

Ion PGM™ Checklist — Sequencing

Pub. no. MAN0009127 Rev. 1.0

Follow the steps of this checklist if you are an experienced user of the Ion PGM™ Sequencing 200 Kit v2 (Catalog no. 4482006) with the Ion Personal Genome Machine® (PGM™) System.

For complete kit and protocol information, including precautions and safety information, refer to the *Ion PGM™ Sequencing 200 Kit v2 User Guide* (Pub. no. MAN0007273), available on the **Ion Community** website.

Before starting

Thaw Sequencing Primer on ice.

Add Control Ion Sphere Particles (IPs) to the enriched, template-positive ISPs

1. Ensure that enriched, template-positive ISPs (prepared using the Ion PGM™ Template OT2 200 Kit) are in a 0.2-mL non-polystyrene PCR tube.
2. Vortex Control ISPs and centrifuge for 2 seconds.
3. Add 5 µL of Control ISPs to the appropriate volume of enriched, template-positive ISPs for your chip type:

Ion 316™ Chip v2/Ion 318™ Chip v2	Ion 314™ Chip v2
Entire volume of prepared ISPs	Half the volume of prepared ISPs (transfer and store half of the volume)

Anneal the Sequencing Primer

1. Mix the tube containing ISPs and controls by thoroughly pipetting up and down. Centrifuge for 2 minutes at 15,500 × g.
2. Carefully remove the supernatant, leaving the following volume in the tube:

Volume	Ion 316™ Chip v2/Ion 318™ Chip v2	Ion 314™ Chip v2
Remaining in tube	15 µL	3 µL

3. Add Sequencing Primer and confirm the total volume (adjust the volume with Annealing Buffer, if necessary):

Volume	Ion 316™ Chip v2/Ion 318™ Chip v2	Ion 314™ Chip v2
Sequencing Primer	12 µL	3 µL
Total volume	27 µL	6 µL

4. Pipet the sample up and down thoroughly to disrupt the pellet.
5. Program a thermal cycler (95°C for 2 minutes, then 37°C for 2 minutes), insert tube in cycler, and run the program.

Chip Check

1. Remove a new chip from its package and label it. Press **Run** on the Ion PGM™ Sequencer touchscreen.
2. When prompted, ground yourself. Replace the old chip in the chip clamp with the new chip. (*Do not wear gloves.*)
3. When prompted, scan the new chip package barcode or press **Change** to manually enter barcode. Press **Chip Check**.
4. Check for leaks on the chip.
- If there is a leak, press **Abort** and proceed to Troubleshooting in the User Guide.
 - If the chip passes, press **Next**.
 - If the chip fails, re-seat the chip, and press **Calibrate**. If the chip still fails, restart with a new chip.
5. Remove the new chip, re-insert the old chip in the socket, close the clamp, and empty the waste bottle.

Bind Sequencing Polymerase to the ISPs

1. Remove ISPs from the thermal cycler and add polymerase according to the following table:

Volume	Ion 316™ Chip v2/Ion 318™ Chip v2	Ion 314™ Chip v2
Ion PGM™ Sequencing 200 v2 Polymerase	3 µL	1 µL
Total volume	30 µL	7 µL

2. Pipet the ISPs up and down, and incubate at room temperature for 5 minutes.

Load the chip

1. Tilt the new chip 45 degrees. Insert pipette tip into the loading port; remove and discard as much liquid as possible.
2. Place the chip *upside-down* in the Ion MiniFuge bucket; transfer the bucket to the MiniFuge (*chip tab pointing in*).
3. Centrifuge for 5 seconds. Remove the chip from the bucket and wipe off any liquid on the bucket.
4. Put the chip back in the bucket and place the bucket on a flat, stable surface.
5. Following polymerase incubation, collect the entire sample into the pipette tip and insert the tip into chip loading port.
6. *Dial down the pipette* to gently and slowly deposit the ISPs, leaving ~0.5 µL of sample in the pipette tip.
7. Remove and discard any displaced liquid from the other port of the chip.
8. Transfer the chip in the bucket to the MiniFuge (*chip tab in*). Centrifuge for 30 seconds.
9. Set the pipette volume according to the following table:

Volume	Ion 316™ Chip v2/Ion 318™ Chip v2	Ion 314™ Chip v2
Pipette setting	30 µL	5 µL

10. Tilt the chip 45 degrees, and slowly pipet the sample in and out of the chip *three times*.
11. Centrifuge the chip in the Minifuge (*with the chip tab out*) for 30 seconds.
12. Repeat step 10 one more time, then centrifuge for 30 seconds (*chip tab in*).
13. Tilt the chip and slowly remove liquid by dialing the pipette.
14. Discard the liquid. If some liquid remains in the chip, perform a 5-second quick spin (*chip tab out*).
15. If necessary, lightly tap the point of the chip tab against the benchtop, and remove and discard any collected liquid.

Select the Planned Run and perform the run

1. Press **Browse** on the sequencer touchscreen and select the name of the plan you created. Press **Next**.
2. Confirm that the run settings are correct or make any necessary changes. Press **Next**.
3. Load and clamp the chip, then press **Next**.
4. Visually inspect the chip in the clamp for leaks before closing the cover.
- If the chip passes, press **Next**.
 - If it fails, press **Abort**, re-seat chip, and press **Calibrate**.
 - If the chip continues to fail calibration, contact Technical Support.

For Research Use Only. Not for use in diagnostic procedures.

Limited Product Warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at www.lifetechnologies.com/termsandconditions. If you have any questions, please contact Life Technologies at www.lifetechnologies.com/support.

© 2013 Life Technologies Corporation. All rights reserved. The trademarks mentioned herein are the property of Life Technologies Corporation and/or its affiliate(s) or their respective owners.

DISCLAIMER

LIFE TECHNOLOGIES CORPORATION AND/OR ITS AFFILIATE(S) DISCLAIM ALL WARRANTIES WITH RESPECT TO THIS DOCUMENT, EXPRESSED OR IMPLIED, INCLUDING BUT NOT LIMITED TO THOSE OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, OR NON-INFRINGEMENT. TO THE EXTENT ALLOWED BY LAW, IN NO EVENT SHALL LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) BE LIABLE, WHETHER IN CONTRACT, TORT, WARRANTY, OR UNDER ANY STATUTE OR ON ANY OTHER BASIS FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING BUT NOT LIMITED TO THE USE THEREOF.

Headquarters

5791 Van Allen Way | Carlsbad, CA 92008 USA | Phone +1 760 603 7200 | Toll Free in USA 800 955 6288

For support visit iontorrent.com/support or email ionsupport@lifetech.com

lifetechnologies.com

4 September 2013

