



Biomek **NGenius** Cloud Services

SECURITY AND PRIVACY FAQs



An unwavering
commitment to
safety and
security.

ACCELERATING
answers.

 **BECKMAN
COULTER**
Life Sciences



Beckman **NGeniusS** Cloud Services

At Beckman Coulter, we develop technologies to advance and optimize the laboratory. To achieve this, we uphold core values that define our responsibility to those we serve. Among them: an unwavering commitment to safety and security. Therefore, we believe in continuously improving to address the ever-evolving privacy and cybersecurity landscape.

We have implemented privacy and security by design in our development processes. The source of these processes and protections are derived from multiple regulatory frameworks (i.e. FDA, HIPAA, GDPR) and internationally recognized cybersecurity best practices (i.e. ISO 27001), and are based upon the 'CIA' triad of Confidentiality, Integrity and Availability.

Security and Privacy FAQs

Q: What is the general architecture of the Beckman NGeniusS Cloud Services?

A: Beckman Coulter Life Sciences uses a Software as a Service (SaaS) model to deliver Beckman NGeniusS Cloud Services. This means that the infrastructure components are shared across the solution, but your lab data is segmented and protected so that only you, and the team members that you set up, have access to your lab data. In addition, Beckman Coulter has dedicated, standalone cloud platforms across the globe to serve different geographies. For example, a separate cloud platform is hosted in the European Union to serve EU-based countries.

Q: How does Beckman Coulter Life Sciences secure connections to the Beckman NGeniusS Cloud Services?

A: We secure web browser connections to the Beckman NGeniusS Cloud Services with HTTPS and end-to-end encryption via Transport Layer Security (TLS) 1.2. In addition, a Web Application Firewall using Open Web Application Security Project (OWASP) rule sets is configured to prevent attacks against the Beckman NGeniusS Cloud Services. Data

transfer between Beckman NGeniusS instruments and Beckman NGeniusS Cloud Services utilizes certificate-based mutual SSL authentication and is encrypted end-to-end via Transport Layer Security (TLS) 1.2.

Q: Can the Beckman NGeniusS Next Generation Library Prep System be installed on a segmented network?

A: Yes. The Beckman NGeniusS Next Generation Library Prep System can be installed on a network segment that is separate from your enterprise systems. Further details of the connectivity requirements can be found in the product documentation.

Q: Does the Beckman NGeniusS Next Generation Library Prep System require inbound connections to my corporate network?

A: No. Communication is initiated by the Beckman NGeniusS Next Generation Library Prep System to the Beckman NGeniusS Cloud Services. An inbound connection to your network is not required. Further details of the connectivity requirements can be found in the product documentation.



Q: Does Biomek NGenius Cloud Services enforce strong passwords?

A: Yes. Biomek NGenius Cloud Services enforces strong passwords, including complexity and length requirements. Further details can be found in the product documentation.

Q: Can I install my own security software on the Biomek NGenius Next Generation Library Prep System?

A: No. The Biomek NGenius Next Generation Library Prep System is locked down with application allow-listing software to ensure that only pre-approved programs can be run. This security control prevents the installation of any additional software and cannot be disabled.

Q: How does Beckman Coulter Life Sciences manage product security for the Biomek NGenius instrument and Biomek NGenius Cloud Services?

A: We have a dedicated product security team that is involved in product development from concept through production release. Reviews are conducted by the product security team at key checkpoints in the development process to ensure security standards are met or exceeded. In addition, we contract with third-party penetration testing firms to get an objective view of our system performance from the security perspective.

Q: Does Beckman Coulter Life Sciences have an information security policy?

A: Yes. We have a security policy that governs the implementation of information security to ensure the confidentiality, integrity, and availability of data in the care of Beckman Coulter Life Sciences.

Q: What areas does Beckman Coulter's information security policy cover?

A: We consider security to be an essential foundation for our products and services. The information security policy covers areas such as access control and account management, supplier security, physical security, people security, system development, security education, incident management and response, and business continuity.

Q: How does the Biomek NGenius system protect customer data?

A: Biomek NGenius Next Generation Library Prep System and Cloud Services protect your data by:

- Encrypting data at rest and in transit
- Ensuring high availability with fault-tolerant architecture
- Ensuring business continuity by storing encrypted data backups
- Managing security risks with a formal Information Security Management System (ISMS)
- Conducting internal and external penetration tests of our systems

Q: Does Biomek NGenius have a privacy policy?

A: Yes. Biomek NGenius has a privacy policy that covers the types of information collected, the usage of this information, and under what circumstances the information is shared. The privacy policy is presented to each user of the system when they are invited to use the Cloud Services and when the privacy policy is updated.

Continued on next page



Security and Privacy FAQs (continued)

Q: Does Beckman Coulter Life Sciences have an ISO 27001 certification?

A: Yes. The Beckman Coulter Life Sciences Information Security Management System (ISMS) is certified for the ISO 27001 standard. The Biomek NGenius Next Generation Library Prep System and Biomek NGenius Cloud Services are in scope for the ISO 27001 certification and are developed and operated in accordance with the ISMS. In addition, we annually conduct both internal and external audits ensure continued compliance with the ISO 27001 standard. Certification details are available by contacting your sales representative.

Q: Is Beckman Coulter Life Sciences GDPR compliant?

A: Biomek NGenius is built to support customer compliance with GDPR. For our EU customers, the Biomek NGenius Cloud Services are hosted in the EU.