

FLASH-seq protocol with SEQURNA (Cat. No. 9028)

In Brief

SEQURNA is a synthetic thermostable RNase inhibitor for single-cell RNA-sequencing (scRNAseq), yielding single-cell libraries of equal or superior quality compared to ubiquitously used protein-based recombinant RNase inhibitors (RRIs).

SEQURNA provides additional unique improvements in reproducibility and throughput, enables new experimental workflows including retained RNase inhibition throughout heat cycles, and can reduce the need for dry-ice transport.

The following protocol was developed by Khven et al. 2025¹ in the Picelli lab (Institute of Molecular and Clinical Ophthalmology, Basel) and is a modification of the original FLASH-seq protocol by Hahaut et al. 2022².

Important Information

- **No additional RNase inhibitor should be included in the RT mix.** SEQURNA in cell lysis buffer remains active throughout cell lysis (by heating) and the reverse transcription (RT) reaction. Adding additional RNase inhibitor in the RT mix may decrease performance of the protocol.
- **Dimethylformamide (DMF) is toxic** and should be handled under a fume hood, or in accordance to local safety regulations.
- Reagent mixes should be prepared shortly before use.
- Mix thoroughly each mix before dispensing. For higher accuracy use liquid handling robots and/or nano-dispensers whenever possible. In FLASH-seq, the Picelli lab has used the I.DOT (Dispensix) for all the dispensing steps and the Fluent 780 liquid handling robot (Tecan) for sample cleanup, reagent transfers and pooling.
- The protocol described below is meant to be carried out in 384-well plates. There is no need to use mineral oil to prevent evaporation.
- The protocol below uses 1 U/μl of SEQURNA inhibitor. This represents the final concentration in the RT-PCR step. The range 0.75–1 U/μl SEQURNA in the RT-PCR step has been shown by benchmarking to be the most suitable one for the FLASH-seq and results in markedly improved gene detection compared to the use of standard RRIs¹. Note that the optimal SEQURNA concentration may differ for other single-cell RNA-seq protocols, even if using the same SMART-seq chemistry, due to volumetric changes, buffer conditions, and enzymes used in the protocols.
- Always use low-binding (LoBind) plates and tubes (especially for long-term storage) to prevent RNA/cDNA/ amplified DNA from sticking to plastic.

- Keep 384-well PCR plates on a thermoconductive cooling module or rack placed on ice (e.g., SEQblock™ PCR 384-well (SQ-PCR-384, Genovis AB)). This maintains optimal low temperatures for samples and reaction mixtures throughout setup, improves handling convenience, and prevents condensation from forming on the exterior of the plates. Water droplets can otherwise introduce lime, salt, or mineral deposits into the thermocycler heat block over time, which gradually impairs heat transfer efficiency and compromises thermal cycling performance of the thermocycler.

Oligonucleotide Sequences (5' to 3')

Standard FLASH-Seq RT-PCR

Smart dT30VN: 5'-/5Biosg/AAGCAGTGGTATCAACGCAGAGTACT30VN (desalted or HPLC)

FS TSO: 5'- /5Biosg/AAGCAGTGGTATCAACGCAGAGTACrGrGrG (desalted or HPLC)

1. Preparation of Lysis Mix (15 min.)

1.1 Prepare the following lysis buffer mix

Reagent	Conc. in lysis buffer	μl per reaction	384-well plate (422.4 rxns)
Triton-X100 (10% v/v)	0.2%	0.020	8.448
dNTP mix (25 mM each)	6 mM	0.240	101.376
SMART dT30VN (100 μM)	1.8 mM	0.018	7.603
SEQURNA (50 mass units/μl)	5 U/μl*	0.100*	42.240*
Betaine (5 M)	0.2 M	0.200	84.480
Nuclease-free water	-	0.422	178.253
Total	-	1 μl	422.4 μl

* This amount results in 1 mass unit/μl SEQURNA in the downstream RT-PCR reaction.

Table 1. Reagent preparation for lysis buffer: Volumes for 384-well plates

IMPORTANT! NO DTT is used in this version of FLASH-seq. Including DTT will decrease performance of the protocol.

1.2 Add 1 μl Lysis Mix to each well of a 384-well plate.

1.3 Seal the plate with a PCR seal and quickly spin it down to collect the Lysis Mix to the bottom.

1.4 Proceed immediately to the next step or store the plate at -20 °C long-term. Plates that are going to be used on the same day can be stored in the fridge or kept on ice.

SAFE STOPPING POINT – Plates containing lysis buffer can be stored for >6 months at -20 °C.

NOTICE

R&D Systems™ now includes the legacy brands Novus Biologicals™, Tocris Bioscience™, and ProteinSimple™.

Bio-Techne Spatial now includes the legacy brands Lunaphore™ and Advanced Cell Diagnostics™.

Bio-Techne Diagnostics now includes Asuragen®, Bionostics™, Clinica™, RNA Medical®, and R&D Systems™ Clinical Controls.

2. Sample Collection (10 min.)

- 2.1 Sort single cells into 384-well plates containing 1 µl of Lysis Mix.
- 2.2 Seal the plate with an aluminum seal. If processing multiple plates at once, keep each plate on dry ice until ready to transfer them all to -80 °C for long-term storage.
- 2.3 Even if proceeding with the protocol immediately after sorting, it is advisable to put the plate on dry ice for 5 minutes, followed by heat denaturation (Cell Lysis step), as freeze-thawing facilitates cell lysis.

SAFE STOPPING POINT – Sorted cells in lysis buffer can be stored for >6 months at -80 °C. Longer storage might lead to lower yield or increased presence of shorter cDNA fragments.

3. Cell Lysis (3 min.)

- 3.1 Remove the plates from the -80 °C freezer and check that the aluminum seal is still intact. If damaged or not sticking to the plate anymore, wait a few minutes for the plate to partially thaw, remove the damaged foil and replace it with a new one.
- 3.2 Place the plate in a thermocycler with a heated lid and incubate for 3 minutes at 72 °C, followed by a 4 °C hold step.
- 3.3 Spin down any condensation droplets that may have formed during the incubation and return the plate to a cooling rack on ice (e.g. SEQblock™ PCR 384-well (SQ-PCR-384, Genovis AB)). Proceed quickly to the next step. If it is not ready with the RT-PCR mix, keep the plate on the cooling rack on ice at all times.

4. RT-PCR Reaction (3h 30 min.)

- 4.1 While the plate is in the thermocycler (Step 3, Cell Lysis), prepare the following RT-PCR Mix:

Reagent	Conc. in RT	µl per reaction	384-well plate (422.4 rxns)
MgCl ₂ (1 M)	9.2 mM	0.046	19.430
Betaine (5 M)	1 M	0.800	337.920
Nuclease-free water	-	0.422	178.253
dCTP (100 mM)	1.8 mM	0.090	38.016
Maxima H- RT (200 U/µl)	2 U/ul	0.050	21.120
KAPA HiFi HotStart ReadyMix (2x)	1x	2.500	1056.000
FS TSO (100 µM)	1.84 µM	0.092	38.861
Total volume (µl)		4.000	1689.600
Total	-	4 µl	1689.600 µl

Table 2. Reagent preparation for reverse transcription reaction: Volumes for 384-well plates

IMPORTANT! NO additional RNase inhibitor should be included in the RT mix. Adding additional RNase inhibitor in this step will decrease performance of the protocol. **NO DTT** is used in this version of FLASH-seq. Including DTT will decrease performance of the protocol.

NOTICE

R&D Systems™ now includes the legacy brands Novus Biologicals™, Tocris Bioscience™, and ProteinSimple™.

Bio-Techne Spatial now includes the legacy brands Lunaphore™ and Advanced Cell Diagnostics™.

Bio-Techne Diagnostics now includes Asuragen®, Bionostics™, Clinica™, RNA Medical®, and R&D Systems™ Clinical Controls.

- 4.2 Add 4 µl of the RT-PCR Mix into each well of the 384-well plate.
- 4.3 Seal the plate with a PCR seal, gently vortex and spin down to collect the liquid at the bottom.
- 4.4 Place it in a thermocycler with heated lid and start the following RT-PCR program:

Condition	Step	Temperature	Time	Cycles
RT		50 °C	60 min	1x
PCR	Initial Denaturation	98 °C	3 min	1x
	Denaturation	98 °C	20 sec	18-21x*
	Annealing	67 °C	20 sec	
	Elongation	72 °C	5 min	
		15 °C	Hold	

* Adjust the number of cycles according to the cell type used. We recommend 18-19 cycles for HEK 293T cells and 21 cycles for hPBMC.

Table 3. Thermocycling conditions for reverse transcription/PCR

SAFE STOPPING POINT – Amplified cDNA before purification can be stored for several months at -20 °C.

5. cDNA Purification (30 min.)

- 5.1 You can either use AMPure XP beads, SPRI beads, or prepare your own in-house solution of SeraMag beads containing 18% w/v PEG to reduce costs. A detailed protocol for making your own magnetic bead solution is described in Picelli S, Methods Mol Biol. 2019:1979:25-44⁵.
- 5.2 Bring out the magnetic bead working solution from the +4 °C (fridge) storage and equilibrate at room temperature for 15 min. Vortex the bead solution.
- 5.3 We recommend adding extra nuclease-free water to each sample, to increase the volume, simplify the handling and improve recovery rate from the 384-well plates. We generally add 10 µl of nuclease-free water to 5 µl of amplified cDNA.
- 5.4 Add a 0.8× volume ratio of magnetic bead working solution to each well (i.e., 12 µl beads for each 15 µl cDNA). Mix thoroughly by pipetting or vortexing.
- 5.5 Incubate the plate off the magnetic stand for 5 min. at room temperature.
- 5.6 Place the plate on the magnetic stand and leave it for 5 min. or until the solution appears clear.
- 5.7 Remove the supernatant without disturbing the beads.
- 5.8 Performing an ethanol wash is not necessary. We do not recommend it when working in 384-well plates and with liquid handling robots, to avoid cDNA losses.

NOTICE

R&D Systems™ now includes the legacy brands Novus Biologicals™, Tocris Bioscience™, and ProteinSimple™.

Bio-Techne Spatial now includes the legacy brands Lunaphore™ and Advanced Cell Diagnostics™.

Bio-Techne Diagnostics now includes Asuragen®, Bionostics™, Clinika™, RNA Medical®, and R&D Systems™ Clinical Controls.

5.9 Remove the plate from the magnetic stand, add 15 μ l of nuclease-free water and mix well by pipetting or vortexing to resuspend the beads. Do not let the bead pellet dry, as it will decrease the final cDNA yield.

5.10 Incubate for 2 min off the magnetic stand.

5.11 Place the plate back on the magnetic stand and incubate for 2 min or until the solution appears clear.

5.12 Remove 14 μ l of the supernatant, containing purified cDNA, and transfer it to a new plate.

SAFE STOPPING POINT – Amplified and purified cDNA can be stored for several months at -20°C . We recommend using LoBind plates to avoid material losses upon long-term storage.

6. cDNA Quality Control Check (45 min.)

6.1 Check the cDNA quality on Agilent Bioanalyzer High Sensitivity DNA chip, or similar equipment. Follow the instructions described in the user manual. A high-quality cDNA library is characterized by a low proportion of fragments <500 bp, absence of residual primers (ca. 100 bp) and an average cDNA size of 2.0–2.5 Kb. A representative Bioanalyzer trace of successfully amplified FLASH-seq cDNA from a HEK cell (19 cycles) using SEQRNA is shown in Figure 1.

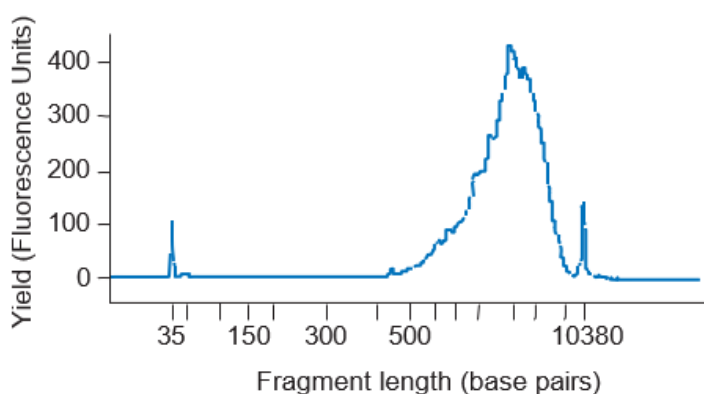


Figure 1. Example of cDNA trace from a HEK cell (19 PCR cycles), using an Agilent Bioanalyzer High Sensitivity DNA Analysis chip (Agilent). x-axis: cDNA yield (fluorescence units), y-axis: fragment length (base pairs). Data obtained from Picelli et al. 2025.³

7. cDNA Quantification (15 min.)

7.1 Allow the Quant-iT PicoGreen reagent to warm to room temperature before opening the vial. PicoGreen is light sensitive; while thawing, wrap in aluminum foil.

7.2 Prepare a 1 $\times$ working solution of TE using 20 $\times$ TE (supplied) and nuclease-free water.

7.3 Prepare a 1:800 dilution of PicoGreen solution and always use a plastic vessel (tubes, Falcon, etc.). Do not use glass as PicoGreen may adsorb to glass.

NOTICE

R&D Systems™ now includes the legacy brands Novus Biologicals™, Tocris Bioscience™, and ProteinSimple™.

Bio-Techne Spatial now includes the legacy brands Lunaphore™ and Advanced Cell Diagnostics™.

Bio-Techne Diagnostics now includes Asuragen®, Bionostics™, Clinika™, RNA Medical®, and R&D Systems™ Clinical Controls.

7.4 Prepare the standard curve using Lambda DNA standard (supplied at a concentration of 100 ng/μl, with the PicoGreen kit) and 1x TE in 8 tubes, as below. The stock tubes can be used multiple times, keep any leftover in the fridge at +4 °C between experiments.

7.5 Vortex well and spin down the DNA standards before every use. Not vortexing thoroughly the standards is going to negatively affect the standard curve and your readings. Serial dilutions should be prepared as shown in the table below.

Tube no.	Contents	Concentration	Final volume
1	90 μl TE + 10 μl Lambda DNA stock	10 ng/μl	100 μl
2	50 μl from Tube 1 + 50 μl TE	5 ng/μl	100 μl
3	50 μl from Tube 2 + 50 μl TE	2.5 ng/μl	100 μl
4	50 μl from Tube 3 + 50 μl TE	1.25 ng/μl	100 μl
5	50 μl from Tube 4 + 50 μl TE	0.625 ng/μl	100 μl
6	50 μl from Tube 5 + 50 μl TE	0.3125 ng/μl	100 μl
7	50 μl from Tube 6 + 50 μl TE	0.15625 ng/μl	100 μl
8	TE only	blank	-

7.6 Prepare the PicoGreen solution by pipetting 0.25 μl of PicoGreen dye + 99.5 μl of 1X TE for each sample. Vortex to mix.

7.7 Pipette 1 μl of each of the 7 standards + 1 Blank into a black, flat-bottom Nunc™ F96 MicroWell™ Plate. Place the standards on one column.

7.8 Pipet 1 μl of your samples into the center of each well of the Nunc™ F96 MicroWell™ Polystyrene Plate.

7.9 Add 99 μl of PicoGreen + TE mix into every well. There is no need to mix.

7.10 Cover the plate with the provided plastic (transparent) lid to prevent possible contamination.

8. Plate Normalization (10 min.)

8.1 Prepare a normalization plate by adding 1 μl of purified cDNA and nuclease-free water to a final concentration of 200 pg/μl.

SAFE STOPPING POINT – Normalized cDNA can be stored for several months at -20°C. LoBind plates must be used to avoid material losses upon long-term storage.

9. Tagmentation and Indexing PCR (1 h)

9.1 Please note that the Tn5 transposase amount indicated below is a suggested starting point for tagmenting 200 pg/μl cDNA. Optimization might be necessary, depending on the specific activity of each batch of Tn5.

NOTICE

R&D Systems™ now includes the legacy brands Novus Biologicals™, Tocris Bioscience™, and ProteinSimple™.

Bio-Techne Spatial now includes the legacy brands Lunaphore™ and Advanced Cell Diagnostics™.

Bio-Techne Diagnostics now includes Asuragen®, Bionostics™, Cliniqa™, RNA Medical®, and R&D Systems™ Clinical Controls.

9.2 Prepare the Tagmentation Mix as described below:

Reagent	Final Concentration	Volume (μl)
TAPS-Mg Buffer, pH=7.3 (5x)	10 mM TAPS, 5 mM MgCl ₂	0.800
Dimethylformamide (DMF) (100%)	20%	0.800
Tn5 transposase (2 μM working dil.)	62.5 nmol	0.125
Nuclease-free water	-	2.275
Total volume (μl)		3.000

IMPORTANT! Dimethylformamide (DMF) is toxic and should be handled under the hood, or according to local safety regulations.

9.3 Dispense 3 μl of Tagmentation Mix in a new 384-well plate.

9.4 Add 1 μl of normalized cDNA (200 pg/μl) to each well containing the Tagmentation Mix.

9.5 Seal the plate, vortex, spin down, and carry out the tagmentation reaction: 55 °C for 8 min, 4 °C hold. Upon completion proceed immediately to the next step.

9.6 Add 1 μl of 0.2% SDS to each well. Seal the plate, vortex, spin down and incubate for 5 min at room temperature. Do not put the plate back on ice.

9.7 Add 2 μl of prediluted N7xx + S5xx Index Adaptors (5 μM each).

9.8 Add 3 μl of Enrichment PCR Mix to each well:

Reagent	Final Concentration	Volume (μl)
KAPA HiFi enzyme (1 U/μl)	0.02 U/μl	0.200
KAPA HiFi Buffer (5x)	1x	2.000
dNTPs (10 mM)	300 mM	0.300
Nuclease-free water	-	0.500
Total volume (μl)		3.000

9.9 Seal the plate, vortex, spin down, and place it in a thermocycler and carry out the Enrichment PCR Reaction. Adjust the number of PCR cycles according to the number of processed cells.

Step		Temperature	Time	Cycles
Gap filling		72° C	3 min	1x
Enrichment PCR	Initial denaturation	98° C	30 sec	1x
	Denaturation	98° C	10 sec	12x
	Annealing	55° C	30 sec	
	Elongation	72° C	30 sec	
		15° C	hold	

SAFE STOPPING POINT - The final unpurified sequencing library can be stored for several months at -20 °C.

NOTICE

R&D Systems™ now includes the legacy brands Novus Biologicals™, Tocris Bioscience™, and ProteinSimple™.

Bio-Techne Spatial now includes the legacy brands Lunaphore™ and Advanced Cell Diagnostics™.

Bio-Techne Diagnostics now includes Asuragen®, Bionostics™, Cliniqa™, RNA Medical®, and R&D Systems™ Clinical Controls.

10. Library Cleanup and Quantification (45 min.)

- 10.1 Take an aliquot from each sample for the final library cleanup (i.e., 5 µl) and transfer it to a 1.5 ml Eppendorf tube. The rest of the library can be stored long-term at -20°C.
- 10.2 Remove the Sera-Mag SpeedBeads™ working solution (alternatively: AMPure XP beads or SPRI beads) from the +4 °C storage and equilibrate it at room temperature for 15 min.
- 10.3 Add Sera-Mag SpeedBeads™ working solution to a final ratio of 0.8× and mix well to homogenization. Use a different ratio if the goal is to recover longer (i.e., lower ratio) or shorter (i.e., higher ratio) fragments.
- 10.4 Incubate the tube off the magnetic stand for 5 min at room temperature.
- 10.5 Place the tube on the magnetic stand and leave it for 5 min or until the solution appears clear.
- 10.6 Remove the supernatant without disturbing the beads.
- 10.7 Recommended: wash the pellet with 1 ml of 80% v/v ethanol. Incubate for 30 sec without removing the tube from the magnetic stand.
- 10.8 Remove any trace of ethanol and let the bead pellet dry for 2 min. Do not cap the tube or remove it from the magnetic stand during this time. Do not completely dry the beads.
- 10.9 Remove the tube from the magnetic stand, add 50 µl of nuclease-free water and mix well by pipetting or vortexing to resuspend the beads.
- 10.10 Incubate for 2 min off the magnetic stand.
- 10.11 Place the tube back on the magnetic stand and incubate for 2 min or until the solution appears clear.
- 10.12 Remove 49 µl of the supernatant and transfer it to a new 1.5-ml LoBind tube. Store the cDNA in a -20 °C freezer long-term or until ready for sequencing.
- 10.13 Use a Qubit fluorometer or a similar fluorimetric assay to quantify the library. Library yield can vary from 1 to 100 ng/µl depending on the number of cells being pooled as well as PCR cycles used.
- 10.14 Check the final library size on the Agilent Bioanalyzer, or similar equipment. Follow the instructions as described in the High Sensitivity DNA chip user manual.
- 10.15 Use the average size indicated on the Bioanalyzer and the concentration reported after Qubit measurement to determine the exact molarity required for sequencing. Example of trace shown in Figure 2.

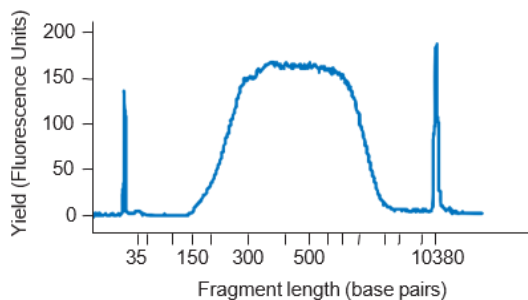


Figure 2. Example of trace of amplified library from a single HEK293 cell (19 cycles). Data obtained from Picelli *et al.* 2025.³

SAFE STOPPING POINT – The final purified sequencing library can be stored for several months at -20°C.

NOTICE

R&D Systems™ now includes the legacy brands Novus Biologicals™, Tocris Bioscience™, and ProteinSimple™.

Bio-Techne Spatial now includes the legacy brands Lunaphore™ and Advanced Cell Diagnostics™.

Bio-Techne Diagnostics now includes Asuragen®, Bionostics™, Cliniqa™, RNA Medical®, and R&D Systems™ Clinical Controls.

11. Pooling and Sequencing

- 11.1 The purified library can be sequenced on any Illumina, Element Bio, or equivalent NGS sequencer. Follow the specifications reported for each instrument.

References

1. **Khven** *et al.* (2025) [SEQURNA enhances FLASH-seq gene detection while eliminating DTT dependence](#). bioRxiv.12.12.693841; *Benchmarking study demonstrating superior performance of SEQURNA Thermostable RNase inhibitor in FLASH-seq.*
2. **Hahaut** *et al.* (2022) [Fast and highly sensitive full-length single-cell RNA sequencing using FLASH-seq](#). Nat Biotechnol **40**, 1447. *Original FLASH-seq publication.*
3. **Picelli S.** (2025) [FLASH-seq with SEQURNA RNase Inhibitor](#). Protocols.io. *For FLASH-seq with SEQURNA, maintained by the Picelli lab.*
4. **Noble** *et al.* (2024). [Introducing synthetic thermostable RNase inhibitors to single-cell RNA-seq](#). Nat Commun **15**, 8373 *Original publication of the SEQURNA Thermostable RNase inhibitor.*
5. **Picelli S.** (2019) [Full-Length Single-Cell RNA Sequencing with Smart-seq2](#). Methods Mol Biol. **1979**, 25. *Protocol including description of in-house preparation of SeraMag beads.*

CONTACT US customerservice.rnd@bio-techne.com / support.rnd@bio-techne.com

NOTICE

R&D Systems™ now includes the legacy brands Novus Biologicals™, Tocris Bioscience™, and ProteinSimple™.

Bio-Techne Spatial now includes the legacy brands Lunaphore™ and Advanced Cell Diagnostics™.

Bio-Techne Diagnostics now includes Asuragen®, Bionostics™, Cliniqa™, RNA Medical®, and R&D Systems™ Clinical Controls.