

Product Information

SEQPLEX™-I WGA Kit

Whole Genome Amplification, DNA Amplification

SeqXi

Storage temperature –20 °C

Product Description

The SeqPlex™-I DNA Amplification Kit for whole genome amplification (WGA) is designed to facilitate Illumina next-generation sequencing (NGS) from extremely small quantities or from degraded/highly fragmented DNA. The SeqPlex™-I WGA kit provides a ligation-free workflow for amplification of such samples for direct input onto Illumina Next Generation Sequencing (NGS) flow cells. The SeqPlex™-I process is comprised of three steps: Pre-amplification/Library Synthesis, Amplification 1 and Amplification 2. See Process Workflow chart on page 2.

First, template DNA is replicated using primers comprised of semi-degenerate 3'- ends with universal 5'- ends. As polymerization proceeds, displaced single strands serve as new templates for additional primer annealing and extension, producing random, overlapping replicons flanked by a universal primer (5') and primer complement (3') sequence.

Next, products from pre-amplification/library synthesis are amplified by single primer PCR via the universal end sequence. These amplification products typically range from 200 to 500+ base pairs. Amplicons from ChIP and/or degraded DNA, such as Formalin Fixed Paraffin Embedded (FFPE), are typically shorter and dependent upon length of the starting DNA.

Last, single primer amplicons from amplification 1 are converted to dual Illumina primer PCR products ready for purification, quantification, and Illumina NGS. Amplification 1 amplicons are amplifiable with Illumina adapter-containing primers ending with 5'- ...GCTCTTCCGATCT-3' such as:

AATGATACGGCGACCACCGAGATCTACAC [i5 index]
ACACTCTTCCCTACACGACGCTCTTCCGATCT

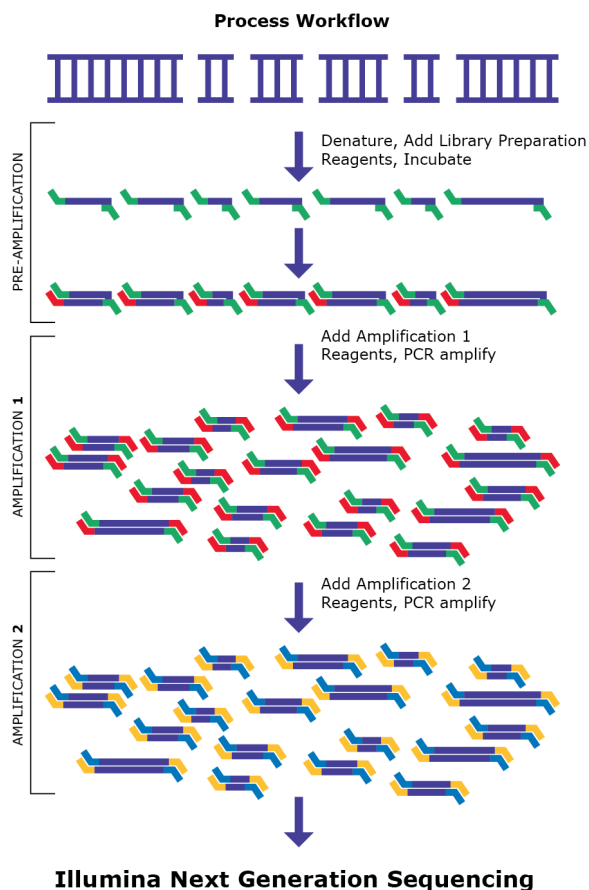
CAAGCAGAAGACGGCATAACGAGAT [i7 index]
GTGACTGGAGTTCAGACGTGTGCTCTTCCGATCT

The example index sequences below are ready to be added to the adapter sequences as is, without reverse complement of the i7 indexes.

Examples of i5 indexes	Examples of i7 indexes
AP100 – AGCGCTAG	AP100 – CCGCGGTT
D501 – TATAGCCT	D701 – CGAGTAAT
D502 – ATAGAGGC	D702 – TCTCCGGA
D503 – CCTATCCT	D703 – AATGAGCG
D504 – GGCTCTGA	D704 – AATGAGCG
D505 – AGGCGAAG	D705 – TTCTGAAT

Note: Entering i5 and i7 index sequences into the Illumina sample sheet is machine specific. The table above shows examples of both index adapters that can be used in many combinations, in addition to the set included in the kit as AP100.

A 20 µL Amplification 2 reaction will produce > 100 ng of amplified double-stranded DNA when starting with 100 pg to 5 ng of high-quality DNA. Higher input quantities and higher quality DNA template generally result in increased yields. For damaged DNA, such as from FFPE, 1-50 ng input DNA is recommended. Reaction volumes can be scaled up or down to accommodate preparation of desired quantities of final product. Pre-amplification/Library synthesis reactions may be scaled to as low as 5 µL. Sufficient reagents have been supplied for the number of indicated kit reactions, at Pre-amplification/Library Synthesis, Amplification 1 and Amplification 2 volumes of 15, 30 and 20 µL respectively. Scaling up will reduce the overall number of reactions that the kit can accommodate. Use the same library synthesis volumes for all samples to be compared.



The SeqPlex-I WGA workflow consists of three phases.

Pre-Amplification

- DNA is replicated using primers composed of a semi-degenerate 3'-end and a universal 5'-end that amplify overlapping fragments with comprehensive genome coverage.
- Input DNA may be intact or fragmented.
- Compatible with low-quantity ChIP or FFPE samples.

Amplification 1

- Pre-Amplification products are amplified by single primer PCR via the proprietary universal end sequence.
- Cycle numbers are variable, depending on quantity and quality of starting sample; amplification is monitored by SYBR green and stopped after signal plateau.

Amplification 2

- Amplification 1 products are converted to dual Illumina-adapted products using adapter primers.
- Adapter primers should be custom ordered with indexes for multiplex capability.
- Amplicon sizes average about 300 bp and are ideal for NGS.
- Following Amplification 2, samples are ready for clean up, quantification, and Illumina NGS.

Reagents Provided

Reagent	Reagent No.	24 RXN	96 RXN	384 RXN
10X Library Synthesis Solution for SeqXi	LP200	36 µL	144 µL	576 µL
10X Library Synthesis Buffer for SeqXi	LB200	36 µL	144 µL	576 µL
Library Preparation Enzyme for SeqXi	EO600	14.4 µL	57.6 µL	230.4 µL
2X Amplification 1 Mix for Seq-I	A8112	345.6 µL	1382.4 µL	5529.6 µL
Amplification 1 Polymerase for Seq-I	SP500	14.4 µL	57.6 µL	230.4 µL
5X Amplification 2 Mix for Seq-I	BA400	86.4 µL	345.6 µL	1382.4 µL
Amplification 2 Enzyme for Seq-I	BA500	12 µL	48 µL	192 µL
Dual Index Adapter Primers for Seq-I	AP100	48 µL	192 µL	768 µL
Water, Molecular Biology Reagent	W4502	5 mL	10 mL	20 mL

Materials and Reagents Required

(But not provided)

- Thermal cycler & qPCR instrument
- Sample DNA
- Spectrophotometer, NanoDrop or similar
- 0.2 mL Thin-Walled PCR tubes, strip or plates
- Pipettors
- Pipette tips with aerosol barriers
- SYBR® Green I, (Cat. No. S9430) (optional)
- GenElute™ PCR Clean-up Kit, (Cat. No. NA1020)
- Additional index primers if multiplexing samples (Cat. No. OLIGO)

Precautions and Disclaimer

The SeqPlex™-I DNA Amplification Kit for whole genome amplification (WGA) is for research use only. Not for drug, household, or other uses. Please consult the Safety Data Sheet for information regarding hazards and safe handling practices.

Storage/Stability

All components should be stored at $-20\text{ }^{\circ}\text{C}$. When thawed for use, components should be kept on ice. Dissolve any precipitate in these solutions by briefly heating at $37\text{ }^{\circ}\text{C}$, with thorough mixing. Stability of the EO600, SP500 and BA500 will be affected if stored above $-20\text{ }^{\circ}\text{C}$ or allowed to remain for long periods at temperatures over $4\text{ }^{\circ}\text{C}$.

Procedure

The following procedure has been used successfully to amplify and sequence from small quantities, such as single cell and/or degraded/fragmented DNA, such as that from ChIP-isolated or FFPE tissues.

Note: Final yield after amplification varies significantly depending upon the quality of starting DNA. In most cases, $> 100\text{ ng}$ can be expected. If larger quantities are needed, reactions can be scaled up to accommodate this need, but will limit the total reactions in the kit.

This procedure was developed using the specific reagents provided with or recommended for use with this kit. Substitutions may result in suboptimal results.

Pre-amplification/Library Synthesis

Most DNA has some level of fragmentation, such as DNA from many ChIP or FFPE samples, which are ready for input. Most DNA samples do not require additional fragmentation before inputting into this method. If needed, DNA may be fragmented using sonication, cavitation, enzymatic or thermal fragmentation.

1. Thaw the 10X Library Synthesis Solution for SeqXi (LP200), 10X Library Synthesis Buffer for SeqXi (LB200) and water (W4502). Mix thoroughly before use.

Combine 100 pg to 10 ng of high-quality DNA, $1\text{-}50\text{ ng}$ FFPE DNA or a single cell with LP200 at the following single-reaction scale:

1.5 μL of 10X Library Synthesis Solution for SeqXI (LP200)

X μL DNA (for example: 1 ng)

Y μL water (W4502)

12.9 μL Total reaction volume

2. Mix thoroughly, centrifuge briefly, and incubate in a thermal cycler programmed for:
 - 95 $^{\circ}\text{C}$ for 2 minutes
 - 16 $^{\circ}\text{C}$ for 5 minutes
3. Add 1.5 μL of 10X Library Synthesis Buffer for SeqXi (LB200) and 0.6 μL of Library Preparation Enzyme for SeqXi (EO600) to the sample.
 - 1.5 μL of 10X Library Synthesis Buffer for SeqXi (LB200)
 - 0.6 μL of Library Preparation Enzyme (EO600)
 - 12.9 μL from previous step

15 μL Total reaction volume

Cap tube and mix thoroughly. Centrifuge briefly and immediately proceed to next step. (For multiple reactions, a master mix comprised of LB200 and EO600 may be prepared. Add 2.1 μL of the master mix to each sample).

4. Place reaction(s) in a thermal cycler and incubate:
 - 16 $^{\circ}\text{C}$ for 10 minutes
 - 25 $^{\circ}\text{C}$ for 10 minutes
 - 37 $^{\circ}\text{C}$ for 10 minutes
 - 16 $^{\circ}\text{C}$ for 10 minutes
 - 25 $^{\circ}\text{C}$ for 10 minutes
 - 37 $^{\circ}\text{C}$ for 10 minutes
 - 42 $^{\circ}\text{C}$ for 5 minutes
 - 95 $^{\circ}\text{C}$ for 1 minutes
 - 4 $^{\circ}\text{C}$ Hold

Remove reaction(s) from thermal cycler and centrifuge briefly. Amplification 1 may be started immediately or store Pre-Amplification product at $-20\text{ }^{\circ}\text{C}$ for up to three days.

Caution—Experienced WGA and WTA users:

Several components found in the SeqPlex RNA Amplification kit (SEQR), SeqPlex Enhanced DNA Amplification kit (SEQXE), GenomePlex WGA kits, Transplex® WTA1 kit and Complete Whole Transcriptome Amplification Kit (WTA2) are similarly named. Though generally analogous in function, they are not interchangeable.

Amplification 1

5. Add the following reagents to the 15 μL of Pre-amplification/Library Synthesis product from Step 4. (For multiple reactions, a master mix comprised of reagents listed below may be prepared. Add 15/15.6 μL +/- SYBR[®] of the master mix to each reaction):

14.4 μL 2X Amplification 1 Mix
for Seq-I (A8112)

0.6 μL Amplification 1 Polymerase
for Seq-I (SP500)

0.6 μL SYBR[®] Green I (S9430) diluted
1/1000* (recommended)

15.0 μL Pre-amp/Lib. Synthesis from Step 4
- μL Instrument Specific Reference Dye
(optional, if needed add to final total)

**30/30.6 μL (-/+ SYBR[®])
Total reaction volume**

* - For the best representation, real-time PCR with addition of freshly diluted SYBR[®] Green I to the amplification reaction is strongly recommended to enable monitoring of the reaction progress. SYBR[®] Green I (S9430) must be diluted 1,000-fold (1/1000) in water and 0.6 μL used per 30 μL of amplification 1 reaction for a total of 30.6 μL to avoid inhibiting the amplification reaction. SYBR[®] Green I formulations other than S9430 have not been tested and are not recommended.

Optimal results are achieved by not proceeding past the amplification "plateau" (see **Figure 1**). The optimal number of amplification cycles varies with starting DNA template quantity and quality.

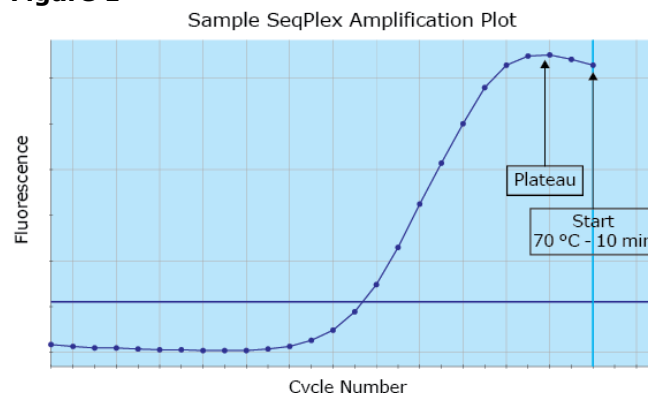
If amplification is performed without adding SYBR[®] Green I, 17-22 cycles will usually give good results with 0.1-1.0 ng of high-quality DNA. Low quality DNA may require higher input quantities and/or more cycles. If input amounts are near 10 pg or lower, as many as 29 cycles may be required to reach amplification plateau.

DNA Input Amount and Typical # of Cycles to Reach Plateau Reference

Input Quantity	# of Cycles
10 pg	25
100 pg	22
1 ng	17
10 ng	14

Note: If more than 29 cycles are required to achieve plateau, subsequent NGS results may be unsatisfactory. Consult the Troubleshooting Guide.

Figure 1



6. Cap tube(s) and mix thoroughly. Centrifuge briefly and cycle in a real-time thermal cycler:

Initial Denaturation:

94 °C for 2 minutes

Cycle until plateau: (< 25)

94 °C Denature for 15 seconds

70 °C Anneal/Extend for 5 minutes
(read fluorescence)

After cycling:

70 °C for 10 minutes

4 °C Hold

After cycling is complete, remove reaction(s) and centrifuge briefly. Amplification 2 may be started immediately or store Amplification 1 product at -20 °C for up to three days.

Caution—Experienced WGA and WTA users:

- SeqPlex[™]-I uses a 2X Amplification Mix for Amplification 1.
- SYBR[®] Green I (S9340) is recommended to monitor the amplification.
- Annealing/Extension temperature is 70 °C.
- A 10-min 70 °C post-amp 1 hold is required.

Amplification 2

7. Add the following reagents to a new tube or prepare a master mix for multiple reactions (Add 18 μL of the master mix to each tube). You will use 2 μL of the Amplification 1 product from Step 6 per reaction:

3.6 μL 5X Amplification 2 Mix for Seq-I (BA400)

0.5 μL Amplification 2 Enzyme for Seq-I (BA500)

2.0 μL Dual Index Adapter Primers for Seq-I (AP100)

11.5 μL Water, Molecular Biology Reagent (W4502)

0.4 μL SYBR[®] Green I (S9430) diluted 1/1000*

2.0 μL Amplification 1 product from Step 6

- μL Instrument Specific Reference Dye (optional, if needed add to final total)

20.0 μL Total reaction volume

Note: The dual index adapter primers (AP100; see table on page 1 for i5 and i7 index sequences) provided in this kit will only work for one set of samples. If pooling samples for sequencing is required, the user must provide additional index primer sets. See example index primer sequences page 1. Using index primers at a concentration of 0.2 μM to 2 μM each is recommended. When used as instructed, the final concentrations of the provided dual index adapter primers (AP100) are 0.2 μM . Lower primer concentration results in higher single stranded hetero duplexes while higher concentrations produce more fully duplex product. Higher concentrations also produce more adapter dimers.

* - For the best representation, real-time PCR with addition of freshly diluted SYBR[®] Green I to the amplification reaction is strongly recommended to enable monitoring of the reaction progress. SYBR[®] Green I (S9430) must be diluted 1,000-fold (1/1000) in water and 0.4 μL used per 20 μL of amplification 2 reaction to avoid inhibiting the amplification reaction. SYBR[®] Green I formulations other than S9430 have not been tested and are not recommended.

Optimal results are achieved by not proceeding past the amplification "plateau". Proceeding past 1-2 cycles after "plateau" may increase bias and decrease amplicon representation. The optimal number of amplification cycles varies with starting DNA template quantity and quality.

If amplification is performed without adding SYBR[®] Green I, 8-12 cycles will usually yield sufficient product. Low quality DNA may require higher input quantities and/or more cycles as observed in the amplification 1 cycling.

8. Cap tube(s) and mix thoroughly. Centrifuge briefly and cycle in a real-time thermal cycler:

Initial Incubation/Denaturation:

37 °C for 5 minutes

94 °C for 2 minutes

Two Cycles:

94 °C for 30 seconds

60 °C for 5 minutes

70 °C for 1 minute

Cycle until plateau: (< 12)

94 °C Denature for 15 seconds

70 °C Anneal/Extend for 1 minute (read fluorescence)

After cycling:

70 °C for 10 minutes

4 °C Hold

Note: The extended incubation at 70 °C after cycling is essential for maximal duplex synthesis.

After cycling is complete, remove reaction(s) and centrifuge briefly. Reactions can be purified now or stored at -20 °C for up to three days.

Caution—Experienced WGA and WTA users:

- SeqPlex[™]-I uses a 5X Amplification Mix for Amplification 2.
- SYBR[®] Green I (S9340) is recommended to monitor the amplification.
- Annealing/Extension temperature is 70 °C.
- A 10-minute 70 °C post-amp 2 hold is required.

Purification

The final Amplification 2 reaction(s) can be purified using the GenElute™ PCR Clean-Up Kit (NA1020). Follow the Technical Bulletin for the kit. Elute in 50 µL nuclease-free water, instead of the Elution Solution provided in the kit. Alternatively, the reactions can be purified with SPRI magnetic beads. In order to minimize sequencing of primer dimers, use a clean up method that removes fragments less than 200 bp in length. The purified reaction(s) can be used immediately or stored at –20 °C for several weeks.

Yield Determination

The amplification product is generally a mixture of single- and double- stranded amplicons. Yield will vary depending on the quality of starting DNA. Concentration can be measured by:

- Traditional A₂₆₀ absorbance can be used to measure concentration. One A₂₆₀ unit is equivalent to 50 ng/µL dsDNA.
- Fluorescent dyes such as PicoGreen®. Duplex-specific dyes may underestimate the actual DNA yield.
- Quantitative PCR-based library quantification methods are recommended, as quantification is independent of strandedness.

Optional: Amplification quality may be assessed by capillary electrophoresis. Typically, 1 µL of crude or purified amplification product is sufficient for capillary chips, such as those for Agilent's Bioanalyzer or TapeStation. The amplification product is generally a mixture of single and double stranded amplicons.

Sequencing

The SeqPlex-I generated library is now ready to enter Illumina Next-Generation Sequencing (NGS). The first nine bases of each read should be disregarded, as these bases result from the pre-amplification/library synthesis primers used in that section of the kit.

Frequently Asked Questions

1. Is SeqPlex™-I WGA Kit compatible with microarrays and qPCR?
Yes, libraries made using SeqPlex™-I WGA Kit may be used in these applications like genomic DNA or existing GenomePlex products.
2. Are there advantages to SeqPlex™-I over GenomePlex?
SeqPlex™-I Pre-Amplification primers have been designed to target more frequently than existing GenomePlex WGA primers and therefore may provide the advantage of superior genome coverage in some regions.

3. Will reducing cycles during amplification improve representation?
No, you need to reach "plateau" for optimum representation. Proceeding past 1-2 cycles will not negate the reactions, but best representation is achieved around "plateau". Insufficient cycling leads to a significant reduction in representation/coverage.
4. Will SeqPlex™-I libraries require special NGS sequencing protocols?
No, SeqPlex™-I libraries fit directly into NGS workflows. Sequencing instrument operators should be notified of running SeqPlex™-I DNA and to expect a slight signal from any remaining primers in the IVC plots.

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 at [SigmaAldrich.com/techservice](https://www.sigmaaldrich.com/techservice).

Standard Warranty

The applicable warranty for the products listed in this publication may be found at [SigmaAldrich.com/terms](https://www.sigmaaldrich.com/terms).

Contact Information

For the location of the office nearest you, go to [SigmaAldrich.com/offices](https://www.sigmaaldrich.com/offices).

SeqXi Experienced User Protocol

Pre-Amplification/Library Synthesis

Reagent	Step 1	Step 2
DNA	1 µL to 10 µL	Incubate 95 °C — 2 min 16 °C — 5 min Add reagents in Step 3 to samples
10X Library Synthesis Solution (LP200)	1.5 µL	
Water (W4502)	10.4 µL to 1.4 µL	
Total	12.9 µL	
Reagent	Step 3	Step 4
Reaction from Step 2	(12.9 µL)	PCR 16 °C — 10 min } 25 °C — 10 min } 2 cycles 37 °C — 10 min } 42 °C — 5 min 95 °C — 1 min 4 °C — hold
10X Library Synthesis Buffer (LB200)	1.5 µL	
Library Preparation Enzyme (EO600)	0.6 µL	
Total	15 µL	

Amplification 1

Reagent	Step 5	Step 6
2X Amplification 1 Mix for Seq-I (A8112)	14.4 µL	PCR 94 °C — 2 min 94 °C — 15 sec } 70 °C — 5 min } < 25 cycles (monitor) 70 °C — 10 min 4 °C — hold
Amplification 1 Polymerase for Seq-I (SP500)	0.6 µL	
SYBR® Green (diluted 1:1000) <i>(S9430, not included in kit, but recommended for monitoring cycling)</i>	(0.6 µL)	
Pre-Amplification/Library Synthesis product from Step 4	15 µL	
Total	30.0 (30.6) µL	

Amplification 2

Reagent	Step 7	Step 8
5X Amplification 2 Mix for Seq-I (BA400)	3.6 µL	PCR 37 °C — 5 min 94 °C — 2 min 94 °C — 30 sec } 60 °C — 5 min } 2 cycles 70 °C — 1 min } 94 °C — 15 sec } 70 °C — 1 min } < 12 cycles (monitor) 70 °C — 10 min 4 °C — hold
Amplification 2 Enzyme for Seq-I (BA500)	0.5 µL	
Dual Index Adapter Primers for Seq-I (AP100; see page 1 for i5 and i7 index sequences)	2 µL	
Water (W4502)	11.5 or 11.9 µL	
SYBR® Green (diluted 1:1000) <i>(S9430, not included in kit, but recommended for monitoring cycling)</i>	(0.4 µL)	
Amplification 1 Reaction (undiluted) from Step 6	2 µL	
Total	20 µL	

Troubleshooting Guide

Observation	Cause	Recommended Solution
No product was detected after amplification	Incorrect Annealing/Extension temperature or time	Perform Amp 1 reaction again with 70 °C Annealing/Extension for 5 minutes each cycle.
	Too few PCR cycles during Amplification	Perform reaction again with more cycles (up to 29) and monitor amplification with SYBR® Green in a real-time thermal cycler.
	Starting DNA was insufficient or too severely degraded	Perform reaction again with more starting DNA.
	DNA may be single stranded	Use a non-duplex-dependent endpoint method (such as qPCR) to detect DNA amplification.
No amplification curve was seen during real-time PCR (qPCR) monitoring	SYBR® Green was not added	SYBR® Green I, Cat. No. S9430, is not included with SeqPlex™-I reagents, but must be added to monitor real-time PCR.
	A reference dye specific to the qPCR instrument may be required	Add instrument specific reference dye. If adding reference dye is not possible, excess cycles should be completed to ensure complete cycling.
Presence of enriched small product(s), particularly 176 bp, after Amplification 2	Primer dimers	Decrease concentration of index adapter primers used in amplification 2. It is recommended to use primers at a final concentration of 0.2 uM to 2 uM; high concentrations increase risk of primer dimer formation.
		Increase template concentration; primer dimer formation decreases with additional template. If primer dimer cannot be eliminated by decreasing adapter primer concentration or increasing template concentration, primers can be removed using a left side size selection or double size selection with SPRI beads.
Poor representation after NGS	Input DNA not fragmented enough	Verify size and fragment if necessary prior to starting the SeqPlex procedure.
	Not enough cycles during SeqPlex Amplification	Monitor amplification cycling and cycle until plateau is reached for both Amp1 and Amp 2. If monitoring completely fails, a default 29 cycles may be used for Amp 1 and 12 cycles for Amp 2.
NGS IVC abnormalities	IVC plots show identical sequence (primer) for the first few nucleotides in multiple reads	Optimum instrument cluster calling may be achieved by normalizing the run to a lane that does not contain SeqPlex™-I DNA.

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

MilliporeSigma, 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.
© 2022 Merck KGaA, Darmstadt, Germany and/or its affiliates. All Rights Reserved.