

Quick Start Guide

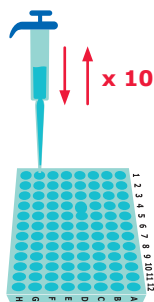
5R-PLEX Kit

MBD6000

Prepare PCR 1 Mix

Water, microbial DNA-free	_____ μL
5X HF Buffer	10 μL
dNTPs	1 μL
5R-PLEX PCR1 Primers mix	0.25 μL
HF DNA Polymerase (add last)	0.5 μL
DNA (Add separately)	_____ μL
Total PCR1 volume	50 μL

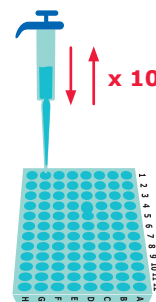
1. Dispense PCR1 mix in each well.
 2. Add _____ μL DNA.
 3. Mix up and down 10 times.
 4. Add 2 μL of diluted 5R-PLEX Positive Control.
 5. Add 6 negative controls (no DNA template).
- Total reaction volume/well 50 μL .



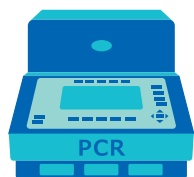
Prepare PCR 2 Mix

Water, microbial DNA-free	_____ μL
5X HF Buffer	10 μL
dNTPs	1 μL
5R-PLEX PCR2 For Primers mix	0.25 μL
HF DNA Polymerase (add last)	0.5 μL
PCR1 product (Add separately)	2–5 μL
Total PCR2 volume	50 μL

1. Dispense PCR2 mix in each well of the 5R-PLEX index plate.
 2. Add 2–5 μL of PCR1 Mix.
 3. Mix up and down ten times.
- Total reaction volume/well 50 μL .



PCR1 Program



98 °C	2 minutes	x 30
98 °C	10 seconds	
62 °C	15 seconds	
72 °C	35 seconds	
72 °C	5 minutes	

PCR2 Program



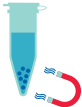
98 °C	2 minutes	x 6
98 °C	10 seconds	
64 °C	15 seconds	
72 °C	25 seconds	
72 °C	5 minutes	

Bead Purification Process

The Bead Purification Process is performed twice, the first time is post PCR2, and the second in the Dual Pooled Library Purification Step. Note step 2 below.

1. Vortex beads 30 seconds

2. Add 42.5 μL beads to 50 μL DNA sample (PCR2 product **OR** the purified Pooled Library)

3. Mix up and down ten times.

4. Place tubes on magnetic stand for 5 minutes.

5. Remove and discard the supernatant.

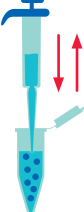
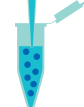
6. Add 200 μL of freshly prepared 80% ethanol.

7. Incubate the tubes on the magnetic stand for 30 seconds

8. Remove and discard the supernatant

Repeat steps 6-8 x 2
9. **IMPORTANT**
After the second wash, remove all traces of the ethanol wash.
 - a. Briefly centrifuge.

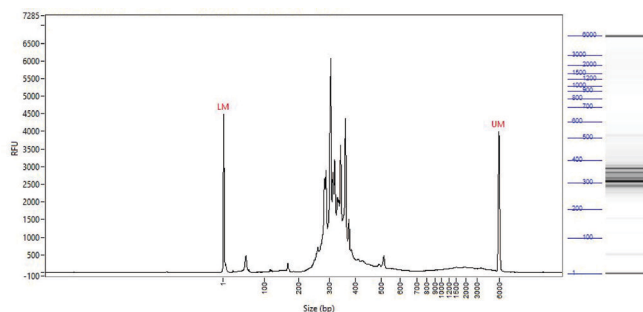
 - b. Return tubes to the magnetic stand and remove the ethanol residual with 10 μL pipette.

 - c. Air-dry for 3–5 minutes.
Do not over-dry the bead.
10. Remove tubes from the magnetic stand and add 20 μL of EB to each tube.

11. Mix up and down 10 times.

12. Incubate at room temperature 2 minutes
13. Place tubes on magnetic stand for 2–5 minutes or until supernatant is clear.

14. Transfer 18 μL of the supernatant (DNA) into a new tube.

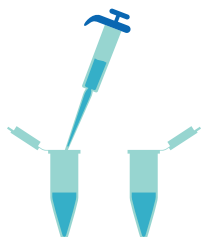
Sample Library QC

Run on a Multiplexed Capillary Electrophoresis (CE) such as Bioanalyzer, TapeStation, or equivalent. and verify amplicons traces in the range size of 250–400 bp.



Quantify and Pool

1. Measure the DNA concentration of the purified libraries with a fluorometric quantification method that uses dsDNA binding dyes, such as Qubit® dsDNA HS Assay Kit or equivalent.
2. Pool equal amount of each sample into a single tube (use the maximum amount possible).
3. Transfer a volume that is equivalent to up to 10 µg DNA of the pooled library in a new tube and continue to library purification.



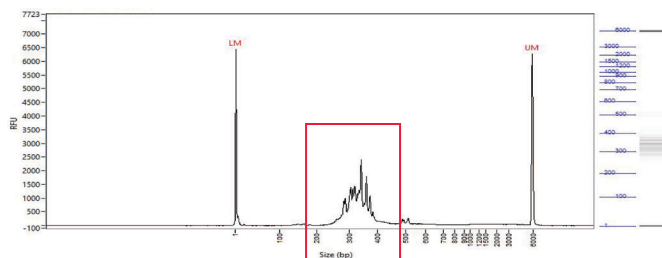
Dual Pooled Library Purification

1. Perform PCR clean up using the GenElute™ PCR Purification Kit (NA1020) or equivalent. Follow the GenElute™ PCR Clean-Up Procedure video at [SigmaAldrich.com/genelute-pcr-cleanup](https://www.sigmaaldrich.com/genelute-pcr-cleanup).
2. Repeat **Bead Purification Process** on the previous page.

Pooled Library QC and Quantification

1. Perform QC for library size: Run on a Multiplexed Capillary Electrophoresis instrument and verify amplicons traces in the range size of 250–400 bp.

IMPORTANT: If primer dimers are observed (~160–180 bp), it is highly recommended to repeat beads purification.



2. Library quantification: Measure the DNA concentration of the purified pooled library using KAPA Library Quantification Kit Illumina® Platforms (Roche, KK4873) (recommended) or Qubit® dsDNA HS Assay Kit.

Prepare the Library for NGS

Follow the Illumina® guide for preparing the library for sequencing of 150 cycles paired-end. The 5R-PLEX single index list can be found in the Appendix A in the Product Information Sheet. The Product Information Sheet can be downloaded from the 5R-PLEX product page at [SigmaAldrich.com](https://www.sigmaaldrich.com).

Bioinformatic Analysis

- Create account and log-in to m-camp.info/microbiome.
- Use the Key-code printed on the Quick Start Guide provided in the kit packaging.
- Upload fastQ in the 5R-PLEX module and follow the instructions in the Product Information Sheet.

Additional Information

The Product Information Sheet and other documents, are available on the product page at [SigmaAldrich.com](https://www.sigmaaldrich.com).

Notice

We provide information and advice to our customers on application technologies and regulatory matters to the best of our knowledge and ability, but without obligation or liability. Existing laws and regulations are to be observed in all cases by our customers. This also applies in respect to any rights of third parties. Our information and advice do not relieve our customers of their own responsibility for checking the suitability of our products for the envisaged purpose.

The information in this document is subject to change without notice and should not be construed as a commitment by the manufacturing or selling entity, or an affiliate. We assume no responsibility for any errors that may appear in this document.

Technical Assistance

Visit the tech service page on our web site at SigmaAldrich.com/TechService.

Terms and Conditions of Sale

Warranty, use restrictions, and other conditions of sale may be found at SigmaAldrich.com/Terms.

Contact Information

For the location of the office nearest you, go to SigmaAldrich.com/Offices.

Merck, and Sigma-Aldrich are trademarks of Merck KGaA, Darmstadt, Germany or its affiliates. All other trademarks are the property of their respective owners. Detailed information on trademarks is available via publicly accessible resources.
© 2023 Merck KGaA, Darmstadt, Germany and/or its affiliates. All Rights Reserved.

The life science business of Merck operates as MilliporeSigma in the U.S. and Canada.