



Biomek **NGenius**

NEXT GENERATION LIBRARY PREP SYSTEM

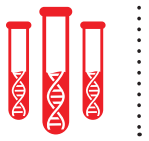
IT'S NOT JUST SMART. IT'S **INGENIOUS.**



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

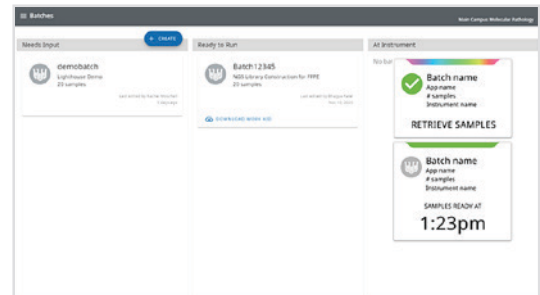
Control user access to applications with a simple authorization process (if required). Ensure that only authorized users have access to non-lab validated applications.



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



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.

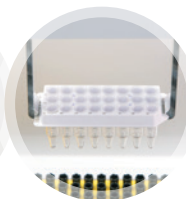


* Requires a Google Chrome or Microsoft Edge enabled computer.

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.

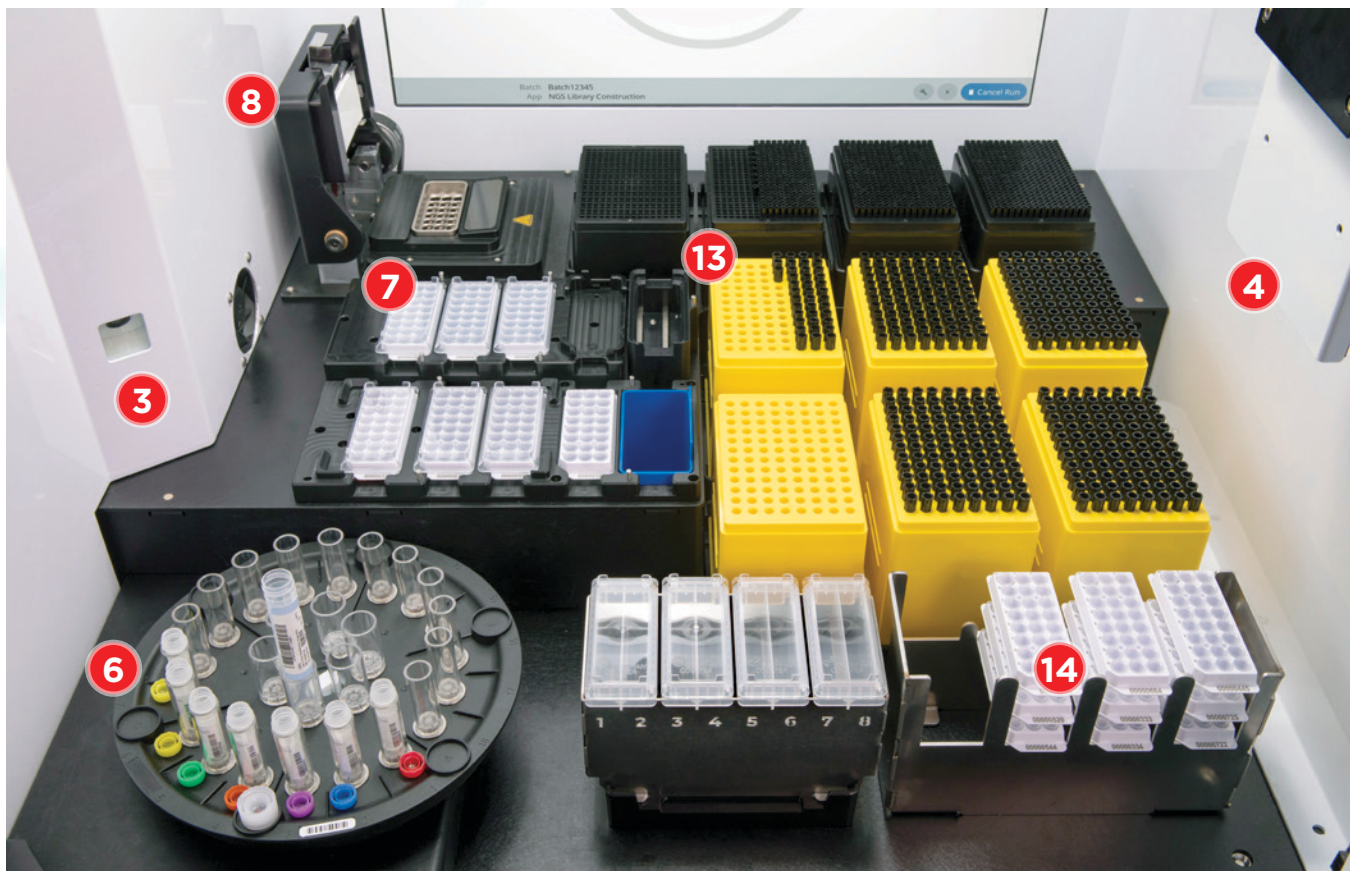


Error reduction

- 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

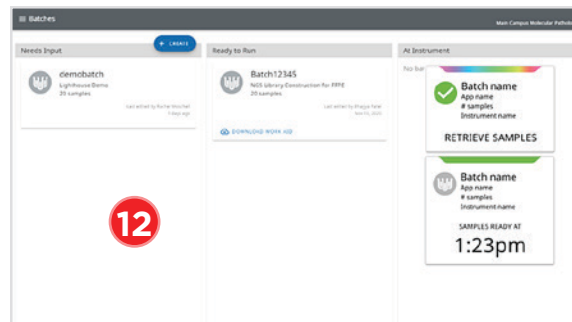
Less hands-on-time

- 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 μ 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.



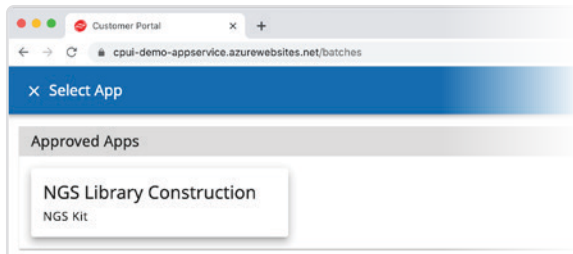
The **Genius** is in its **Simplicity**

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.

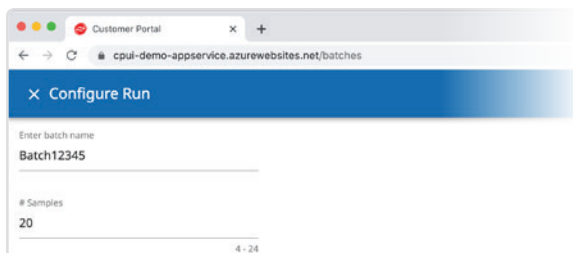
Set up and run a batch in 6 easy steps...

On the Biomek NGenius Portal software:

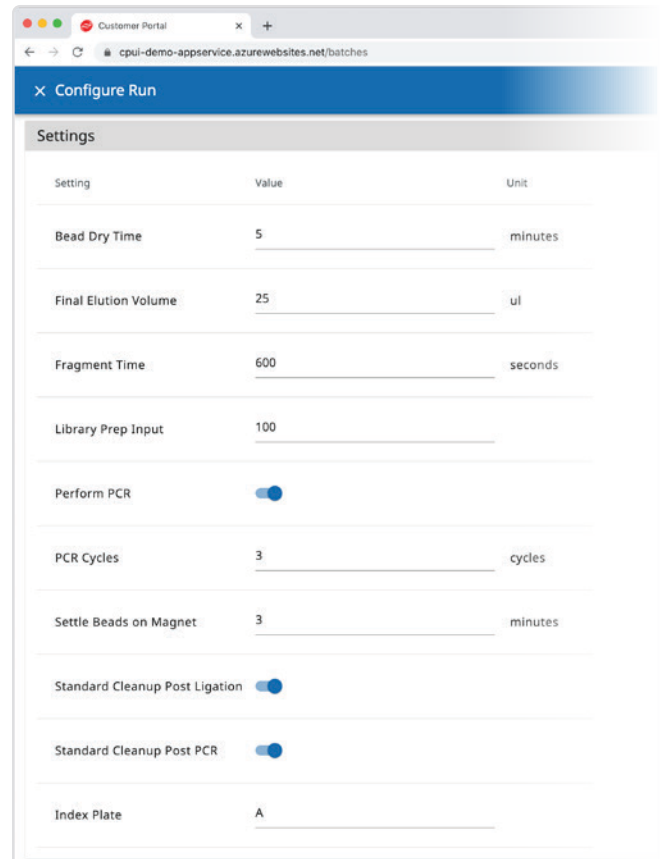
Step 1 Select an application



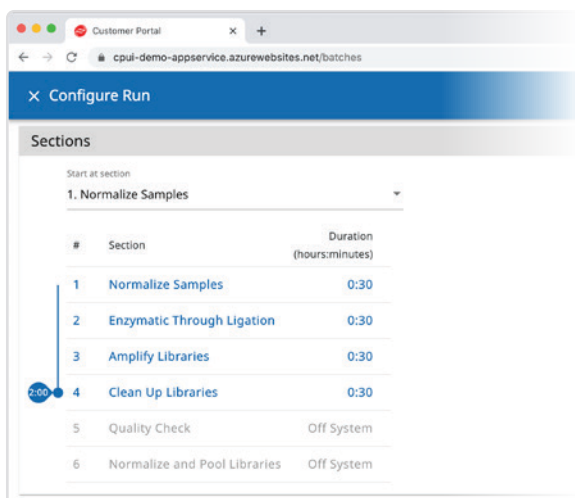
Step 2 Enter batch name and number of samples from 4 to 24



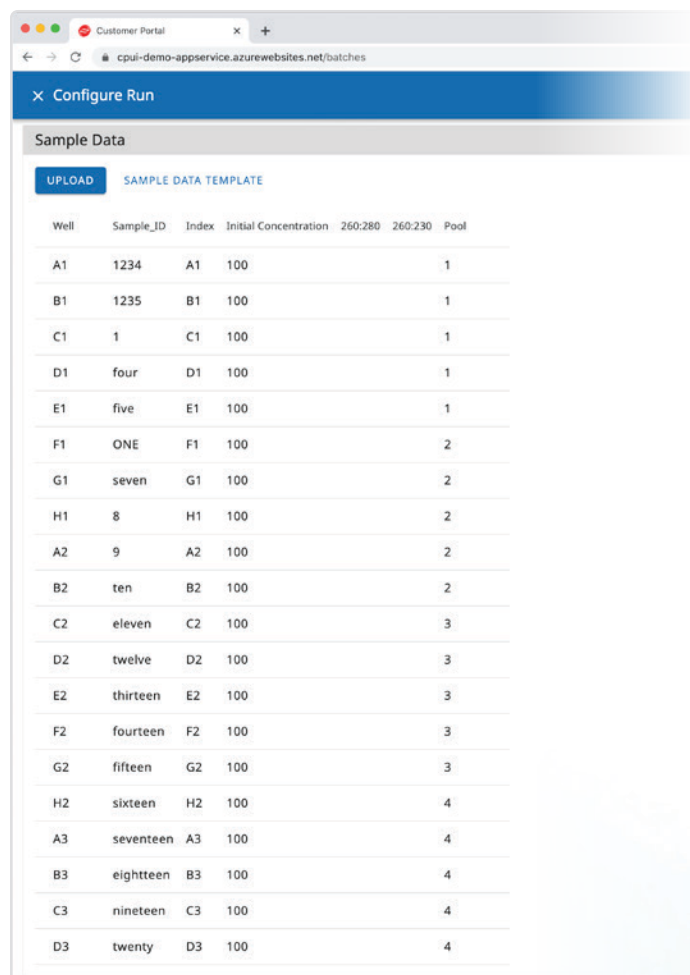
Step 3 Make any necessary run adjustments



Step 4 Select the stop point that fits your schedule

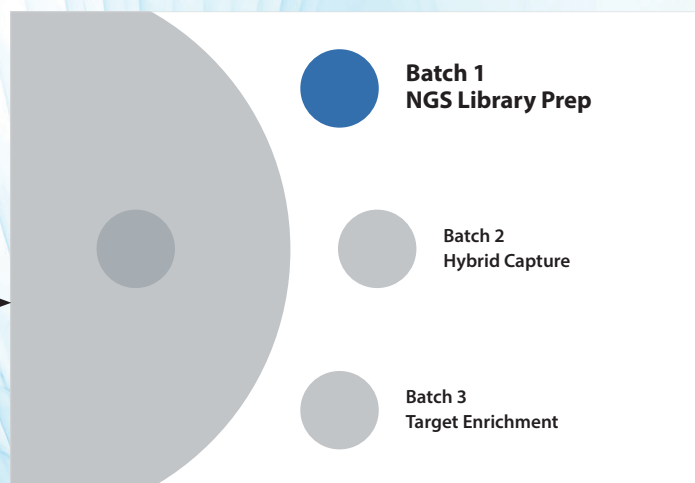


Step 5 Upload your Sample Data and select 'Ready to Run'



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	Biomek NGenius Instrument 1 Year Service Agreement
C81657	Biomek NGenius Instrument IQ Service Agreement
C81658	Biomek NGenius Instrument OQ Service Agreement
C89885	Getting Started with Your Biomek NGenius System 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	Learning to Use Your Biomek NGenius System 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	Using Your Biomek NGenius System - for Administrators 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 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 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

20.12.476.GEN

