



DISCOVERY
in motion.

LEARN VIA YOUR PREFERRED METHOD

WITH “MULTICHANNEL” TRAINING OPTIONS FOR BIOMEK I-SERIES WORKSTATIONS

Pick your preferred channel: in-person (our lab or yours), online or Virtual Reality (VR). The choice is yours when scheduling training for your Biomek instrument(s). Some questions to consider:



Are you interested in a state-of-the-art immersive technology such as Virtual Reality that lets you interact with an instructor and get virtual “hands-on” training on your Biomek instrument?

Our 2-day, instructor-led Virtual Reality training session (VR headset provided) lets you train from almost anywhere you can access the internet. To complement instructor presentations and workbook exercises, you'll be able to “physically” interact with a virtual Biomek workstation as if it were right in front of you.



Do you prefer to learn at your own speed? Are you already comfortable using videos and other digital content to teach yourself something new?

Our e-Learning modules were developed for new and intermediate Biomek users to walk through the basics of liquid handling with Biomek 5 software, and to learn best practices. These self-paced videos are separated into short 10-45 min modules to allow you to learn whenever it is best for you.



Do you prefer to learn by talking to someone in real time? Can you only dedicate two full days to learning?

Our instructor-led virtual classes are only two days long. You will learn the basics of Biomek 5 software while watching and listening to your instructor discuss concepts via a virtual meeting. Virtual training allows you to ask questions during class, and includes 60 days of email support for follow-up questions as you work through assigned hands-on training with your instrument.



Do you prefer in-person instruction and can travel? Are you not quite comfortable operating your instrument on your own?

We provide in-person training at our facility in Indianapolis. This 3-day course starts with new concepts in software simulation before moving to hands-on instrument training to build your confidence using Biomek automation.



Still prefer in-person training but not able to travel to Indianapolis?

We provide customer site training where we come to you! You will start by learning new concepts in simulation before we move to learning on your own instrument.

Ready to take your Biomek to the next level?

Talk to your local Sales Representative or email us: AutomationLearningCenter@beckman.com



TRAINING CLASS DESCRIPTIONS



Online & Virtual Learning Options

Self-paced Biomek i-Series Method Programming Training for Beginners

- Multichannel and Span-8 Pods (C73060)
- Span-8 Pod (C73061)
- Multichannel Pod (C73062)

Available in our Learning Center, these self-paced e-Learning modules will introduce the software and cover the use and configuration of all the Liquid Handling Steps as well as the Project Editors and how to use them during method programming. Available as Multichannel & Span-8 Curriculum, Multichannel Only Curriculum, and Span-8 Only Curriculum.

Instructor-led online Biomek i-Series Method Programming Training for Beginners (C73056)

This 2-day class taken from your home or office covers material using a mix of presentation and workbook exercises and allows for hands-on exercises to be performed after training is complete on your instrument. Method development skills will gradually build as we learn how to configure and use the Liquid Handling Steps tab as well as how to use the Project Editors. 60 days of email support are provided to guide you through the hands-on exercises.

Instructor-led Virtual Reality (VR) for Beginners (C73056)

During this virtual 2-day training session—which you can take almost anywhere with internet access—we'll use a mix of presentation and workbook exercises, and you'll be able to interact with the instrument and instructor as if they were right in front of you.



In-Person Options

Biomek i-Series Method Programming Training for Beginners at Beckman Indianapolis Site (B87510)

This 3-day class covers material using a mix of presentation, workbook exercises, and hands-on exercises to help you gradually learn how to configure and use the Liquid Handling Steps tab as well as how to use the Project Editors. The last half of day 3 includes an opportunity to discuss topics such as file management, revision history, and common troubleshooting techniques, as well as a hands-on opportunity to perform framing of the instrument.

Biomek i-Series Method Programming Training for Beginners at Customer Site (C02855)

This course is presented in your own lab and covers the same content that would be given at the Indianapolis site or in a virtual class. As time permits, device integrations on your system may be included.

Biomek i-Series Software Training for Experienced Users at Beckman Indianapolis Site (B86572)

This Biomek software training is intended for those who have previous Biomek experience and have been recommended for the class. The course will cover advanced method programming topics such as: Selection as Text, Expressions, Arrays, Global Variables, Run Program, Procedures, If/Then, Worklists and Data Sets.

Contact us at AutomationLearningCenter@beckman.com to learn more.



© 2023 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.

For Beckman Coulter's worldwide office locations and phone numbers, please visit Contact Us at [beckman.com](https://www.beckman.com)
2023-AMER-EN-100952-V1