



# Biomek NGenius

NEXT GENERATION LIBRARY PREP SYSTEM

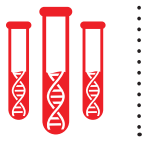
IT'S NOT JUST SMART. IT'S **INGENIOUS.**



You asked for an accurate, easy to use, cost-effective, NGS sample prep system with built-in error reduction and the flexibility to run an array of library preparation kits from multiple vendors—without requiring in-depth programming knowledge, and now it's here.

ACCELERATING  
*answers.*

 **BECKMAN  
COULTER**  
*Life Sciences*



# Introducing Biomek NGeniusS System

## The **Genius** is in its Total Workflow Solution



### 1. Flexible application selection

As part of our comprehensive service agreement, the Biomek NGeniusS System provides a wide range of complimentary demonstrated methods from a variety of reagent manufacturers, including:

#### DNA Sequencing

- Whole genome
- Targeted re-sequencing (Amplicon)

#### RNA Sequencing

- mRNA
- Total RNA
- Targeted RNA (Amplicon)



### 2. Application control

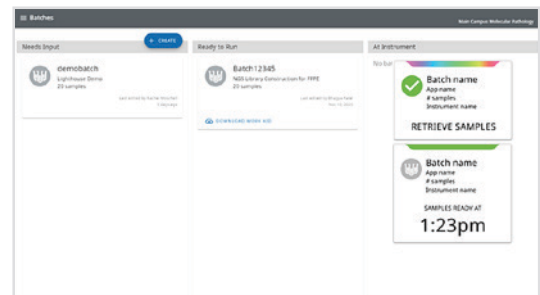
Control user access to applications with a simple authorization process (if required). Ensure that only authorized users have access to non-lab validated applications.



### 3. Work-from-anywhere batch set up\*

Using the **Biomek NGeniusS Portal software**, set up your batch run when and where it's most convenient for you.

Whether working from home, office, or benchtop, you can program batch sizes from 4-24 samples.



### 4. Reagent preparation made easy

Once you select the application and enter the number of samples, the Biomek NGeniusS software creates a **Work Aid**, which includes:

- Reagent list needed for the run
- Volume requirements for each reagent
- Instructions on how to prepare and load reagents

The Work Aid saves time and effort while ensuring reproducibility during reagent preparation.

REAGENT PREPARATION	REAGENT	REQUIRED
1. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
2. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
3. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
4. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
5. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
6. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
7. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
8. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
9. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
10. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
11. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
12. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
13. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
14. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
15. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
16. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
17. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
18. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
19. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
20. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
21. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
22. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
23. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
24. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
25. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
26. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
27. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
28. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
29. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
30. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
31. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
32. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
33. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
34. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
35. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
36. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
37. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
38. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
39. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
40. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
41. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
42. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
43. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
44. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
45. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
46. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
47. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
48. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
49. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
50. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
51. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
52. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
53. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
54. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
55. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
56. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
57. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
58. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
59. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
60. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
61. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
62. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
63. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
64. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
65. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
66. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
67. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
68. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
69. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
70. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
71. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
72. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
73. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
74. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
75. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
76. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
77. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
78. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
79. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
80. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
81. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
82. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
83. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
84. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
85. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
86. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
87. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
88. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
89. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
90. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
91. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
92. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
93. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
94. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
95. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
96. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
97. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
98. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
99. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL
100. Pipette 100 µL of each reagent into the designated wells.	100 µL	100 µL



## 5. Error-reduced instrument setup

Sophisticated optical analytics combined with the head-up display give you real-time feedback on labware placement, virtually eliminating loading errors.

Now you can allow less-experienced operators to run complex applications with confidence.



## 6. Remote system monitoring\*

The 360° multicolored, pattern-encoded light bar shows current instrument status.

Not within visual range? The Biomek NGenius Portal software allows you to monitor the system status from any computer.



## 7. Simplified sample processing

The **Multi-functional Reaction Vessel (RV)** combines three consumables into one, reducing time spent on labware inventory management:

- Sample input plate
- Thermal cycling plate
- Clean-up plate

Track your samples—with up to **100 million unique RV barcodes**—throughout the application run.

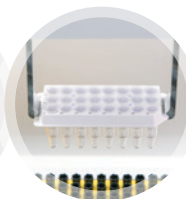


\* Requires a Google Chrome or Microsoft Edge enabled computer.

# Biomek NGenius System

## The latest in a long line of Smart Automation Technology

Inspired by Beckman Coulter's 30-plus years of laboratory automation and Smart Automation Technology, Biomek NGenius System represents an evolutionary leap in NGS liquid-handling systems with its best-in-class performance, error-reduction technology, ease-of-use, reduced hands-on-time, and flexibility.

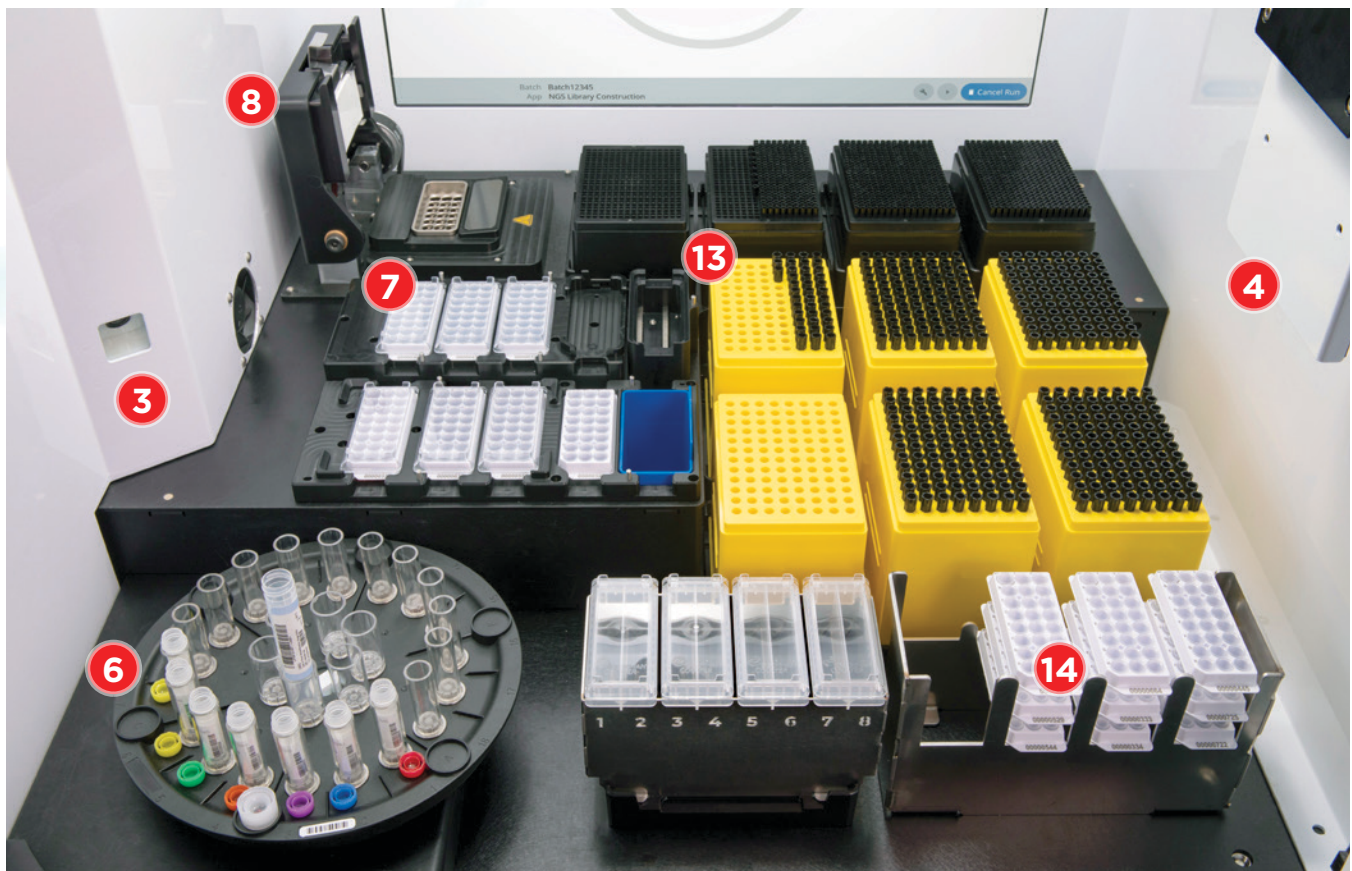


### Error reduction

- 1 Dynamic DeckOptix** system provides software-guided labware placement and confirmation
- 2 Head-up Display** ensures mission critical information is in sight
- 3 Reagent Carousel Auto ID** confirms the correct reagents are on board for the samples being processed
- 4 Multi-channel Liquid Level Sensing** aids in the detection of insufficient reagents, alerting the operator before it becomes a processing issue
- 5 Sample Tracking** eliminates the worry of mixing up samples when running in pre- and post-PCR lab environments

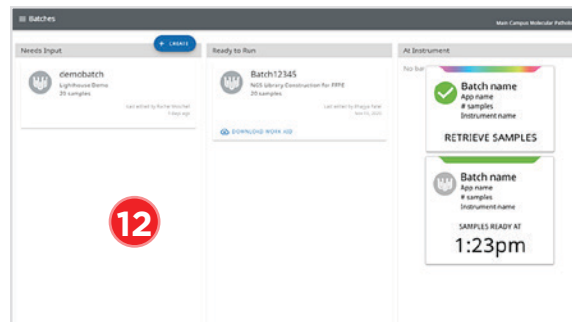
### Less hands-on-time

- 6 Reagent Carousel** reduces the need to manually transfer reagents and automates most reagent kits designed for manual use
- 7 Temperature-controlled Reagent Storage** protects reagents from thermal degradation and allows you to run end-to-end chemistries
- 8 Integrated Thermal Cycler** extends walkaway time
- 9 Multi-functional Pipetting Pod**
  - **Pipetting Range** - 1 to 1000  $\mu$ L
  - **Selective Tip Type** - ensures the right tip for the right volume without manual intervention
  - **Select Tip** - 1 to 8 tips for pipetting efficiency
  - **Labware Transport** - enables longer walkaway time
- 10 UV-C Light** standard on the system



## Easier to use

- 11 **Software Navigation Dial** makes system programming as simple as turn and select
- 12 **Biomek NGenius Portal Software\*** allows setup and monitoring of batch runs from wherever it's most convenient for you
- 13 **Partial Tip Box Use** eliminates manual tip transfers and/or wasting partially used tip boxes
- 14 **Multi-functional Reaction Vessel** provides 3-in-1 functionality and less inventory management
- 15 **Multicolor Status Light** provides 360° line of sight system status



## Flexible design

- ✓ **Dynamic Sample Batch Size** accommodates batch sizes of 4 to 24 samples
- ✓ **Complimentary Application Library** is designed to evolve with your changing chemistry needs
- ✓ **Network Enabled Scalability** enables you to seamlessly set-up and run multiple instruments from your Biomek NGenius portal.
- ✓ **Continuing Application Development** ensures your investment is protected for the foreseeable future

\* Requires a Google Chrome or Microsoft Edge enabled computer.



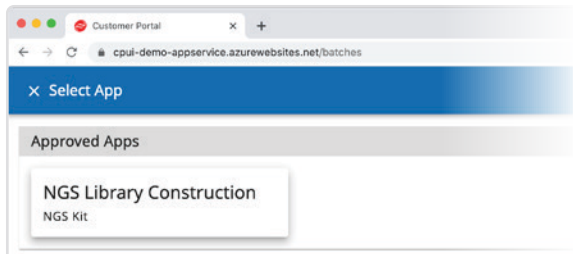
# The **Genius** is in its **Simplicity**

The unique instrument design and software functionality of the **Biomek NGenius** System not only make it easy to operate, but advanced programming skills are not necessary to set up an application and complete a run.

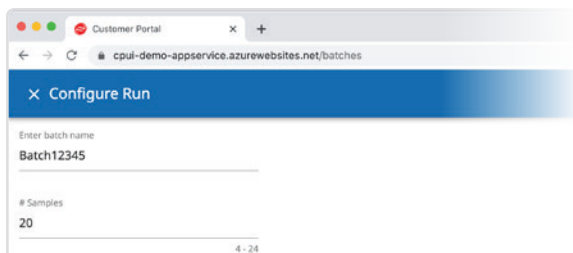
Set up and run a batch in 6 easy steps...

On the Biomek NGenius Portal software:

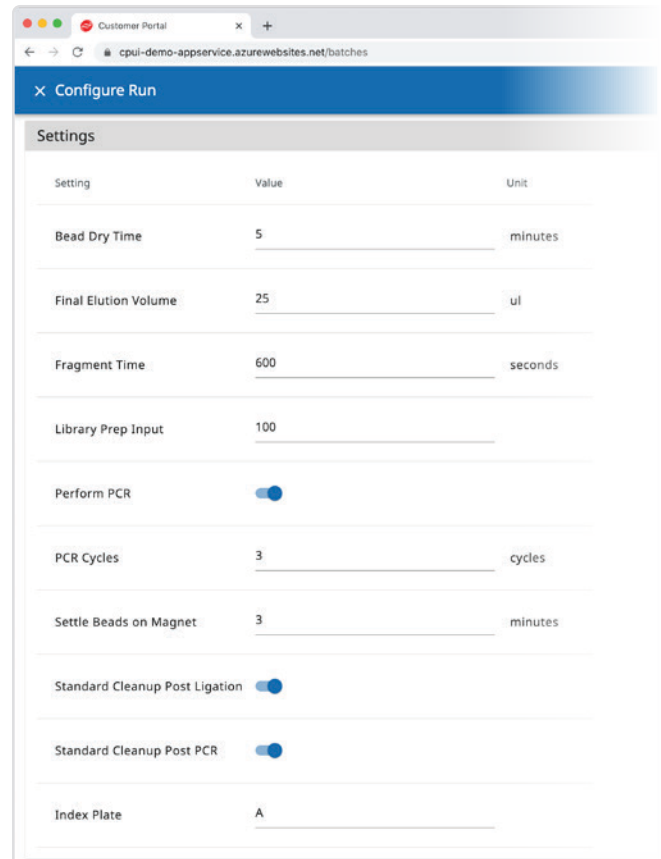
**Step 1** Select an application



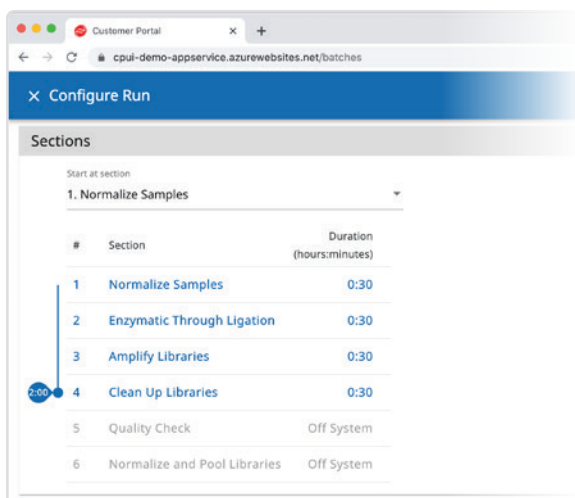
**Step 2** Enter batch name and number of samples from 4 to 24



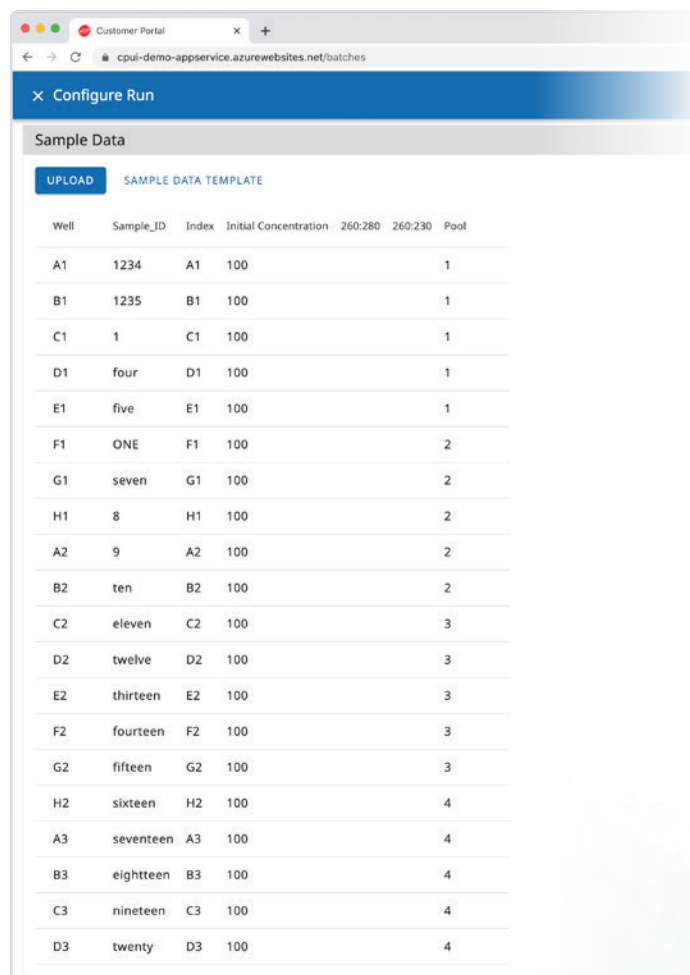
**Step 3** Make any necessary run adjustments



**Step 4** Select the stop point that fits your schedule

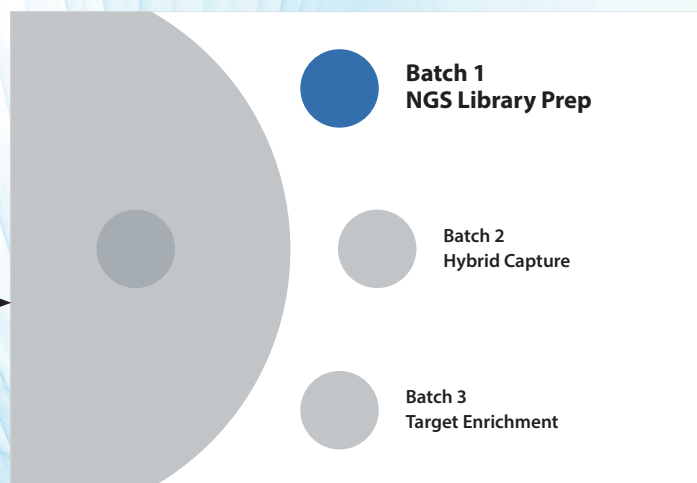
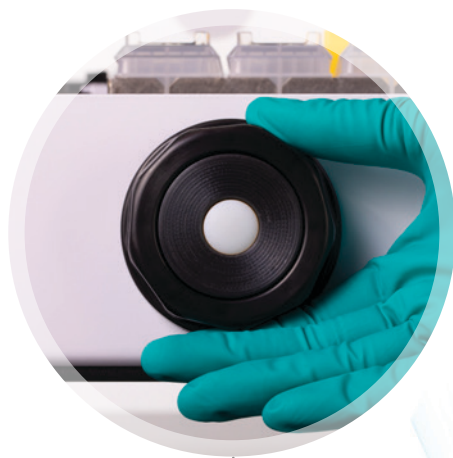


**Step 5** Upload your Sample Data and select 'Ready to Run'



On the instrument:

**Step 6** Use the dial on the instrument to select the batch and you are ready to run



# Biomek NGenius System

## Small footprint, unimaginable power and performance

Part Number	Description	Power Requirements	Weight	Dimensions (W x D X H)
C62703	Biomek NGenius System	10 A, 100-240 V, 50/60 Hz	135 lb (61.23 kg)	73.2 cm x 78.5 cm x 128.5 cm (with door open)

## Biomek NGenius Consumables

### The most critical point of contact

The Biomek NGenius System is optimized to perform as a complete system, which is why only Biomek consumables from Beckman Coulter Life Sciences are validated and approved for use.

Part Number	Description	Quantity
C62705	Biomek Reaction Vessel, 24 well	64/Box
C62706	Biomek Lid, 24 well	96/Box
C62707	Biomek Bulk Reservoir, 20 mL/section	32/Box
C59585	Case, Tips, 1025 µL, Conductive, Filtered, 480 Tips	5 Racks/Box (96 per rack)
C62712	Case, Tips, 70 µL, Conductive, Filtered, 3840 Tips	10 Racks/Box (384 per rack)
C70672	Case, Biomek 1025 µL Tip Box, Empty, 5 Racks	5 Racks/Box
C70673	Case, Biomek 70 µL Tip Box, Empty, 10 Racks	10 Racks/Box
C70665	Biomek Seal Plate, 24 well	20/Box

Part Number	Description
C81659	<b>Biomek NGenius Instrument 1 Year Service Agreement</b>
C81657	<b>Biomek NGenius Instrument IQ Service Agreement</b>
C81658	<b>Biomek NGenius Instrument OQ Service Agreement</b>
C89885	<b>Getting Started with Your Biomek NGenius System</b> This training is for administrators to be introduced to the hardware and consumables as well as understand instrument activation and account management. This training should take place prior to instrument installation if an existing administrator is not already trained and available.
C89886	<b>Learning to Use Your Biomek NGenius System</b> This training is for the administrator to better understand how App Setup and testing can be performed. It also demonstrates how Batches, Work Aids, and Batch Reports are used within the Biomek NGenius Customer Portal.
C89887	<b>Using Your Biomek NGenius System - for Administrators</b> This training is for operators to be introduced to the hardware and consumables used on the system as well as understanding how Batches, Work Aids, and Batch Reports are used within the Biomek NGenius Customer Portal.

Contains Mercury Dispose per State, Local, and National Laws  
For Canada customers: Visit [mybeckman.ca](http://mybeckman.ca) for disposal and recycling information for this product

Contient du mercure Éliminer selon les lois d'État, locales et nationales Pour les clients au Canada: Visitez [mybeckman.ca](http://mybeckman.ca) pour l'élimination et le recyclage de l'information pour ce produit



© 2022 Beckman Coulter, Inc. All rights reserved. Beckman Coulter, the stylized logo, and the Beckman Coulter product and service marks mentioned herein (including Biomek and Biomek NGenius) are trademarks or registered trademarks of Beckman Coulter, Inc. in the United States and other countries.

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

Each app template for use in the Biomek NGenius Next Generation Library Prep System has been demonstrated but has not been validated by Beckman Coulter. Beckman Coulter makes no warranties of any kind whatsoever express or implied, with respect to app templates, including but not limited to warranties of fitness for a particular purpose or merchantability or that app templates are non-infringing. All warranties are expressly disclaimed. Your use of any app template is solely at your own risk, without recourse to Beckman Coulter.

The app templates have been demonstrated for use on the Biomek NGenius system for the chemistry kit version and the release date shown at the time when the app is selected and created, but has not been validated by Beckman Coulter for use in the diagnosis of disease or other clinical conditions.

Biomek NGenius Next Generation Library Prep Systems and demonstrated applications are not intended or validated for use in the diagnosis of disease or other conditions.

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

20.12.476.GEN

