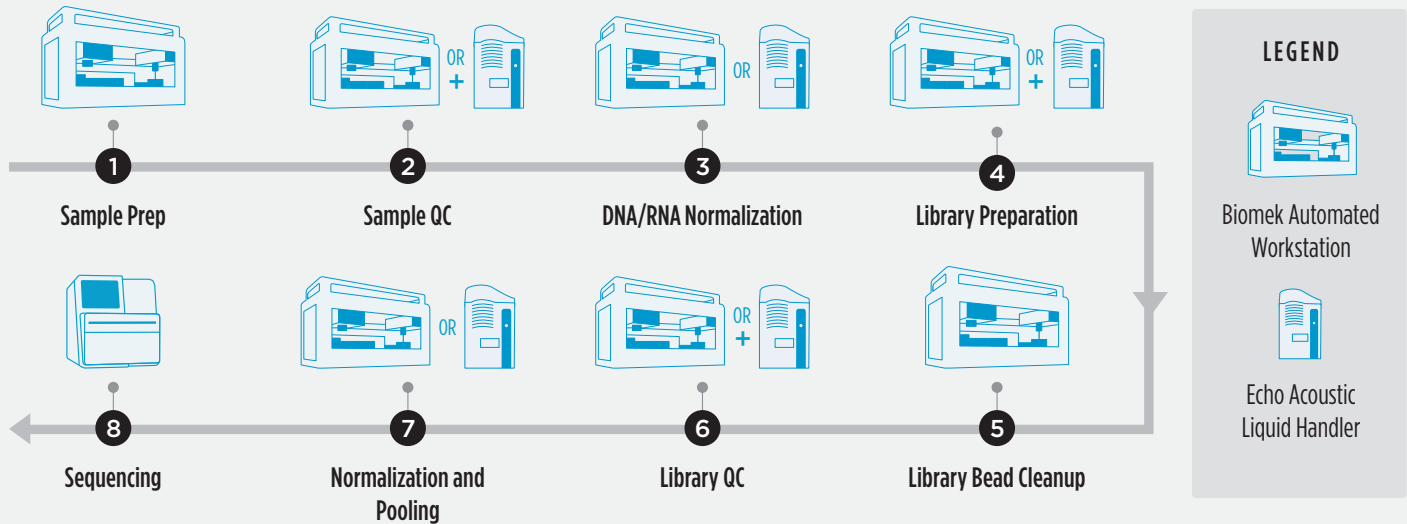


Generate Accurate Data Faster by Automating Your NGS Workflow



1 Sample Prep

Use a Biomek Workstation and our bead-based reagents to automate DNA and RNA extraction.

2 Sample QC

Use either a Biomek Workstation or a combination of Biomek Workstation and Echo LH, depending on your workflow needs.

3 DNA/RNA Normalization

Use either a Biomek Workstation or an Echo LH depending on your workflow needs.

4 Library Preparation

For standard volumes, use a Biomek Workstation or a combination of an Echo LH and Biomek Workstation (i.e., reduce reaction volumes of precious reagents using the Echo LH while doing larger volume transfers on the Biomek Workstation).

5 Library Bead Cleanup

Use a Biomek Workstation for 96- and 384-well bead cleanups.

6 Library QC

Use either a Biomek Workstation, or an Echo LH and Biomek Workstation depending on your workflow needs.

7 Normalization and Pooling

Use a Biomek Workstation for preparing the dilution plates and pooling of the libraries. Alternatively, use an Echo LH for simultaneous normalization and pooling, for up to 97% faster processing time.

8 Sequencing

Load the pooled libraries onto a flow cell of a sequencer.

The method/application illustrated is for demonstration only, and is not validated by Beckman Coulter. Beckman Coulter makes no warranties of any kind whatsoever express or implied, with respect to this protocol, including but not limited to warranties of fitness for a particular purpose or merchantability or that the protocol is non-infringing. All warranties are expressly disclaimed. Your use of the method is solely at your own risk, without recourse to Beckman Coulter. Not intended or validated for use in the diagnosis of disease or other conditions.

Products identified are not for use in diagnostics procedures.

© 2022 Beckman Coulter, Inc. All rights reserved. Beckman Coulter, the stylized logo, and the Beckman Coulter product and service marks mentioned herein are trademarks or registered trademarks of Beckman Coulter, Inc. in the United States and other countries. ECHO is a trademark or registered trademark of Labcyte Inc. in the United States and other countries.

All other product names and brands are properties of their respective owners.

For Beckman Coulter's worldwide office locations and phone numbers, please visit Contact Us at [beckman.com](https://www.beckman.com)

21.08.2274.AUTO