

Ion Chef System

Simple, reproducible, automated workflows for Ion Torrent sequencing

The Ion Chef™ System simplifies the workflow for the Ion GeneStudio™ S5 series, Ion Proton™, and Ion PGM™ systems, providing a walk-away, reproducible solution for automated Ion AmpliSeq™ library preparation and automated template preparation and chip loading.

Ion Chef System

- Automates library generation, equalization, and pooling for Ion AmpliSeq™ 1- and 2-pool designs for up to 8 samples in a single instrument run
- Simplifies template preparation and chip loading using a single automated process for loading of up to two Ion Torrent™ semiconductor sequencing chips per instrument run
- Helps reduce sources of variability for users at all experience levels
- Helps save time and labor with just 15 minutes of hands-on time for setup prior to start of automated workflows
- Supports the Ion GeneStudio S5 series, Ion PGM, and Ion Proton systems, and all Ion Torrent semiconductor chips and sequencing chemistries

Simple, automated workflow

The Ion Chef System simplifies the Ion Torrent semiconductor sequencing workflow by integrating several processes into a single instrument for maximum efficiency and throughput (Figure 1). The entire process is automated, whether starting from DNA samples and obtaining equalized Ion AmpliSeq libraries or starting from prepared libraries and obtaining a final output of up to two sequencing-ready chips. The instrument has an intuitive touch-screen interface, making setup easy for users at any experience level.



Reduced variability

The Ion Chef System helps reduce sources of variability for users at any experience level by greatly minimizing pipetting errors, providing reproducible chip loading, and enabling sample tracking and system checks with the onboard tracking system.

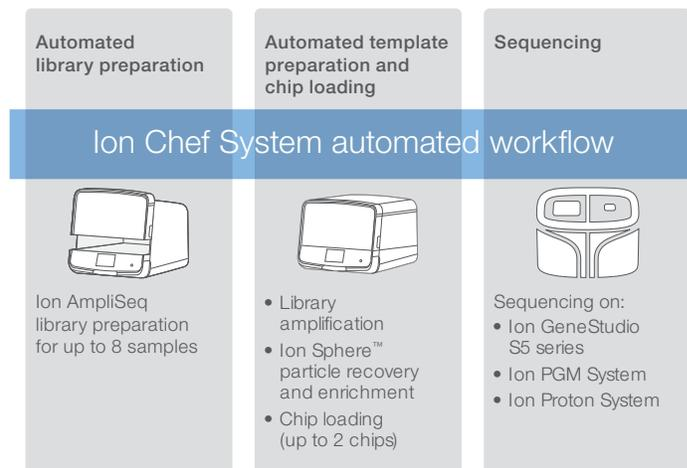


Figure 1. The Ion Chef System automated workflow includes less than 45 minutes of hands-on time. The Ion Chef System fully automates library preparation and template preparation and produces sequencing-ready Ion Torrent semiconductor chips.

Equipped with barcode-reading capabilities, the tracking system enables identification of samples, onboard reagents, and chips, and integrates with Torrent Suite™ Software and/or a LIMS system. For added assurance, the tracking system also performs setup checks, including scanning the instrument deck for positioning of reagents and consumables, calibration checks, and instrument diagnostics. Combined with the simple and robust Ion Chef System workflow, the tracking system helps reduce user-introduced errors and variability for maximum accuracy and efficiency.

Time and labor savings

The Ion Chef System enables time and labor savings, with just 15 minutes of hands-on time for setup. Simply select the run parameters using the intuitive touch-screen interface and load the library samples, reagents, consumables, and chips onto the deck. A single run of the Ion Chef System will produce one pooled Ion AmpliSeq library for up to 8 samples, or up to two loaded chips for sequencing on the Ion GeneStudio S5 series, Ion PGM System, or Ion Proton System, in a single-day workflow. The minimal hands-on time and walk-away automation of the Ion Chef System enables overnight runs, helping to maximize the productivity of your lab.

Support for Ion GeneStudio S5 series, Ion PGM, and Ion Proton systems

The Ion Chef System supports all Ion Torrent semiconductor sequencing chips and provides flexibility for most project types and sizes.

Ion Chef System specifications	
Platform compatibility	Ion GeneStudio S5 series Ion PGM System Ion Proton System
Chip compatibility	Ion 510™ Chip (200, 400 bases) Ion 520™ Chip (200, 400, 600 bases) Ion 530™ Chip (200, 400, 600 bases) Ion 540™ Chip (200 bases) Ion 550™ Chip (200 bases) Ion 314™ Chip (200, 400 bases) Ion 316™ Chip (200, 400 bases) Ion 318™ Chip (200, 400 bases) Ion PI™ Chip (200 bases)
Ion AmpliSeq panel compatibility	Supports all Ion AmpliSeq 1- and 2-pool panel designs. Reverse transcription for Ion AmpliSeq RNA panels must be done prior to loading cDNA on the Ion Chef instrument.
Library compatibility	Supports template preparation for a broad range of DNA and RNA libraries, including Ion AmpliSeq libraries
Setup time	15 minutes
Weight	150 lb
Dimensions	22.1 x 27.6 x 28.1 in (56.1 x 70 x 71.4 cm)
Power requirements	Voltage: 100 V (min) to 240 V (max) Current: 14 A (max) Frequency: 50/60 Hz Power draw: 1,350 W
Operating environment	Temperature: 20–25°C Humidity: 40–60% (noncondensing)

Ordering information

Product	Cat. No.
Ion Chef System	4484177
Ion AmpliSeq Kit for Chef DL8	A29024
Ion 510 & Ion 520 & Ion 530 Kit – Chef	A34019
Ion 540 Kit-Chef	A30011
Ion 550 Kit-Chef	A34541
Ion 510 Chip Kit	A34292
Ion 520 Chip Kit	A27762
Ion 530 Chip Kit	A27764
Ion 540 Chip Kit	A27766
Ion 550 Chip Kit	A34538
Ion PGM Hi-Q View Chef Kit	A29902
Ion 314 Chip Kit v2 BC	4488144
Ion 316 Chip Kit v2 BC	4488149
Ion 318 Chip Kit v2 BC	4488150
Ion PI Hi-Q Chef Kit	A27198
Ion PI Chip v3, 8-pack	A26771

Find out more at thermofisher.com/ionchef