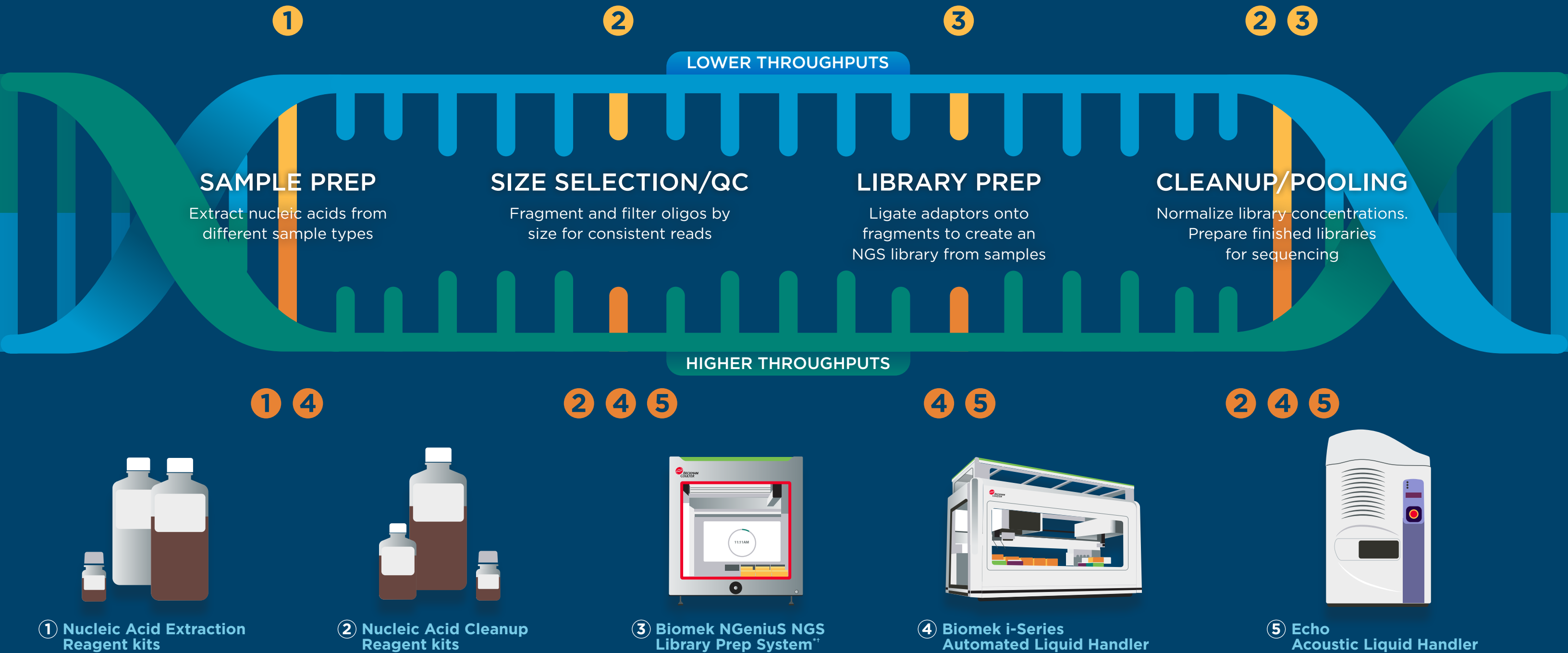


# NGS Library Prep

Next Generation Sequencing utilization is increasing at a rapid pace. As the cost of sequencing decreases, NGS is becoming a common technique for knowledge discovery and disease research. No matter the throughput of your laboratory, NGS library prep takes time and is costly. For lower throughputs, automating with Biomek NGenius allows for a completely automated, walk-away library prep solution. At higher throughputs, combining the flexibility of a Biomek i-Series with the speed of an Echo Acoustic Liquid Handler will save time and reagent costs. No matter your desired throughput, Beckman Coulter Life Sciences has an efficient solution.



\*Product in development. Performance characteristics have not been validated.  
\*\*Normalization may require additional equipment. Biomek NGenius pooling capability not available at launch.