



Reimagine low-throughput library prep.

BioQule™ NGS System

EASY TO USE, COMPACT SOLUTION FOR AUTOMATING LOW THROUGHPUT LIBRARY PREP

Reimagine low-throughput library prep with the BioQule™ NGS System and simplify your library prep and quantitation with walk-away automation. This low-cost, open, benchtop instrument delivers libraries ready to load into your sequencer in only 15 minutes of hands-on time (HOT), allowing you to focus on what really matters.

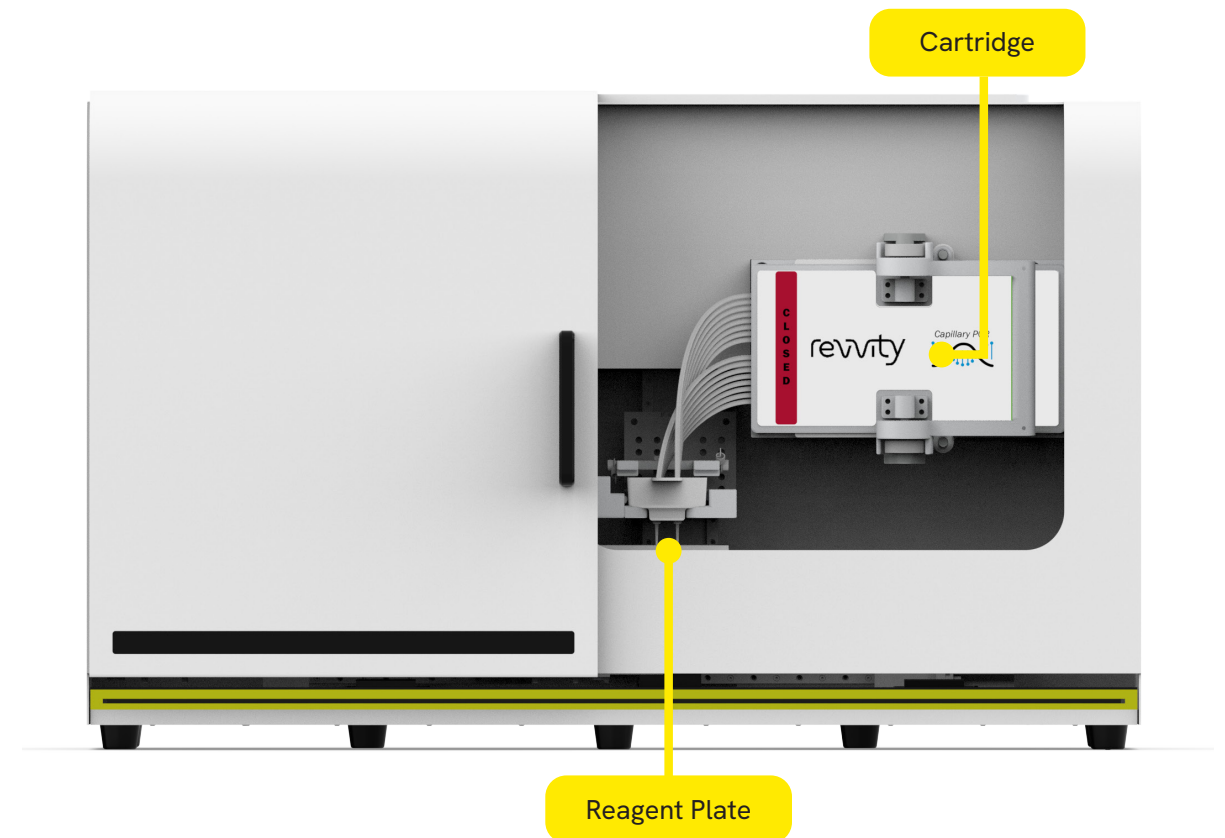
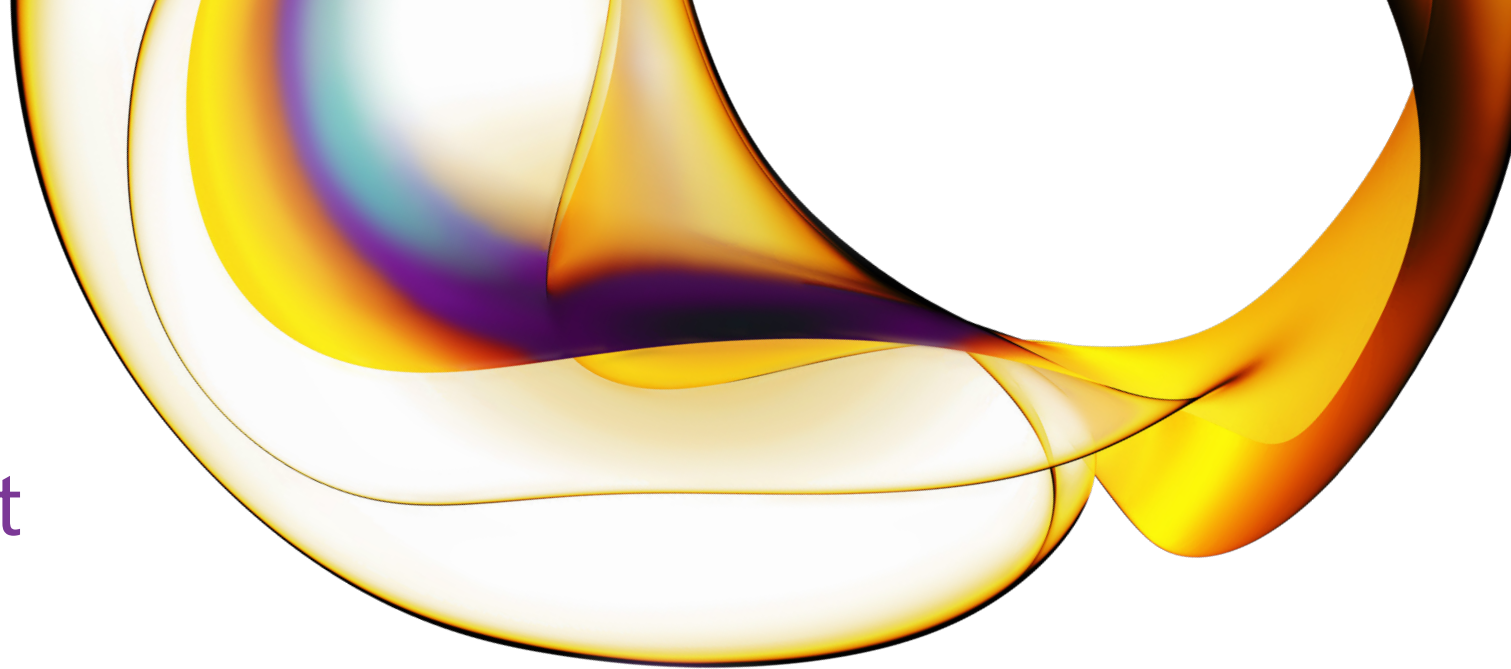
Currently, automated methods are available for NEXTFLEX®, Illumina® and Roche® KAPA® kits.

- Full walkaway automation (80% reduction in HOT)
- Easy to use (only 4 touch points)
- Minimize human error (>99% success rate)
- Open system able to automate different applications



Learn how to automate your low-throughput library prep.

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

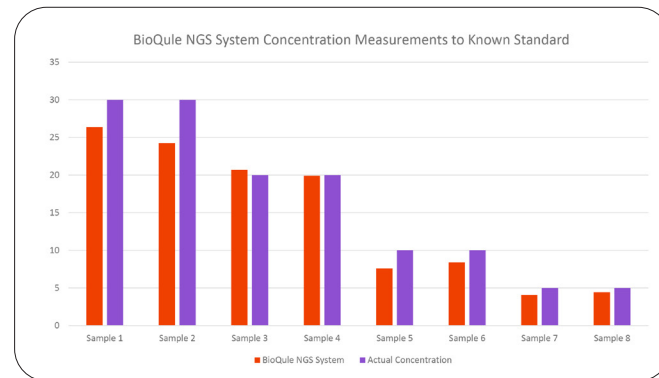


Easy NGS library preparation

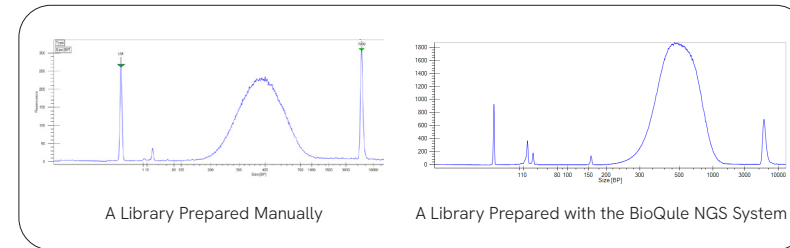
The BioQule NGS System has been specifically designed so no automation experience is required to generate NGS libraries. Simply pipette your samples, beads, and two reagents into the plate, place the cartridge and plate in the instrument, close the door and walk-away. The BioQule NGS System will construct AND quantitate up to 8 libraries while you to focus on what really matters.

Library preparation and quality control in the same system

The BioQule NGS System is equipped with automated fluorescence-based optical quantitation to measure the concentration of each library. You will no longer need to move your samples from one instrument to another to perform concentration analysis on your libraries.



The BioQule NGS System constructs libraries with size distributions that are equivalent to those prepared manually.



Simplified reagent and cartridge preparation

Each run of the BioQule NGS System uses one reagent plate and one cartridge. The reagent plate comes pre-plated with most reagents required for NGS library preparation. Once one run is complete, simply change out the plate and cartridge and you can construct another 8 NGS libraries.



BioQule NGS System Compatible Kits*

- NEXTFLEX® Rapid DNA-Seq kit 2.0
- NEXTFLEX® Rapid XP V2 DNA-Seq kit

Other automated methods for the BioQule NGS System*

- Illumina® DNA Prep protocol
- Roche® KAPA® HyperPrep™ protocol

** More kits and automated methods are in development. Please inquire.

Specifications



ELECTRICAL REQUIREMENTS:
100-240VAC at 50-60Hz



WEIGHT:
45 lbs

revvity