SPECIFICATION SHEET Ion PGM System

## The Ion PGM System

High-quality, accessible sequencing

The Ion PGM™ System makes affordable, high-quality next-generation sequencing accessible to human disease researchers and microbiologists around the world. The Ion PGM System is a reliable sequencing platform that combines simple sample preparation and data analysis solutions with flexible chip output for ultimate project flexibility.



| Ion PGM System performance specifications |                                   |  |   |   |  |  |  |
|---|-----------------------------------|--|---|---|--|--|--|
|   |                                   | lon 314™ Chip v2 or<br>Ion 314 Chip v2 BC  | lon 316™ Chip v2 or<br>Ion 316 Chip v2 BC | Ion 318™ Chip v2 or<br>Ion 318 Chip v2 BC |  |  |  |
| Output*                                   | 200 base<br>400 base <sup>†</sup> | 30-50 Mb<br>60-100 Mb  | 300-600 Mb<br>600 Mb-1Gb                  | 600 Mb-1Gb<br>1.2-2 Gb                    |  |  |  |
| Reads                                     |                                   | 400-550 thousand   | 2–3 million                               | 4–5.5 million                             |  |  |  |
| Run time                                  | 200 base<br>400 base              | 2.3 hr<br>3.7 hr   | 3.0 hr<br>4.9 hr                          | 4.4 hr<br>7.3 hr                          |  |  |  |
| Research areas                            |                                   | Cancer research, inherited disease research, microbial genomics, stem cell research, agriculture, epigenomics, metagenomics, forensic science, and ancient DNA genomics  |   |   |  |  |  |
| Key applications                          |                                   | Targeted DNA sequencing, copy number analysis, targeted RNA sequencing, small RNA sequencing, de novo microbial sequencing, bacterial typing research, viral typing research, ChIP sequencing, methylation analysis, SNP verification, and genotyping by sequencing                |   |   |  |  |  |
| Target selection solutions                |                                   | Ion AmpliSeq <sup>™</sup> technology, Ion TargetSeq <sup>™</sup> kits (3 custom enrichment kits and 1 multiblocker kit)  |   |   |  |  |  |
| Library solutions                         |                                   | lon AmpliSeq™ Library Kits, Ion Xpress™ Plus Fragment Library Kit, Ion Total RNA-Seq Kit v2, Ion Library Equalizer™ Kit, and 384 barcodes supported by Torrent Suite™ Software   |   |   |  |  |  |
| Data analysis solutions                   |                                   | Torrent Suite Software processes signals, calls bases, aligns sequences, and performs basic variant calling as the primary analysis software for the lon PGM™ Sequencer.   |   |   |  |  |  |
|   |                                   | Optimized data analysis workflows have been developed for targeted DNA sequencing, targeted RNA sequencing, microbial <i>de novo</i> sequencing, microbial typing research, and other research applications using third-party software solutions and freely downloadable plug-ins. |   |   |  |  |  |
|   |                                   | Ion Reporter™ Software performs controlled analysis, annotation, and reporting of variants along with application-specific workflows to support single, trio, and paired tumor–normal analysis.  |   |   |  |  |  |



## **ion**torrent

| Ion PGM Sequencer specifications             |  |  |  |  |  |  |
|--|--|--|--|--|--|--|
| Working environment<br>(for indoor use only) | Temperature:<br>68-77°F (20-25°C)<br>Humidity: 40-60%,<br>noncondensing<br>Altitude: <6,500 ft (2,000 m) | Clearances: 12 in. (30.5 cm) in rear 4 in. (10 cm) on left side 4 in. (10 cm) on right side 4 in. (10 cm) from front edge of bench to sequencer beze 36 in. (90 cm) aisle in front of bench for operator access 8 in. (20 cm) from front edge of bench to the conical tube |  |  |  |  |
| Gas supply                                   | Connection: 0.25 in. push-to-connect fitting   | Pressure: 35-45 psi  | Composition: nitrogen (grade 4.5, 99.995% or better) |  |  |  |
| Other connections                            | Ethernet: 1 GigE   | USB: 2x USB 2.0  |  |  |  |  |
| Power  | Voltage: 100 V (min) to<br>240 V (max)<br>Current: 9 A (max)   | Frequency: 50/60 Hz<br>Power draw: 200–300 W   |  |  |  |  |
| Dimensions                                   | 24 x 20 x 21 in./61 x 51 x 53 cm (W x D x H)   |  |  |  |  |  |
| Weight                                       | Crated for shipment:<br>85 lb/39 kg  | Free-standing: 65 lb/30kg  | 1  |  |  |  |

| Ion PGM Torrent Server specifications                                  |  |  |  |  |  |  |
|--|--|--|--|--|--|--|
| Product configuration  | A single free-standing tower computer appliance, included with the purchase of the Ion PGM System. Includes Torrent Suite Software with all necessary software components to deliver signal processing, base calling, read alignment, and variant calling. |  |  |  |  |  |
| Processor  | Dual 8-core 2.6 GHz CPUs   |  |  |  |  |  |
| Memory   | 64 GB RAM  |  |  |  |  |  |
| Storage (approx.)  | 18 TB  |  |  |  |  |  |
| Operating System   | Ubuntu™  |  |  |  |  |  |
| Dimensions (approx.)   | 9 x 28 x 43 in./22 x 71 x 17 cm (W x D x H)  |  |  |  |  |  |
| Weight (approx.)   | 121 lb/55 kg   |  |  |  |  |  |
| Power  | Voltage: 100 V (min) to 240 V (max)<br>Current: 12 A (max)   | Frequency: 50/60 Hz<br>Power draw: 1,100 W |  |  |  |  |
| Data formats Industry-standard FASTQ, SFF, BAM, and VCF format outputs |  | VCF format outputs                         |  |  |  |  |

## **Ordering information**

| Product  | Cat. No. |
|--|----------|
| Ion PGM System (includes Ion PGM Sequencer and Ion PGM Torrent Server) | 4462921  |
| Ion Chef System  | 4484177  |
| Ion OneTouch 2 System  | 4474779  |

## Find out more at thermofisher.com/pgm



 $<sup>\</sup>dagger$  400 base pair chemistry is not compatible with the Ion Chef  $^{\!\scriptscriptstyle{\text{TM}}}$  Instrument at this time.

