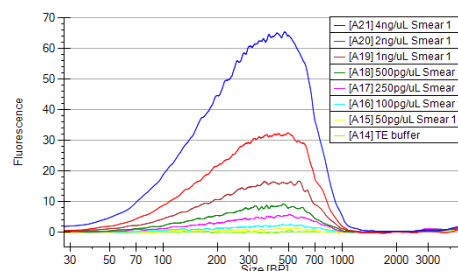
By automating the tedious, labor intensive chip priming, sample loading, and analysis steps using microfluidic technology, the LabChip GX Touch 24 Nucleic Acid Analyzer is the only nucleic acid analysis solution with a completely walk-away workflow, greatly reducing required hands-on time.

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

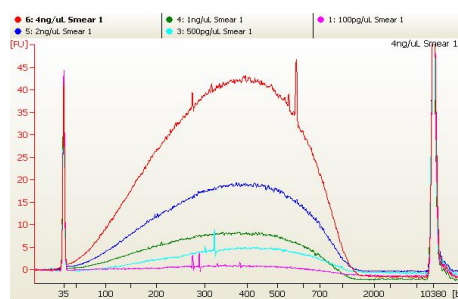
Because the chips are reusable and offer flexible throughput and affordable reagent pricing, the cost of analysis is lower using the LabChip GX Touch 24 Nucleic Acid Analyzer than other low throughput options.

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 PerkinElmer sales representative to find out how you can get more, without spending more.



LabChip trace



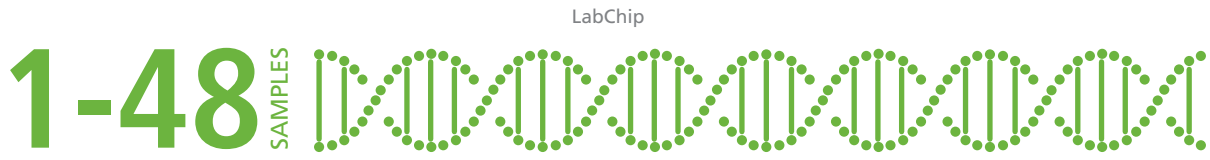
Bioanalyzer trace

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

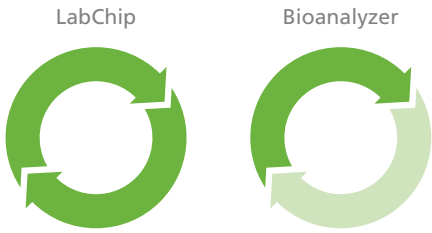
For research use only. Not for use in diagnostic procedures.



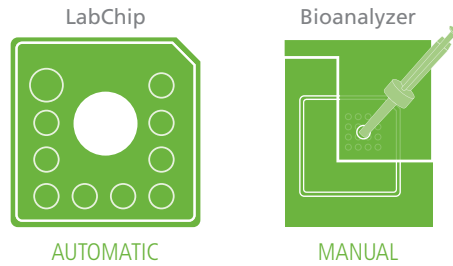
THROUGHPUT



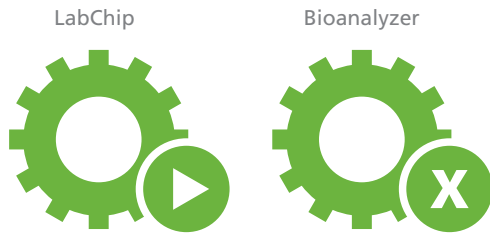
CHIP REUSABILITY



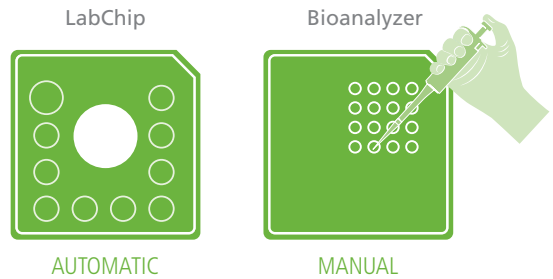
CHIP PRIMING



AUTOMATED SOLUTION



SAMPLE LOADING



COST PER SAMPLE
(12 SAMPLES)



PerkinElmer, Inc.
 940 Winter Street
 Waltham, MA 02451 USA
 P: (800) 762-4000 or
 (+1) 203-925-4602
www.perkinelmer.com



For a complete listing of our global offices, visit www.perkinelmer.com/ContactUs

Copyright © 2017, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

AG011709_20A_FL Printed in USA