

## USER GUIDE

# Universal Plus™ mRNA-Seq with NuQuant®

**REF** 0520, 0520B, 0520C, 0520D, 0521, 0521B, 0521C, 0521D, 0522,  
0522B, 0522C, 0522D

**Publication Number: M01485**

**Revision: v7**



## Limited License

Buyer acknowledges and agrees that the sale by Tecan Genomics Inc. (“Tecan Genomics”), and the purchase and/or use by Buyer, of a Tecan Genomics product (“Product”) is subject to all applicable limited licenses, end user license agreements, notices, terms, conditions and/or use restrictions listed (by full text, URL or hyperlink) (i) on any packaging, label or insert associated with the Product, (ii) in any instructions, warnings, user manuals or other documentation accompanying, or provided by Tecan Genomics regarding or for use in connection with the Product or related service (“Documentation”), or (iii) on [www.tecan.com](http://www.tecan.com).

## Regulatory Compliance

Buyer acknowledges that the Product does not have United States Food and Drug Administration (“FDA”) or equivalent non-U.S. regulatory agency approval (“Approval”). Accordingly, Buyer acknowledges that use of the Product will be limited to RESEARCH USE ONLY and is NOT FOR USE IN DIAGNOSTIC PROCEDURES, THERAPEUTIC PROCEDURES OR COMMERCIAL SERVICES. Buyer also acknowledges that the Product is not to be administered to humans or animals. The Product should be used by qualified professionals in strict accordance with the Documentation. Unless otherwise expressly stated by Tecan Genomics in writing, no claim or representation is made or intended by Tecan Genomics (i) as to any diagnostic or clinical use of the Product; (ii) that the Product has any Approval for use in any diagnostic or other procedure, or for any other use requiring compliance with any law, regulation or governmental policy concerning medical devices, laboratory tests or the like (collectively, “Regulatory Laws”); (iii) that the Product will satisfy the requirements of the FDA or any other regulatory agency; or (iv) that the Product or its performance is suitable or has been validated for clinical or diagnostic use, for safety and effectiveness, for any specific use or application, or for importation into Buyer’s jurisdiction. Buyer agrees that if it elects to use the Product for a purpose that would subject Buyer, its customers or the Product to the application of Regulatory Laws or any other law, regulation or governmental policy, Buyer shall be solely responsible for obtaining any required Approvals and otherwise ensuring that the importation of the Product into Buyer’s jurisdiction and Buyer’s use of the Product complies with all such laws, regulations and policies.

## Patents, Licensing and Trademarks

© Tecan Genomics, Inc. All rights reserved. The Tecan Genomics family of products and methods of their use are covered by several issued U.S. and International patents and pending applications ([www.tecan.com/patents](http://www.tecan.com/patents)). Trademarks and registered trademarks of Tecan Genomics are listed on [www.tecan.com/trademarks](http://www.tecan.com/trademarks). Other marks appearing in Documentation are marks of their respective owners.

The purchase of the Product conveys to the Buyer the limited, non-exclusive, non-transferable right (without the right to modify, reverse engineer, resell, repackage or further sublicense) under these patent applications and any patents issuing from these patent applications to use the Product and associated methods SOLELY FOR RESEARCH PURPOSES and in strict accordance with the intended use described in the Documentation. No license to make or sell products by use of the Product is granted to the Buyer whether expressly, by implication, by estoppels or otherwise. In particular, the purchase of the Product does not include or carry any right or license to use, develop or otherwise exploit the Product commercially, and no rights are conveyed to the Buyer to use the Product or its components for purposes including commercial services or clinical diagnostic procedures or therapeutic procedures.

For information on purchasing a license to use this Product and associated Tecan Genomics patents for uses other than RESEARCH PURPOSES, please contact Tecan Global Intellectual Property ([intellectual.property@tecan.com](mailto:intellectual.property@tecan.com)).

## Warranty

EXCEPT AS EXPRESSLY STATED HEREIN, TECAN GENOMICS DISCLAIMS, ALL REMEDIES AND WARRANTIES, EXPRESS, STATUTORY, IMPLIED, OR OTHERWISE, INCLUDING, BUT NOT LIMITED TO, ANY WARRANTIES OF MERCHANTABILITY, SATISFACTORY QUALITY, NONINFRINGEMENT, OR FITNESS FOR A PARTICULAR PURPOSE, OR REGARDING RESULTS OBTAINED THROUGH THE USE OF THE PRODUCT OR ITS COMPONENTS (INCLUDING, WITHOUT LIMITATION, ANY CLAIM OF INACCURATE, INVALID OR INCOMPLETE RESULTS), Tecan Genomics hereby warrants that the Product meets the performance standards described in the Company’s product and technical literature for a period of six months from the date of shipment, provided that the Product is handled and stored according to published instructions, and that the Product is not altered or misused. If the Product fails to meet these performance standards, Tecan Genomics will replace the Product free of charge or issue a credit for the purchase price. Tecan Genomics’ liability under this warranty shall not exceed the purchase price of the Product. TECAN GENOMICS SHALL ASSUME NO LIABILITY FOR DIRECT, INDIRECT, CONSEQUENTIAL OR INCIDENTAL DAMAGES ARISING











FROM THE USE, RESULTS OF USE OR INABILITY TO USE THE PRODUCT. Tecan Genomics reserves the right to change, alter or modify the Product to enhance its performance and design. Except as expressly set forth herein, no right to modify, reverse engineer, distribute, offer to sell or sell the Product is conveyed or implied by Buyer's purchase of the Product.

### Indemnification of Tecan Genomics

Buyer agrees to defend Tecan Genomics against any claim that arises in connection with Buyer's use of the Product (a) that is not for RESEARCH purposes, or (b) that diverges from the intended use described in the Documentation, or (c) that is in violation of any applicable law or regulation. If Buyer is a U.S. state, city, town or other municipality, or a public university, college or other not-for-profit institution chartered under the laws of a U.S. state, this section shall apply to the maximum extent permitted by applicable law. If Buyer is an agency of the U.S. Government; this section shall apply to the maximum extent permitted by the Federal Tort Claims Act, 28 USC 2671, et seq.

### Miscellaneous

This document will be interpreted in accordance with the laws of the State of New York without giving regards to conflict of law principles requiring the application of other laws, and in no event by the United Nations Convention on Contracts for the International Sale of Goods. The parties hereby consent to personal jurisdiction in the state and federal courts in New York in any suit arising out of this Confirmation. The invalidity in whole or in part of any provision hereof shall not affect the validity of any other provision. In case of difference between the terms and conditions in this document and those contained in any Tecan Genomics Terms and Conditions or those under the heading "Limited License" above, the ones in the present document shall prevail.

SYMBOL	MEANING
	Catalog Number
	Consult instructions for use
	Contains sufficient for <n> tests
	Warning
	Important
	Expiration Date
	Temperature limitation
	Optional stopping point
	Note
	Manufacturer

# Contents

<b>I. Introduction .....</b>	<b>1</b>
A. Overview .....	1
B. Storage and Stability .....	1
C. Warnings and Precautions .....	2
D. Before You Start.....	2
<b>II. Components.....</b>	<b>3</b>
A. Reagents Provided.....	3
B. Additional Equipment, Reagents and Labware .....	5
<b>III. Planning the Experiment .....</b>	<b>7</b>
A. Workflow and Time Required .....	7
B. Input RNA Requirements .....	8
C. Programming the Thermal Cycler.....	8
D. Working with the 24- and 96-Plex Adaptor Plates .....	10
E. Bead Purifications .....	10
F. NuQuant .....	11
G. Sequencing Recommendations and Guidelines.....	12
H. Data Analysis.....	13
I. Amplified Library Storage.....	13
<b>IV. Protocol .....</b>	<b>14</b>
Protocol Notes.....	14
A. Sample Preparation .....	15
B. Poly(A) Selection .....	15
C. RNA Elution .....	16
D. RNA Fragmentation .....	17
E. First Strand cDNA Synthesis.....	18
F. Second Strand cDNA Synthesis.....	18
G. cDNA Purification.....	19
H. End Repair.....	20
I. Adaptor Ligation .....	20
J. Strand Selection.....	21
K. Strand Selection Purification .....	22
L. Library Amplification A.....	23
M. Probe Binding (optional AnyDeplete workflow).....	24
N. Targeted Depletion (optional AnyDeplete workflow).....	24
O. Library Amplification B (optional AnyDeplete workflow).....	25
P. Amplified Library Purification .....	26
Q. Quantitative and Qualitative Assessment of the Library.....	27
<b>V. Technical Support .....</b>	<b>29</b>
<b>VI. Appendix .....</b>	<b>30</b>
A. Barcode Sequences and Guidelines for Multiplex Experiments .....	30
B. Library Amplification qPCR Optimization for Standard Workflow .....	31
C. Library Amplification qPCR Optimization for AnyDeplete Workflow .....	33
D. Frequently Asked Questions (FAQs).....	34

# I. Introduction

## A. Overview

### Intended Use

Universal Plus mRNA-Seq with NuQuant is an end-to-end solution for generation of stranded RNA-Seq libraries derived from poly(A) selected RNA. Universal Plus mRNA-Seq with NuQuant is intended for Research Use Only and not for use in diagnostic procedures.

### Features

Universal Plus mRNA-Seq with NuQuant is a streamlined solution for construction of mRNA-Seq libraries from total RNA. This kit is compatible with high quality total RNA obtained from a broad range of tissues or cell lines. This product includes all the reagents necessary for library construction, including pre-plated Unique Dual Index (UDI) adaptors with **DimerFree**<sup>®</sup> technology for efficient library construction without adaptor dimers. Tecan's novel library quantitation method, **NuQuant**<sup>®</sup>, is also integrated providing time- and cost-savings when measuring molarity for library pooling prior to multiplex sequencing. Finally, optional integration of **AnyDeplete**<sup>®</sup> is available with this kit, enabling an integrated solution for depletion of unwanted transcript sequences, providing more efficient use of sequencing resources.

### Specifications

<b>Input type:</b>	High-quality total RNA
<b>Input amount:</b>	10 ng - 1 µg
<b># Reactions available:</b>	24, 96
<b>Barcodes available:</b>	Up to 384 UDI
<b>Sequencing platforms:</b>	Illumina NGS

## B. Storage and Stability

Universal Plus mRNA-Seq is provided as two shipments. The core library preparation kit and optional AnyDeplete module are shipped on dry ice while the mRNA Selection Module is shipped at 4 °C. Both packages should be unpacked and inspected immediately upon receipt.



**Important:** This product contains components with multiple storage temperature requirements. All shipments should be unpacked immediately upon receipt and stored as directed below.



- The mRNA Selection Module (Part No. 0408) containing the Oligo(dT) Beads, mRNA Elution Buffer, mRNA Wash Buffer, and mRNA Binding Buffer should be stored at 2–8 °C.



- K562 Control RNA (RNA<sub>VER</sub> 1) should be stored at –80 °C (provided only in 24 reaction kits).



- All remaining components should be stored at –20 °C in a freezer without a defrost cycle.




This product has been tested to perform to specifications after as many as six freeze/thaw cycles. Kits handled and stored according to the above guidelines are warranted to perform to specifications for 6 months from the date of shipment. Expiration dates listed on the kit label may be longer than the warranty period. Do not use kits that have passed the expiration date.

# I. Introduction

## C. Warnings and Precautions

1. Unpack and inspect the kits immediately upon receiving. In case of severe kit package damage, no dry ice left in the package or ice pack melted, and/or missing components, please contact Tecan NGS Technical Support immediately. Please provide Tecan Genomics with the kit(s) and/or component(s) part number, and lot number. Do not use damaged components.
2. Follow your institution's safety procedures for working with chemicals and handling of biological samples. Follow good laboratory practices and safety guidelines. Wear lab coats, disposable latex gloves and protective glasses where necessary. Changing gloves between handling samples is recommended to avoid contamination of sample or reagents.
3. Consult your institution's environmental waste personnel on proper disposal of unused reagents. Check state and local regulations as they may differ from federal disposal regulations. This material may exhibit characteristics of hazardous waste requiring specific disposal requirements. Institutions should check their country hazardous waste disposal requirements.
4. If appropriate, an SDS for this product is available on the Tecan Genomics website at [www.nugen.com/products/universal-plus-mrna-seq-library-preparation-nuquant](http://www.nugen.com/products/universal-plus-mrna-seq-library-preparation-nuquant)

## D. Before You Start

 Please review this User Guide before using this kit for the first time, including the “Kit Components”, “Planning the Experiment”, “Overview”, “Protocol” and “FAQ” sections. For more information, visit the Universal Plus mRNA-Seq page at [www.nugen.com/products/universal-plus-mrna-seq-library-preparation-nuquant](http://www.nugen.com/products/universal-plus-mrna-seq-library-preparation-nuquant).

New to NGS? Contact Tecan NGS Technical Support at [techserv-gn@tecan.com](mailto:techserv-gn@tecan.com) for tips and tricks on getting started.

## II. Components

### A. Reagents Provided



This kit is provided in a 24 (-24) or 96 (-A01) reaction format. A01 fill sizes may be used in manual or automation workflows.

Universal Plus mRNA-Seq with NuQuant (Part Nos. 0520, 0520B, 0520C and 0520D) is a bundle of:

- Part No. 0361, Universal Plus RNA-Seq with NuQuant
- Part No. 0408, mRNA Selection Module
- Part No. S02480 (0520, 0521, 0522), S02690 (0520B, 0521B, 0522B), 30185200 (0520C, 0521C, 0522C), or 30185201 (0520D, 0521D, 0522D) Unique Dual Index adaptor plate

Universal Plus mRNA-Seq with NuQuant and AnyDeplete (Part Nos. 0521, 0521B, 0521C, 0521D, 0522, 0522B, 0522C and 0522D) also includes:

- Part No. 0359, AnyDeplete Module
- AnyDeplete Probe Mix

**Table 1. Universal Plus RNA-Seq with NuQuant (Part No. 0361)**

COMPONENT	0361-24 PART NUMBER	0361-A01 PART NUMBER	VIAL LABEL	VIAL NUMBER
2X Fragmentation Buffer	S02621	S02423	Clear	FB1 VER 1
Actinomycin D	S02350	S02379	Brown	—
First Strand Buffer Mix	S02612	S02415	Blue	A2 VER 14
First Strand Enzyme Mix	S02613	S02416	Blue	A3 VER 6
Second Strand Buffer Mix	S02333	S02380 (2)	Yellow	B1 VER 11
Second Strand Enzyme Mix	S02614	S02417	Yellow	B2 VER 4
End Repair Buffer Mix	S01708	S02418	Blue	ER1 VER 7
End Repair Enzyme Mix	S01706	S02419	Blue	ER2 VER 4
End Repair Enhancer	S02615	S02420	Blue	ER3 VER 2
Ligation Buffer Mix	S01847	S01689	Yellow	L1 VER 4
Ligation Enzyme Mix	S02616	S02421	Yellow	L3 VER 4
Strand Selection Buffer Mix I	S02617	S02393 (2)	Purple	SS1
Strand Selection Enzyme Mix I	S02618	S02385	Purple	SS2
Strand Selection Enzyme Mix II	S02619	S02386	Purple	SS4
Amplification Reagent I	S02620	S02627	Red	AR1 VER 1
Amplification Reagent II	S02607	S02610	Red	AR2 VER 1

## II. Components

**Table 1 Universal Plus RNA-Seq with NuQuant (Part No. 0361), continued**

COMPONENT	0361-24 PART NUMBER	0361-A01 PART NUMBER	VIAL LABEL	VIAL NUMBER
Amplification Enzyme Mix	S02628	S02629	Red	AR3 VER 1
K562 Control RNA	S02289	—	Clear	RNA VER 1
Nuclease-free Water	S01113	—	Green	D1
Universal Plus UDI 24-Plex Adaptor Plate	S02622	—	—	—
Universal Plus UDI 96-Plex Adaptor Plate (Part No. 0520, 0521, 0522 for A01- fill only)	—	S02480	—	—
Universal Plus UDI-B 96-Plex Adaptor Plate (Part No. 0520B, 0521B and 0522B for A01-fill only)	—	S02690	—	—
Universal Plus UDI-C 96-Plex Adaptor Plate (Part No. 0520C, 0521C and 0522C for A01-fill only)	—	30185200	—	—
Universal Plus UDI-D 96-Plex Adaptor Plate (Part No. 0520D, 0521D and 0522D for A01-fill only)	—	30185201	—	—
NuQuant Buffer	S02516	S02517	Clear	—
NuQuant Standard	S02512	S02512	Clear	—

**Table 2. mRNA Selection Module (Part No. 0408)**

COMPONENT	0408-24 PART NUMBER	0408-A01 PART NUMBER	VIAL LABEL	VIAL NUMBER
Oligo dT Beads	S02623	S02248	Clear	—
mRNA Elution Buffer	S02624	S02424	Clear	—
mRNA Wash Buffer	S02625	S02425	Clear	—
mRNA Binding Buffer	S02626	S02426	Clear	—



## II. Components

**Table 3. AnyDeplete Module (Part No. 0359)**

COMPONENT	0359-24 PART NUMBER	0359-A01 PART NUMBER	VIAL LABEL	VIAL NUMBER
AnyDeplete Buffer Mix	S02594	S02427	Purple	AD1 VER 1
AnyDeplete Enzyme Mix I	S02595	S02428	Purple	AD2 VER 1
AnyDeplete Enzyme Mix II	S02596	S02392	Purple	AD3 VER 1

**Table 4. AnyDeplete Probe Mix**

COMPONENT	24 REACTIONS	96 REACTIONS	VIAL LABEL	VIAL NUMBER
Universal Plus AnyDeplete Probe Mix – Human Globin	S02364	S02431	Purple	AD14 VER 1

To use Universal Plus mRNA-Seq with NuQuant with custom AnyDeplete probe mixes, contact your local Account Executive or Tecan NGS Technical Support (techserv-gn@tecan.com).



The reagents in Universal Plus mRNA-Seq with NuQuant are similar to reagents in our other kits; however, unless the component part numbers are identical, these reagents do not have exactly the same composition and, therefore, are not interchangeable. Do not exchange or replace one reagent named, for example, A1 with another A1, as it will adversely affect performance.

### B. Additional Equipment, Reagents and Labware

#### Required Materials

- **Equipment**
  - Agilent 2100 Bioanalyzer or 2200 TapeStation Instrument, or other equipment for electrophoretic analysis of nucleic acids
  - Microcentrifuge for individual 1.5 mL and 0.5 mL tubes
  - Microcentrifuge for 0.2 mL tube strips or plates
  - 0.5-10  $\mu$ L pipette, 2-20  $\mu$ L pipette, 20-200  $\mu$ L pipette, 200-1000  $\mu$ L pipette
  - 2-20  $\mu$ L or 5-50  $\mu$ L multichannel pipette and 20-200  $\mu$ L or 20-300  $\mu$ L multichannel pipette for sample mixing
  - Vortexer
  - Thermal cycler with 0.2 mL tube heat block, heated lid, and 100  $\mu$ L reaction capacity
  - Qubit<sup>®</sup> 2.0, 3.0 or 4 Fluorometer (Thermo Fisher Scientific) or appropriate fluorometer and accessories for quantification of fragmented DNA and amplified libraries.

## II. Components

- **Reagents**
  - Agencourt RNAClean XP Beads or AMPure XP Beads (Beckman Coulter, Cat. #A63987 or A63881)
  - Ethanol, Absolute (200 Proof), Molecular Biology Grade (Fisher Scientific Cat. #BP2818), for purification steps
  - Low-EDTA TE Buffer, 1X, pH 8.0 (Fisher Scientific, Cat. #75793), for diluting nucleic acids
  - Nuclease-free water (Fisher Scientific, Cat. #BP2484), for diluting nucleic acids
  - Agilent High Sensitivity DNA Kit (Agilent, Cat. #5067-4626) or equivalent
  - EvaGreen® Dye, 20X in water (Biotium, Cat. #31000) optional; for optimizing Library Amplification with qPCR
- **Supplies and Labware**
  - Barrier (filter) pipette tips, nuclease-free
  - 1.5 mL and 0.5 mL RNase-free microcentrifuge tubes
  - 0.2 mL PCR strip tubes or 0.2 mL thin-wall PCR plates
  - Low-retention microcentrifuge tubes (DNA LoBind Tubes, Eppendorf Cat. # 0030108035 or 0030108051)
  - Magnetic stand for 0.2 mL strip tubes or plates. (Beckman Coulter Cat. #A29164 or A32782; Thermo Fisher Scientific Cat. #12331D, 12027, or 12332D; Promega Cat. #V8351). Other magnetic stands may be used as well, although their performance has not been validated by Tecan.
  - Cleaning solutions such as RNaseZap® RNase Decontamination Solution (Thermo Fisher Scientific, Cat. #AM9780) and DNA OFF™ (MP Biomedicals, Cat. #11QD0500)
  - AlumaSeal II film (Sigma-Aldrich Cat. #Z721530)
  - Disposable gloves
  - Kimwipes
  - Ice bucket

### To Order:

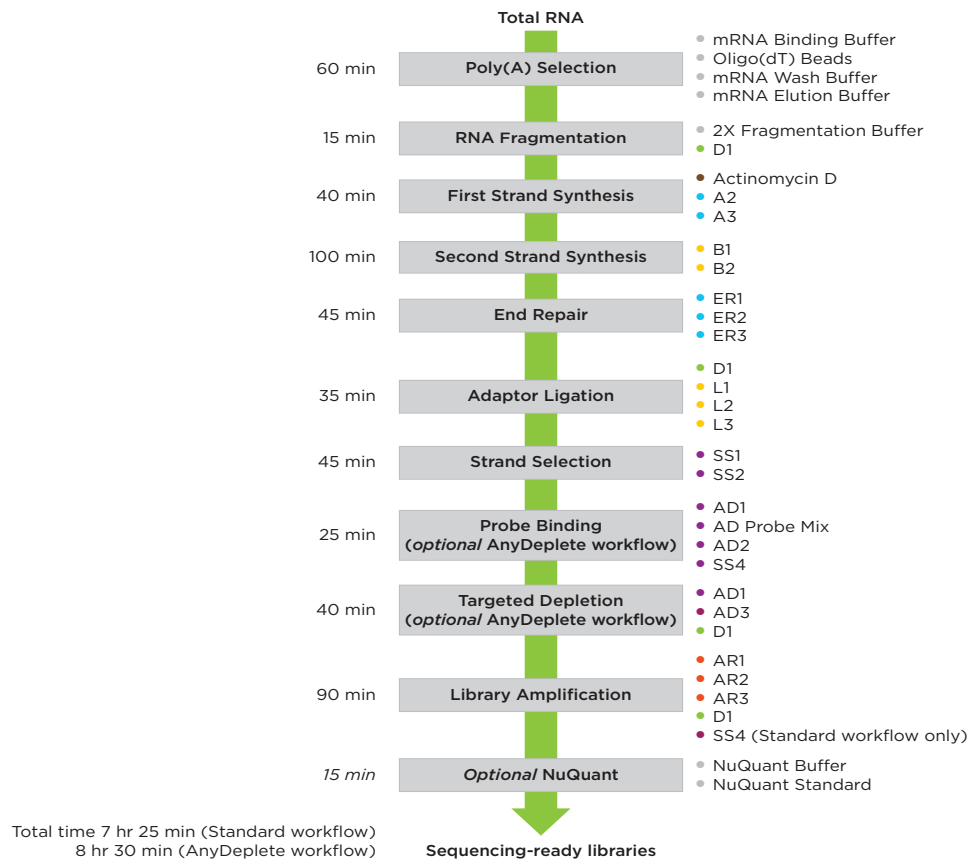
- Agilent, [www.agilent.com](http://www.agilent.com)
- Beckman Coulter, [www.beckmancoulter.com](http://www.beckmancoulter.com)
- Biotium, [www.biotium.com](http://www.biotium.com)
- Eppendorf, [www.eppendorf.com](http://www.eppendorf.com)
- Fisher Scientific, [www.fishersci.com](http://www.fishersci.com)
- MP Biomedicals, [www.mpbio.com](http://www.mpbio.com)
- Qiagen, [www.qiagen.com](http://www.qiagen.com)
- Promega, [www.promega.com](http://www.promega.com)
- Sigma-Aldrich, Inc., [www.sigmaaldrich.com](http://www.sigmaaldrich.com)
- Thermo Fisher Scientific, [www.thermofisher.com](http://www.thermofisher.com)

### III. Planning the Experiment

#### A. Workflow and Time Required

The Universal Plus mRNA-Seq with NuQuant consists of poly(A) RNA selection, RNA fragmentation and double-stranded cDNA generation using a mixture of random and oligo(dT) priming. The library is then constructed by end repairing the cDNA to generate blunt ends, ligation of UDI adaptors, strand selection, and PCR amplification (Figure 1). An optional AnyDeplete step may be performed prior to library amplification to remove unwanted high-abundance transcripts (e.g. globin). The entire workflow can be completed in one day, and yields cDNA libraries ready for either single read or paired-end sequencing on Illumina sequencing platforms.

**Figure 1. Universal Plus mRNA-Seq with NuQuant Workflow.**



### III. Planning the Experiment

#### B. Input RNA Requirements

##### RNA Quantity

Total RNA input must be between 10 ng – 1 µg. We strongly recommend accurate quantification of total RNA to ensure the minimum input requirement is met.

##### RNA Purity

RNA samples must be free of contaminating proteins and other cellular material, organic solvents (including phenol and ethanol) and salts used in many RNA isolation methods. If using an RNA isolation method based on organic solvents, such as TRIzol, we recommend column purification after isolation.

One measure of RNA purity is the ratio of absorbance readings. The A260:A280 ratio for RNA samples should be greater than 1.8 and A260:A230 should be greater than 2.0.

##### RNA Integrity

The Universal Plus mRNA-Seq with NuQuant kit is designed for use with RNA samples of high molecular weight with little or no evidence of degradation. RNA integrity can be determined using the Agilent 2100 Bioanalyzer and the RNA 6000 Nano LabChip or RNA 6000 Pico LabChip.

#### C. Programming the Thermal Cycler

Use a thermal cycler with a heat block designed for 0.2 mL tubes, equipped with a heated lid, and with a capacity of 100 µL reaction volume. Prepare the programs shown in Table 5 following the operating instructions provided by the manufacturer. For thermal cyclers with an adjustable heated lid, set the lid temperature at 100 °C. For thermal cyclers with a fixed-temperature heated lid use the default settings (typically 100–105 °C).

**Table 5. Thermal Cycler Programming**

mRNA SELECTION		VOLUME
<b>Program 1</b> Poly(A) RNA Binding	65 °C – 5 min, hold at 4 °C	110 µL
<b>Program 2</b> RNA Elution	80 °C – 2 min, hold at 25 °C	50 µL
RNA FRAGMENTATION		VOLUME
<b>Program 3a</b> RNA Fragmentation (200 bp insert size)	94 °C – 8 min, hold at 4 °C	20 µL
<b>Program 3b</b> RNA Fragmentation (300 bp insert size)	86 °C – 8 min, hold at 4 °C	20 µL

### III. Planning the Experiment

**Table 5. Thermal Cycler Programming, continued**

cDNA SYNTHESIS		VOLUME
<b>Program 4</b> First Strand Synthesis	25 °C – 5 min, 42 °C – 15 min, 70 °C – 15 min, hold at 4 °C	25 µL
<b>Program 5</b> Second Strand Synthesis	16 °C – 60 min, hold at 4 °C	75 µL
END REPAIR		VOLUME
<b>Program 6</b> End Repair	25 °C – 30 min, 70 °C – 10 min, hold at 4 °C	15 µL
LIGATION		VOLUME
<b>Program 7</b> Adaptor Ligation	25 °C – 30 min, hold at 4 °C	30 µL
STRAND SELECTION		VOLUME
<b>Program 8</b> Strand Selection	72 °C – 10 min, hold at 4 °C	100 µL
AMPLIFICATION (FOR STANDARD WORKFLOW)		VOLUME
<b>Program 9</b> Library Amplification A	37 °C – 10 min, 95 °C – 2 min, 2x(95 °C – 30 s, 60 °C – 90 s), 15x*(95 °C – 30 s, 65 °C – 90 s), 65 °C – 5 min, hold at 4 °C	100 µL
ANYDEPLETE (OPTIONAL)		VOLUME
<b>Program 10</b> Probe Binding	37 °C – 10 min, 95 °C – 2 min, 50 °C – 30 s, 65 °C – 5 min, hold at 4 °C	25 µL
<b>Program 11</b> Targeted Depletion	60 °C – 30 min, 95 °C – 5 min, hold at 4 °C	50 µL
AMPLIFICATION (FOR OPTIONAL ANYDEPLETE WORKFLOW)		VOLUME
<b>Program 12</b> Library Amplification B	95 °C – 2 min, 2x(95 °C – 30 s, 60 °C – 90 s), 17x*(95 °C – 30 s, 65 °C – 90 s), 65 °C – 5 min, hold at 4 °C	100 µL



**Important:** The number of cycles (\*) used for Library Amplification depends on the starting amount and quality of RNA and should be optimized by qPCR. For more information, contact Tecan NGS Technical Support.

### III. Planning the Experiment

#### D. Working with the 24- and 96-Plex Adaptor Plates

The Adaptor Plate included with the 24 and 96 reaction Universal Plus mRNA-Seq kit contains Unique Dual Index (UDI) adaptor mixes with unique eight-base barcodes. Each well contains sufficient volume for preparation of a single library. The Universal Plus mRNA-Seq Adaptor Plates are sealed with a foil seal designed to provide airtight storage.

Thaw adaptor plate on ice and spin down. Do NOT warm above room temperature. Make sure all adaptor mixes are collected at the bottom of the wells and return the adaptor plate to ice after centrifuging. When pipetting the adaptor mixes, puncture the seal for each well you wish to use with a fresh pipet tip, and transfer the appropriate volume of adaptor into your sample. The remaining wells of the plate should remain sealed for use at a later date. Cover used wells with a new foil seal to prevent any remaining adaptor-containing liquid from contaminating future reactions.

For details regarding barcode sequences, please see **Appendix A** on page 30.

#### E. Bead Purifications

##### Agencourt® Beads

Ampure XP or RNA Clean XP Beads (Agencourt beads) are suitable for use with this kit. There are modifications to the Agencourt beads' standard procedure; therefore, you must follow the protocols outlined in this user guide for the use of these beads.

- Remove beads from 4 °C and leave at room temperature for at least 30 minutes prior to use. Cold beads and buffer will result in reduced recovery.
- Prior to use, ensure beads are fully resuspended by vortexing or inverting and tapping the tube.
- If using strip tubes or partial plates, ensure they are firmly placed on the magnetic plate. The use of individual tubes is not advised as they are not very stable on the magnetic plates.
- It is critical to let the beads separate on the magnet for the full time indicated at each step. Removing the binding buffer before the beads have completely separated will impact yields.
- After the binding step has been completed, take care to minimize bead loss when removing the binding buffer. Loss of beads at this step may impact yields. With the samples placed on the magnetic plate, carefully remove the specified quantity of binding buffer from each sample to avoid disturbing the beads.
- Ensure that the 70% ethanol wash is freshly prepared from fresh ethanol stocks. Lower percent ethanol mixes will reduce recovery.
- It is critical that all residual ethanol be removed prior to elution. Therefore, when removing the final ethanol wash, first remove most of the ethanol, then allow the excess to collect at the bottom of the tube before removing the remaining ethanol. This also reduces the required bead air-drying time.
- After drying the beads, inspect each tube carefully and make certain that all the ethanol has evaporated before proceeding with the amplification step.

### III. Planning the Experiment

#### Preparation of EtOH Wash Solution

Prepare a 70% or 80% EtOH wash solution by combining 100% EtOH and nuclease-free water. Make the ethanol mixes fresh, carefully measuring both the ethanol and water with pipettes. This protocol requires 0.8 mL of 70% EtOH solution per sample.



**Important:**

- It is critical that the ethanol solution in the purification steps be prepared fresh on the same day of the experiment from a recently opened stock container.
- Measure both the ethanol and the water components carefully prior to mixing. Failure to do so can result in a higher than anticipated aqueous content, which may reduce amplification yield.

**Table 6. 70% EtOH Wash Recipe**

EtOH SOLUTION	1X REACTION VOLUME*	100% EtOH	NUCLEASE-FREE WATER
70% EtOH	0.8mL	0.56 mL	0.24 mL

\*A minimum of 10% extra volume should be prepared for each sample.

#### F. NuQuant

NuQuant® is a novel method to accurately measure molar concentrations of NGS libraries without the need for separate fragment size analysis. The library molar concentration can be directly measured using fluorometers or standard plate readers, then pooled and quantified for sequencing.

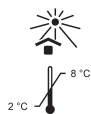
NuQuant is compatible with Qubit 2.0, 3.0 or 4 as well as a wide variety of fluorescent plate readers.

For Qubit-based quantification, an app is required. The apps and installation instructions are available on GitHub: <https://nugentechnologies.github.io/NuQuant/>

#### Preparation of Diluted NuQuant Standard

Universal Plus mRNA-Seq includes a 50X NuQuant Standard stock solution. The fluorescence of this stock corresponds to a 32.2  $\mu$ M library generated with the kit. This stock solution should be diluted prior to use on fluorometers or fluorescent plate readers following the protocol below. The diluted 1X NuQuant Standard corresponds to a 644 nM Universal Plus mRNA-Seq library.

**Note:**



- The 50X and 1X NuQuant Standards should be protected from light.
- NuQuant Standard diluted to 1X for Qubit may be stored at 2–8 °C for up to two months. Do not freeze the 1X NuQuant Standard.

### III. Planning the Experiment

1. Remove concentrated 50X NuQuant Standard stock solution and thaw on ice. Mix by vortexing, spin down and place on ice.
2. Prepare diluted NuQuant Standard in a DNA LoBind tube.
  - A. For use with Qubit, combine 2  $\mu$ L of 50X NuQuant Standard stock solution and 98  $\mu$ L of NuQuant Buffer. Mix thoroughly by vortexing, spin down and store at 2–8 °C.
  - B. For use with fluorescent plate readers, make a fresh dilution of the 50X NuQuant Standard to 1X with Low-EDTA TE Buffer, 1X, pH 8.0. Mix thoroughly by vortexing and spin down.

#### NuQuant Application

NuQuant is compatible with Qubit 2.0, 3.0 or 4 as well as a wide variety of fluorescent plate readers.

For Qubit-based quantification, an app is required. The apps and installation instructions are available on GitHub: <https://nugentechnologies.github.io/NuQuant/>

For other fluorometers, a standard curve must be prepared. For using NuQuant with a fluorescent plate reader, contact Tecan NGS Technical Support at [techserv-gn@tecan.com](mailto:techserv-gn@tecan.com).



**Important:** NuQuant quantitation must be performed prior to storage at –20 °C.

#### G. Sequencing Recommendations and Guidelines

The Universal Plus mRNA-Seq kit produces RNA-Seq libraries compatible with Illumina NGS platforms. These libraries should be sequenced using the Illumina protocol for multiplex sequencing, following the recommendations for the specific sequencer. For sequencing recommendations and questions contact Tecan NGS Technical Support at [techserv-gn@tecan.com](mailto:techserv-gn@tecan.com).

#### Index Read Recommendations

Universal Plus mRNA-Seq with NuQuant libraries contain 8 base Unique Dual Indexes (UDI) for sample multiplexing. Both index 1 (i7) and index 2 (i5) should be sequenced for the detection of “index (barcode) hopping”. These barcodes differ from the sequences used by Illumina and can be found in Appendix A.

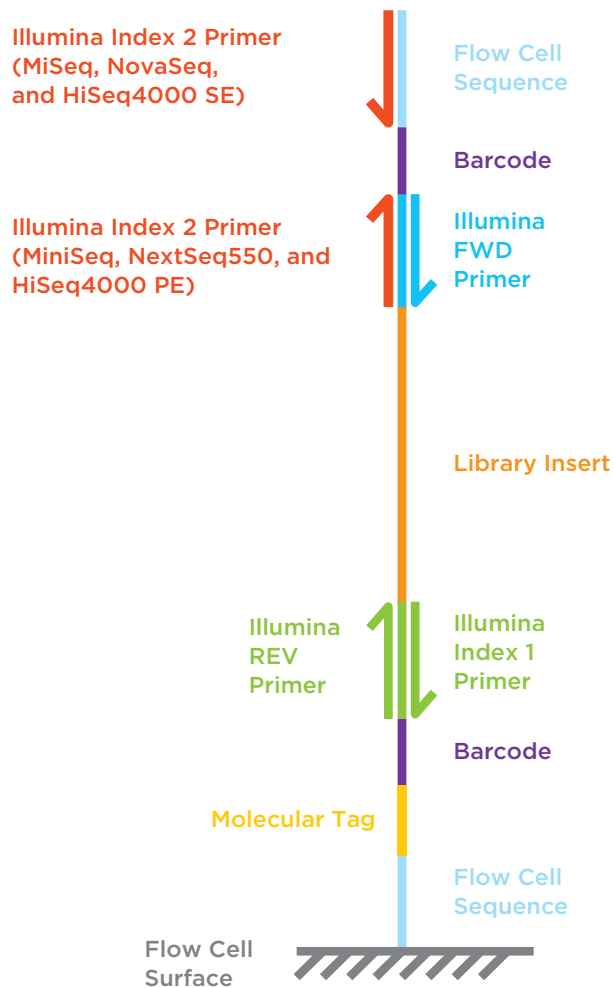
The 24-reaction kit (0520-24, 0521-24 or 0522-24) contains barcodes A01-H03 from the Universal Plus UDI 96-Plex Adaptor Plate (S02480).

The Universal Plus mRNA-Seq with NuQuant kit includes an optional 8 nucleotide (nt) molecular tag which can be used in conjunction with genomic sequence for duplicate read determination. To capture the molecular tag information, use 16 cycles for the index 1 (i7) read.



### III. Planning the Experiment

Figure 2. Universal Plus mRNA-Seq with NuQuant Library Structure.




#### H. Data Analysis

Once the data have been parsed according to sample, additional sample-specific data analysis may be employed according to the requirements of the experiment.



**Note:** The forward read from Universal Plus mRNA-Seq with NuQuant libraries represents the sense strand. This may be opposite to stranded data from other library preparation kits and may require slight modification to the data analysis workflow. Contact Tecan NGS Technical Support for more information.

#### I. Amplified Library Storage

 Amplified libraries may be stored at  $-20^{\circ}\text{C}$ .

### Protocol Notes

#### Controls

- We recommend the routine use of the K562 positive control RNA, provided at 1 µg/µL in the 24 reaction kit, especially the first time a reaction is set up. The use of a positive control RNA will establish a baseline of performance and provide the opportunity to become familiar with the bead purification step. This step may be unfamiliar to many users and can be especially prone to handling variability in using the magnet plate, so a practice run with the plate is highly recommended.
- Routine use of a no template control (NTC) is recommended to monitor the work environment for potential carryover contamination of previous libraries.

#### General Workflow

- Set up no fewer than 8 reactions at a time to ensure that you are not pipetting very small volumes and to ensure sufficient reagent recoveries for the full nominal number of samples from the kit. Making master mixes for fewer than 8 samples at a time may affect reagent recovery volumes.
- Thaw components used in each step and immediately place them on ice. Always keep thawed reagents and reaction tubes on ice unless otherwise instructed. Do not thaw all reagents at once.
- After thawing and mixing buffer mixes, if any precipitate is observed, re-dissolve the precipitate completely prior to use. You may gently warm the buffer mix for 2 minutes at room temperature followed by brief vortexing.
- Keep enzyme mixes on ice after briefly spinning to collect the contents. Do not vortex enzyme mixes nor warm any enzyme, adaptor or primer mixes.
- When preparing master mixes, use the minimal amount of extra material to ensure you are able to run the maximum number of reactions using the components provided in the kit.
- When placing small amounts of reagents into the reaction mix, pipet up and down several times to ensure complete transfer from the pipet tip to the reaction mix.
- When instructed to mix via pipetting, gently aspirate and dispense a volume that is at least half of the total volume of the reaction mix.
- Always allow the thermal cycler to reach the initial incubation temperature prior to placing the tubes or plates in the block.

#### Reagents

- Use the water provided with the kit (green: D1; 24 reaction kit only) or an alternate source of nuclease-free water. We do not recommend the use of DEPC-treated water with this protocol.
- Components and reagents from other Tecan products should not be used with this product.
- Use only fresh ethanol stocks to make ethanol for washes in the purification protocols.
- Make the ethanol mixes fresh, carefully measuring both the ethanol and water with pipettes. Lower concentrations of ethanol in wash solutions will result in loss of yield as the higher aqueous content will dissolve the cDNA and wash it off the beads or column.

## IV. Protocol

This protocol includes workflows for RNA-Seq library construction with and without AnyDeplete.

For **standard workflows**, follow sections **A. Sample Preparation** through **L. Library Amplification A**, then continue to **P. Amplified Library Purification**.

This workflow requires the following components:

Part No. 0408 – mRNA Selection Module

Part No. 0361 – Universal Plus RNA-Seq with NuQuant

For **AnyDeplete workflows**, follow sections **A. Sample Preparation** through **K. Strand Selection Purification**, then continue to **M. Probe Binding (optional AnyDeplete workflow)**.

This workflow requires the following components:

Part No. 0408 – mRNA Selection Module

Part No. 0361 – Universal Plus RNA-Seq with NuQuant

Part No. 0359 – AnyDeplete Module

AnyDeplete Probe Mix

For each section of the protocol, remove reagents from recommended storage conditions listed in “B. Storage and Stability” on page 1. Thaw and place reagents at room temperature or on ice as instructed. After each section, continue immediately to the next section of the protocol unless otherwise directed.

Return all reagents to their appropriate storage conditions promptly after use unless otherwise instructed.

### A. Sample Preparation

1. Remove Nuclease-free water (Green: D1; 24 reaction kit only) from storage and place at room temperature.
2. Aliquot each total RNA input sample (10 ng – 1 µg) into a 0.2 mL strip tube or 96-well plate.
3. Dilute the RNA with D1 to a final volume of 50 µL and place on ice.



**Note:** Keep D1 at room temperature for use in subsequent sections.

### B. Poly(A) Selection

1. Remove mRNA Wash Buffer (Clear) and mRNA Elution Buffer (Clear) from storage and place at room temperature for use in the next section.
2. Remove mRNA Binding Buffer (Clear) and Oligo(dT) Beads (Clear) from storage and place at room temperature at least 10 minutes prior to use. Ensure these reagents have completely reached room temperature before proceeding.

**Table 7. Oligo(dT) Bead Master Mix**

REAGENT	OLIGO(dT) BEADS (CLEAR)	mRNA BINDING BUFFER (CLEAR)
1X REACTION VOLUME	10 $\mu$ L	50 $\mu$ L

3. Prepare Oligo(dT) Bead Master Mix by combining mRNA Binding Buffer and Oligo(dT) Beads in a 0.5 mL capped tube according to the volumes shown in Table 6.
4. Aliquot 60  $\mu$ L of Oligo(dT) Bead Master Mix into a tube for each sample. Add 50  $\mu$ L of total RNA to the Oligo(dT) Bead Master Mix for a total of 110  $\mu$ L. Mix thoroughly by pipetting up and down at least 10 times, taking care when pipetting to minimize foaming.
5. Place the tubes in a pre-warmed thermal cycler programmed to run Program 1 (Poly(A) RNA Binding; see Table 5):  
65 °C – 5 min, hold at 4 °C
6. Remove the tubes from the thermal cycler, spin to collect condensation and place on ice.

### C. RNA Elution

1. Remove 2X Fragmentation Buffer (Clear: FB1) from storage for use in the next section. Thaw at room temperature, mix by vortexing, spin down and place on ice.
2. Mix samples thoroughly by pipetting slowly up and down at least 10 times.
3. Incubate at room temperature for 5 minutes.
4. Repeat steps 2 and 3 for a total of two incubations.
5. Transfer the tubes to the magnet and let stand for 2 minutes to completely clear the solution of beads.
6. Carefully remove the binding buffer and discard it, taking care not to disturb the beads.
7. Remove samples from the magnet.
8. Add 200  $\mu$ L of mRNA Wash Buffer to the beads. Mix thoroughly by slowly pipetting up and down at least 10 times.
9. Transfer the tubes to the magnet and let stand for 2 minutes to completely clear the solution of beads.
10. Carefully remove the wash buffer and discard it, taking care not to disturb the beads.
11. Remove tubes from the magnet.
12. Add 50  $\mu$ L of mRNA Elution Buffer to each sample. Resuspend beads by slowly pipetting up and down at least 10 times.
13. Place the tubes in a pre-warmed thermal cycler programmed to run Program 2 (RNA Elution; see Table 5):  
80 °C – 2 min, hold at 25 °C
14. Remove tubes from the thermal cycler, spin to collect condensation and place on ice.

## IV. Protocol

15. Add 50  $\mu\text{L}$  of mRNA Binding Buffer to bead/elution buffer mix. Mix thoroughly by slowly pipetting up and down at least 10 times.
16. Incubate at room temperature 5 minutes.
17. Transfer the tubes to the magnet and let stand for 2 minutes to completely clear the solution of beads.
18. Carefully remove the buffer mix and discard it, taking care not to disturb the beads.
19. Remove the tubes from the magnet.
20. Add 200  $\mu\text{L}$  of mRNA Wash Buffer to the beads. Mix thoroughly by slowly pipetting up and down at least 10 times.
21. Transfer the tubes to the magnet and let stand for 2 minutes to completely clear the solution of beads.
22. Continue to **D. RNA Fragmentation** with the tubes incubating on the magnet.

### D. RNA Fragmentation

1. Remove Actinomycin D (Brown) from storage for use in the next section. Thaw at room temperature, mix by vortexing, spin down and place at room temperature.

**Table 8. 1X Fragmentation Buffer**

REAGENT	2X FRAGMENTATION BUFFER (CLEAR: FB1 VER 1)	NUCLEASE-FREE WATER (GREEN: D1)
1X REACTION VOLUME	10 $\mu\text{L}$	10 $\mu\text{L}$

2. Prepare 1X Fragmentation Buffer by combining FB1 and D1 in a 0.5 mL capped tube according to the volumes shown in Table 7. Resuspend beads by slowly pipetting up and down at least 10 times.
3. Carefully remove the wash buffer from the samples and discard it, taking care not to disturb the beads.
4. Remove the tubes from the magnet.
5. Add 20  $\mu\text{L}$  of 1X Fragmentation Buffer to each sample. Resuspend beads by slowly pipetting up and down at least 10 times.
6. Place the tubes in a pre-warmed thermal cycler programmed to run Program 3 (RNA Fragmentation; see Table 5).  
  
Program 3A (200 bp insert size)    94 °C – 8 min, hold at 4 °C  
Program 3B (300 bp insert size)    86 °C – 8 min, hold at 4 °C
7. Remove tubes from the thermal cycler, spin to collect condensation and place on ice.
8. Transfer the tubes to the magnet and let stand for at least 2 minutes to completely clear the solution of beads.
9. Transfer 20  $\mu\text{L}$  of fragmented mRNA to a new 0.2 mL tube and place on ice.

### E. First Strand cDNA Synthesis

**Table 9. First Strand Master Mix**

REAGENT	ACTINOMYCIN D (BROWN)	FIRST STRAND BUFFER MIX (BLUE: A2 VER 14)	FIRST STRAND ENZYME MIX (BLUE: A3 VER 6)
1X REACTION VOLUME	1.25 µL	2.75 µL	1 µL

- Spin down the contents of A3 and place on ice.
- Thaw A2 at room temperature. Mix by vortexing, spin down and place on ice.
- Prepare a master mix by combining Actinomycin D, A2 and A3 in a 0.5 mL capped tube according to the volumes shown in Table 8. Mix well by pipetting, spin down, and place on ice.
- Add 5 µL of First Strand Master Mix to each sample tube for a total of 25 µL. Mix well by pipetting, spin down and place on ice.
- Place the tubes in a thermal cycler programmed to run Program 4 (First Strand Synthesis; see Table 5):  
25 °C – 5 min, 42 °C – 15 min, 70 °C – 15 min, hold at 4 °C
- Remove the tubes from the thermal cycler, spin to collect condensation and place on ice.

### F. Second Strand cDNA Synthesis

- Remove Agencourt beads from storage and place on the bench top to reach room temperature for use in the next section.

**Table 10. Second Strand Master Mix**

REAGENT	SECOND STRAND BUFFER MIX (YELLOW: B1 VER 11)	SECOND STRAND ENZYME MIX (YELLOW: B2 VER 4)
1X REACTION VOLUME	48 µL	2 µL

- Spin down the contents of B2 and place on ice.
- Thaw B1 at room temperature. Mix by vortexing, spin down and place on ice.
- Prepare a master mix by combining B1 and B2 in a 0.5 mL capped tube according to the volumes shown in Table 9. Mix well by pipetting, spin down, and place on ice.
- Add 50 µL of Second Strand Master Mix to each sample tube for a total of 75 µL. Mix well by pipetting, spin down and place on ice.
- Place the tubes in a pre-cooled thermal cycler programmed to run Program 5 (Second Strand Synthesis; see Table 5):  
16 °C – 60 min, hold at 4 °C
- Remove the tubes from the thermal cycler and spin to collect condensation.

### G. cDNA Purification

1. Ensure the Agencourt beads and D1 have completely reached room temperature before proceeding.
2. Prepare a 70% ethanol wash solution.

#### Important

- It is critical that the ethanol solution in the purification steps be prepared fresh on the day of the experiment from a recently opened stock container.
- Measure both the ethanol and the water carefully prior to mixing. Failure to do so can result in a higher than anticipated aqueous content, which may reduce sample recovery.

3. Resuspend the beads by inverting and tapping the tube. Ensure that the beads are fully resuspended before adding to the sample. After resuspending, do not spin the beads.
4. At room temperature, add 135  $\mu\text{L}$  (1.8 volumes) of Agencourt beads to each reaction and mix by pipetting 10 times.



**Note:** At this step, the reaction will be 210  $\mu\text{L}$ . Pipet carefully to avoid spilling the sample.

5. Incubate at room temperature for 10 minutes.
6. Transfer the tubes to the magnet and let stand 5 minutes to completely clear the solution of beads.
7. Keeping the tubes on the magnet, carefully remove the binding buffer and discard it.
8. With the tubes still on the magnet, add 200  $\mu\text{L}$  of freshly prepared 70% ethanol and allow to stand for 30 seconds.



**Note:** The beads should not disperse; instead, they will stay on the walls of the tubes. Significant loss of beads at this stage will impact cDNA yields, so ensure beads are not removed with the binding buffer or the washes.

9. Remove the 70% ethanol wash using a pipette.



**Note:** It is critical to remove as much of the ethanol as possible. Use at least 2 pipetting steps and allow excess ethanol to collect at the bottom of the tubes after removing most of the ethanol in the first pipetting step.

10. Air-dry the beads on the magnet for 10 minutes. Inspect each tube carefully to ensure that all the ethanol has evaporated.
11. Remove the tubes from the magnet.
12. Add 11  $\mu\text{L}$  D1 to the dried beads. Mix thoroughly to ensure all beads are resuspended.
13. Incubate at room temperature for 5 minutes.
14. Transfer the tubes to the magnet and let stand for at least 5 minutes to completely clear the solution of beads.
15. Transfer 10  $\mu\text{L}$  of sample to a fresh 0.2 mL tube.



*Optional stopping point: Store samples at  $-20\text{ }^{\circ}\text{C}$ .*

### H. End Repair

**Table 11. End Repair Master Mix**

REAGENT	END REPAIR BUFFER MIX (BLUE: ER1 VER 7)	END REPAIR ENZYME MIX (BLUE: ER2 VER 4)	END REPAIR ENHANCER (BLUE: ER3 VER 2)
1X REACTION VOLUME	4 $\mu$ L	0.5 $\mu$ L	0.5 $\mu$ L

1. Spin down the contents of ER2 and ER3 and place on ice.
2. Thaw ER1 at room temperature. Mix by vortexing, spin down and place on ice.
3. Prepare a master mix by combining ER1, ER2 and ER3 in a 0.5 mL capped tube according to the volumes shown in Table 11. Mix well by pipetting, spin down and place on ice.
4. Add 5  $\mu$ L of End Repair Master Mix to each sample tube for a total of 15  $\mu$ L. Mix well by pipetting, spin down and place on ice.
5. Place the tubes in a thermal cycler programmed to run Program 6 (End Repair; see Table 5):  
25 °C – 30 min, 70 °C – 10 min, hold at 4 °C
6. Remove the tubes from the thermal cycler, spin to collect condensation and place on ice.

### I. Adaptor Ligation

1. Thaw Adaptor Plate on ice, spin down, and return to ice.

**Table 12. Ligation Master Mix**

REAGENT	NUCLEASE-FREE WATER (GREEN: D1)	LIGATION BUFFER MIX (YELLOW: L1 VER 4)	LIGATION ENZYME MIX (YELLOW: L3 VER 4)
1X REACTION VOLUME	4.5 $\mu$ L	6.0 $\mu$ L	1.5 $\mu$ L

2. Spin down L3 and place on ice.
3. Thaw L1 at room temperature. Mix by vortexing, spin down and place on ice.
4. Add the entire 15  $\mu$ L of sample to the appropriate adaptor well, mix well by pipetting, then transfer the entire 18  $\mu$ L of sample to a PCR tube.



**Note:** The L1 Ligation Buffer Mix is very viscous. Pipet this reagent slowly and mix thoroughly.



## IV. Protocol

5. Prepare a master mix by combining D1, L1 and L3 in a 0.5 mL capped tube according to the volumes shown in Table 12 Mix by pipetting slowly, without introducing bubbles, spin down and place on ice. Use the master mix immediately.



**Note:** The Ligation Master Mix is very viscous. Pipet this reagent slowly and mix thoroughly.

6. Add 12  $\mu\text{L}$  of Ligation Master Mix to each sample tube for a total of 30  $\mu\text{L}$ . Mix thoroughly by pipetting slowly and gently, spin down and place on ice. Proceed immediately with the incubation.
7. Place the tubes in a thermal cycler programmed to run Program 7 (Adaptor Ligation; see Table 5):  
25 °C – 30 min, hold at 4 °C
8. Remove the tubes from the thermal cycler, spin to collect condensation and place on ice.

### J. Strand Selection

1. Remove Agencourt beads from storage and place on the bench top to reach room temperature for use in the next section.

**Table 13. Strand Selection Master Mix**

REAGENT	STRAND SELECTION BUFFER MIX I (PURPLE: SS1)	STRAND SELECTION ENZYME MIX I (PURPLE: SS2)
1X REACTION VOLUME	69 $\mu\text{L}$	1 $\mu\text{L}$

2. Thaw SS1 at room temperature. Mix by vortexing, spin down and place on ice.
3. Spin down SS2 and place on ice.
4. Prepare a master mix by combining SS1 and SS2 in a 0.5 mL capped tube according to the volumes shown in Table 13.
5. Add 70  $\mu\text{L}$  of Strand Selection Master Mix to 30  $\mu\text{L}$  of each sample for a total of 100  $\mu\text{L}$ . Mix by pipetting, spin down and place on ice.
6. Place the tubes in a pre-warmed thermal cycler programmed to run Program 8 (Strand Selection; see Table 5):  
72 °C – 10 min, hold at 4 °C
7. Remove the tubes from the thermal cycler and spin to collect condensation.

### K. Strand Selection Purification

1. Ensure the Agencourt beads have completely reached room temperature before proceeding.
2. Resuspend the beads by inverting and tapping the tube. Ensure the beads are fully resuspended before adding to samples. After resuspending, do not spin the beads.
3. Add 80  $\mu\text{L}$  (0.8 volumes) of the bead suspension to the Strand Selection reaction product. Mix thoroughly by pipetting up and down.
4. Incubate at room temperature for 10 minutes.
5. Transfer the tubes to the magnet and let stand 5 minutes to completely clear the solution of beads.
6. Keeping the tubes on the magnet, carefully remove the binding buffer and discard it.



**Note:** The beads should not disperse; instead, they will stay on the walls of the tubes. Significant loss of beads at this stage will impact the amount of DNA carried into subsequent steps of the protocol, so ensure beads are not removed with the binding buffer or the wash.

7. With the tubes still on the magnet, add 200  $\mu\text{L}$  of freshly prepared 70% ethanol and allow to stand for 30 seconds.
8. Remove the 70% ethanol wash using a pipette.



**Note:** It is critical to remove as much of the ethanol as possible. Use at least two pipetting steps and allow excess ethanol to collect at the bottom of the tubes after removing most of the ethanol in the first pipetting step.

9. Air-dry the the beads on the magnet for 10 minutes. Inspect each tube carefully to ensure that all of the ethanol has evaporated. It is critical that all residual ethanol be removed prior to continuing.
10. Remove the tubes from the magnet.
11. Add 16  $\mu\text{L}$  D1 to the dried beads. Mix thoroughly to ensure all beads are resuspended.
12. Transfer the tubes to the magnet and let stand for 3 minutes for the beads to clear the solution.
13. Carefully remove 15  $\mu\text{L}$  of the eluate, ensuring as few beads as possible are carried over, transfer to a fresh set of PCR tubes and place on ice.
14. Continue to **L. Library Amplification A** (for standard workflow) or to **M. Probe Binding** (for optional AnyDeplete workflow).



*Optional stopping point: Store samples at  $-20\text{ }^{\circ}\text{C}$ .*

### L. Library Amplification A



**Note:** The number of cycles used for PCR amplification should be optimized via real-time PCR whenever using a sample for the first time with the kit, or using a new amount of input. See **Appendix B** for a protocol to determine the appropriate number of cycles for your sample.

1. Remove Agencourt beads from storage and place on the bench top to reach room temperature for use in the next section.

**Table 14. Library Amplification A Master Mix**

REAGENT	AMPLIFICATION REAGENT I (RED: AR1 VER 1)	AMPLIFICATION REAGENT II (RED: AR2 VER 1)	AMPLIFICATION ENZYME MIX (RED: AR3 VER 1)	STRAND SELECTION ENZYME MIX II (PURPLE: SS4)	NUCLEASE-FREE WATER (GREEN: D1)
1X REACTION VOLUME	10 µL	8 µL	0.5 µL	1 µL	65.5 µL

2. Thaw AR1 and AR2 at room temperature. Mix by vortexing, spin down and place on ice.
3. Spin down AR3 and SS4 and place on ice.
4. Make a master mix by sequentially combining D1, AR1 and AR2 in an appropriately sized capped tube according to the volumes shown in Table 14. Add AR3 and SS4 at the last moment and mix well by pipetting, taking care to avoid bubbles. Spin down and place on ice.
5. On ice, add 85 µL of Library Amplification A Master Mix to each sample for a total of 100 µL. Mix by pipetting, spin down and place on ice.
6. Place the tubes in a pre-warmed thermal cycler programmed to run Program 9 (Library Amplification A; see Table 5):  
 37 °C – 10 min, 95 °C – 2 min, 2x(95 °C – 30 s, 60 °C – 90 s), 15x\*(95 °C – 30 s, 65 °C – 90 s), 65 °C – 5 min, hold at 4 °C  
 \*The recommended number of PCR cycles given in the example above is based on starting with 100 ng of total RNA. The number of PCR cycles may be decreased or increased based on the requirements for a given sample.
7. Remove the tubes from the thermal cycler and spin to collect condensation.
8. Continue to **P. Amplified Library Purification**.

**M. Probe Binding (optional AnyDeplete workflow)****Table 15. Probe Binding Master Mix**

REAGENT	ANYDEplete BUFFER MIX (PURPLE: AD1 VER 1)	ANYDEplete PROBE MIX (PURPLE)	ANYDEplete ENZYME MIX I (PURPLE: AD2 VER 1)	STRAND SELECTION ENZYME MIX II (PURPLE: SS4)
1X REACTION VOLUME	5 $\mu$ L	4 $\mu$ L	0.5 $\mu$ L	0.5 $\mu$ L

1. Thaw AD1 and AnyDeplete Probe Mix at room temperature. Mix by pipetting, spin down and place on ice.
2. Spin down AD2 and SS4 and place on ice.
3. Prepare a master mix by combining AD1, AnyDeplete Probe Mix, AD2 and SS4 in a 0.5 mL capped tube, according to the volumes shown in Table 15.
4. Add 10  $\mu$ L of Probe Binding Master Mix to each sample for a total of 25  $\mu$ L. Mix by pipetting, spin down and place on ice.
5. Place the tubes in a pre-warmed thermal cycler programmed to run Program 10 (Probe Binding; see Table 5):  
37 °C – 10 min, 95 °C – 2 min, 50 °C – 30 s, 65 °C – 5 min, hold at 4 °C
6. Remove the tubes from the thermal cycler spin to collect condensation and place on ice.

**N. Targeted Depletion (optional AnyDeplete workflow)****Table 16. Targeted Depletion Master Mix**

REAGENT	ANYDEplete BUFFER MIX (PURPLE: AD1 VER 1)	ANYDEplete ENZYME MIX II (PURPLE: AD3 VER 1)	NUCLEASE-FREE WATER (GREEN: D1)
1X REACTION VOLUME	5 $\mu$ L	4 $\mu$ L	16 $\mu$ L

1. Thaw AD1 at room temperature. Mix AD1 by vortexing, spin down and place on ice.
2. Spin down AD3 and place on ice.
3. Prepare a master mix by combining AD1, AD3 and D1 in a 0.5 mL capped tube according the volumes in Table 16. Mix thoroughly by pipetting, spin down and place on ice.
4. Add 25  $\mu$ L of Targeted Depletion Master Mix to each sample for a total of 50  $\mu$ L. Mix by pipetting, spin down and place on ice.

## IV. Protocol

- Place the tubes in a pre-warmed thermal cycler programmed to run Program 11 (Targeted Depletion; see Table 5):  
60 °C – 30 min, 95 °C – 5 min, hold at 4 °C
- Remove the tubes from the thermal cycler, spin to collect condensation and place on ice.

### O. Library Amplification B (optional AnyDeplete workflow)



**Note:** The number of cycles used for PCR amplification should be optimized via real-time PCR whenever using a sample for the first time with the kit, or using a new amount of input. See **Appendix C** for a protocol to determine the appropriate number of cycles for your sample.

- Remove Agencourt beads from storage and place on the bench top to reach room temperature for use in the next section.

**Table 17. Library Amplification B Master Mix**

REAGENT	AMPLIFICATION REAGENT I (RED: AR1 VER 1)	AMPLIFICATION REAGENT II (RED: AR2 VER 1)	AMPLIFICATION ENZYME MIX (RED: AR3 VER 1)	NUCLEASE-FREE WATER (GREEN: D1)
1X REACTION VOLUME	10 µL	8 µL	0.5 µL	31.5 µL

- Thaw AR1 and AR2 at room temperature. Mix by vortexing, spin down and place on ice.
- Spin down AR3 and place on ice.
- Prepare a master mix by sequentially combining D1, AR1 and AR2 in an appropriately sized capped tube according to the volumes shown in Table 17. Add AR3 at the last moment and mix well by pipetting, taking care to avoid bubbles. Spin down and place on ice.
- On ice, add 50 µL of Library Amplification B Master Mix to each sample for a total of 100 µL. Mix by pipetting, spin down and place on ice.
- Place the tubes in a pre-warmed thermal cycler programmed to run Program 12 (Library Amplification B; see Table 5):  
95 °C – 2 min, 2x(95 °C – 30 s, 60 °C – 90 s), 17x\*(95 °C – 30 s, 65 °C – 90 s), 65 °C – 5 min, hold at 4 °C

\*The number of PCR cycles given in the example above is based on starting with a 100 ng of total RNA from whole blood containing approximately 30% globin, and performing globin AnyDeplete. Decrease or increase the cycle number as needed based on the requirements for a given sample.

- Remove the tubes from the thermal cycler and spin to collect condensation.



*Optional stopping point: Store samples at -20 °C.*

### P. Amplified Library Purification

1. Remove NuQuant reagents from storage for use in the next section:
  - Remove diluted NuQuant Standard from storage. Mix thoroughly by vortexing, spin and place at room temperature. Protect from light.
  - Remove NuQuant Buffer from storage and thaw at room temperature. Mix thoroughly by vortexing, spin and place on benchtop.
2. Ensure the Agencourt beads and D1 have completely reached room temperature before proceeding.
3. Resuspend the beads by vortexing. Ensure the beads are fully resuspended before adding to samples. After resuspending, do not spin the beads.
4. At room temperature, add Agencourt beads to 100  $\mu$ L of library amplification product. Mix thoroughly by pipetting, taking care to avoid spilling the sample.<sup>†</sup>

REQUIRED INSERT SIZE	BEAD SUSPENSION AMOUNT
200 bp	100 $\mu$ L
300 bp	70 $\mu$ L

5. Incubate at room temperature for 10 minutes.
6. Transfer the tubes to the magnet and let stand 5 minutes to completely clear the solution of beads.
7. Carefully remove and discard the supernatant, taking care not to disturb the beads.



**Important:** It is critical to remove as much of the supernatant as possible. Use at least two pipetting steps and allow excess buffer to collect at the bottom of the tubes after removing most of the supernatant in the first pipetting step.

8. With the tubes still on the magnet, add 200  $\mu$ L of freshly prepared 70% ethanol and allow to stand for 30 seconds.
9. Remove the 70% ethanol wash using a pipette.
10. Repeat the 70% ethanol wash (steps 8 and 9) one more time, for a total of two washes.



**Important:** It is critical to remove as much of the supernatant as possible. Use at least two pipetting steps and allow excess buffer to collect at the bottom of the tubes after removing most of the supernatant in the first pipetting step.

11. Air dry the beads on the magnet for 10 minutes. Inspect each tube carefully to ensure that all of the ethanol has evaporated. It is critical that all residual ethanol be removed prior to continuing.
12. Remove the tubes from the magnet.
13. Add 31  $\mu$ L D1 to the dried beads. Mix thoroughly to ensure all beads are resuspended.
14. Transfer the tubes to the magnet and let stand for at least 3 minutes for the beads to completely clear the solution.

<sup>†</sup> If tubes have become noticeably warped or plastic integrity compromised, transfer all 200  $\mu$ L of sample to fresh tubes.

- Carefully remove 30  $\mu\text{L}$  of the eluate, ensuring as few beads as possible are carried over; transfer to a fresh set of PCR tubes and place on ice.

### Q. Quantitative and Qualitative Assessment of the Library

#### Library Quantification with NuQuant<sup>†</sup>



**Note:** A detailed protocol for the NuQuant Qubit assay is provided below. For information on using NuQuant with other fluorometers and fluorescent plate readers, please contact Tecan NGS Technical Support.

- Ensure NuQuant Buffer and diluted NuQuant Standard (prepared in section III. F.) have reached room temperature before proceeding.
- Aliquot 200  $\mu\text{L}$  NuQuant Buffer into a new thin-wall, clear, 0.5 mL tube. Label this tube S1 (blank).
- Aliquot 195  $\mu\text{L}$  NuQuant Buffer into thin-wall, clear, 0.5-mL tubes. Prepare one tube for the diluted standard and one tube for each library. Label the diluted standard tube S2.



**Note:** The S2 Standard represents a Universal Plus mRNA-Seq with NuQuant library with a concentration of 16.1 nM.

- Add 5  $\mu\text{L}$  of diluted NuQuant Standard or library to each tube containing NuQuant Buffer from Step 3. Close lids securely, mix thoroughly by vortexing, and spin briefly to collect all liquid into the bottom of the tube.



- The diluted NuQuant Standard may settle over time. Mix thoroughly before aliquoting.
- If standards and libraries are not thoroughly mixed, inaccurate results may occur.

- Measure samples as directed for your specific quantification platform:
  - Qubit 2.0: Select “Univ. Plus” from the Qubit home screen. Follow the on screen prompts to read S1, S2, and samples.



**Note:** For Qubit 2.0, the “Assay Concentration” is displayed after reading each sample. To calculate the stock concentration of each library, select the “Calculate Stock Conc.” button, and change the “Volume of Sample Used” to 5  $\mu\text{L}$ . The correct stock concentration of your library will be displayed.

- Qubit 3.0 and 4: Navigate to the saved location of Tecan NuQuant app and select “Univ. Plus”. Follow the on screen prompts to read S1, S2, and samples.



**Note:** For Qubit 3.0 and 4, enter the original sample volume (5  $\mu\text{L}$ ) after reading standard 2. The result for each sample will be displayed as the “Original Calculated Sample Concentration” and the “Qubit Tube Concentration.”

<sup>†</sup> Note: For quantification of libraries without use of NuQuant, a qPCR-based method is recommended. Contact Tecan NGS Technical Support for more information.

## IV. Protocol

### Library Pooling and Assessment

1. Pool the libraries according to NuQuant concentration. An example using the calculation  $V_1=(C_2*V_2)/C_1$ , where C= concentration and V=volume, is provided in Table 18 below.

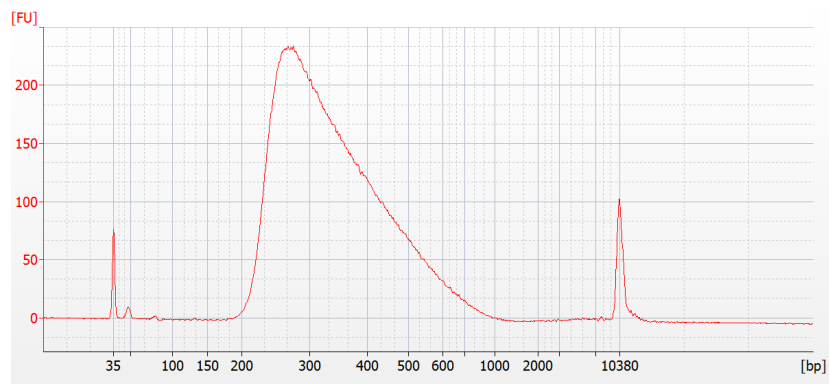
**Table 18. Example Library Pooling Calculation using NuQuant Concentration**

LIBRARY	NUQUANT CONCENTRATION (C <sub>1</sub> )	TARGET CONCENTRATION OF POOL (C <sub>2</sub> )	TARGET VOLUME OF POOL (V <sub>2</sub> )	VOLUME OF LIBRARY TO ADD TO POOL (V <sub>1</sub> )
1	250 nM	10 nM	200 µL	2.0 µL
2	220 nM			2.3 µL
3	280 nM			1.8 µL
4	240 nM			2.1 µL
Volume of low-EDTA or nuclease-free water to add				191.8 µL
Total volume of library pool				200 µL

2. Validate the library pool for sequencer loading following the Illumina guidelines, “Best practices for manually normalizing library concentrations” for your specific sequencer. See **Appendix A** of this guide for guidelines on color balancing and multiplexing of Tecan libraries. A typical fragment distribution for Universal Plus mRNA-Seq with NuQuant libraries prepared with high quality input RNA is shown in Figure 3.

**Important:** As with any library quantitation method, individual operators, lab equipment and instruments will influence the optimal loading quantitation. Validation of the library pool should be determined empirically for each new library type and sequencing platform.

**Figure 3. Fragment distribution when 1 µL of 5 ng/µL library is loaded into a High Sensitivity DNA Chip from 100 ng K562 total RNA starting material.**



3. Prepare libraries for sequencing following the Illumina “Denature and Dilute Libraries Guide” for your specific sequencer.



## V. Technical Support

For help with any of our products, please contact Tecan NGS Technical Support at 650.590.3674 (direct) or 888.654.6544, option 2 (toll-free, U.S. only) or email [techserv-gn@tecan.com](mailto:techserv-gn@tecan.com).

In Europe contact Tecan NGS Technical Support at +31.13.5780215 (Phone) or email [europe-gn@tecan.com](mailto:europe-gn@tecan.com).

In all other locations, contact your Tecan NGS reagent distributor for technical support.

## A. Barcode Sequences and Guidelines for Multiplex Experiments

Unique Dual Index (UDI) barcode sequences for the Universal Plus UDI 96-Plex Adaptor Plate (S02480) are given below. Barcodes in the 24 reaction kits are found in wells A01–H03.

Text versions of all barcode sequences for the Universal Plus UDI 24-Plex Adaptor Plate (S02622) or Universal Plus UDI (S02480), UDI-B (S02690), UDI-C (30185200) and UDI-D (30185201) 96-Plex Adaptor plates can be found at <https://tecangenomics.github.io/>, or contact Tecan NGS Technical Support.

**Table 19. Barcode sequences for Universal Plus UDI 96-Plex Adaptor Plate (Indexes 1-96; S02480)**

PLATE LOCATION	INDEX 1 (I7) SEQUENCE	INDEX 2 (I5) SEQUENCE	PLATE LOCATION	INDEX 1 (I7) SEQUENCE	INDEX 2 (I5) SEQUENCE	PLATE LOCATION	INDEX 1 (I7) SEQUENCE	INDEX 2 (I5) SEQUENCE
A01	CGCTACAT	AACCTACG	A05	AGGTTCTT	TCGAACCT	A09	GCCTTAAC	CCGTTATG
B01	AATCCAGC	GCATCCTA	B05	GAACCTTC	CAAGGTAC	B09	ATTCCGCT	TGTCGACT
C01	CGTCTAAC	CAACGAGT	C05	AAGTCCTC	AGCTACCA	C09	ATCGTGGT	CTCTATCG
D01	AACTCGGA	TGCAAGAC	D05	CCACAACA	CATCCAAG	D09	GCTACAAC	ACTGCTTG
E01	GTCGAGAA	CTTACAGC	E05	ATAACGCC	CTCACCAA	E09	TCTACGCA	CGCCTTAT
F01	ACAACAGC	ACCGACAA	F05	CCGGAATA	TCAGTAGG	F09	CTCCAATC	ATAGGTCC
G01	ATGACAGG	ACATGCCA	G05	CCAAGTAG	GAACGTGA	G09	ACTCTCCA	TGATCACG
H01	GCACACAA	GAGCAATC	H05	AAGGACCA	AGGAACAC	H09	GTCTCATC	CGGATCAA
A02	CTCCTAGT	CCTCATCT	A06	ACGCTTCT	CCTAAGTC	A10	GCCAGAAT	TACTAGCG
B02	TCTTCGAC	TACTGCTC	B06	CTATCCAC	AACGCACA	B10	AATGACGC	TGGACCAT
C02	GACTACGA	TTACCGAC	C06	TGACAACC	GTCAACAG	C10	GTACCACA	GCGCATAT
D02	ACTCCTAC	CCGTAACT	D06	CAGTGCTT	ACACCTCA	D10	ACGATCAG	ATCGCAAC
E02	CTTCCTTC	TTCCAGGT	E06	TCACTCGA	TATGGCAC	E10	TAACGTCG	TCAGCCTT
F02	ACCATCCT	CCATGAAC	F06	CTGACTAC	CGCAATGT	F10	CGCAACTA	CATTGACG
G02	CGTCCATT	TTCCTCCT	G06	GTGATCCA	ACTCAACG	G10	AACACTGG	ACAGGCAT
H02	AACTTGCC	CCAACTTC	H06	ACAGCAAG	GTCTGCAA	H10	CCTGTCAA	AGGTCTGT
A03	GTACACCT	GAGACCAA	A07	TGCTGTGA	CACGATTC	A11	TCCTGGTA	CAGATCCT
B03	ACGAGAAC	ACAGTTCG	B07	CAACACAG	AGAAGCCT	B11	CATCAACC	CTCCTGAA
C03	CGACCTAA	CTAACCTG	C07	CCACATTG	TACTCCAG	C11	AGCAGACA	AGAGGATG
D03	TACATCGG	TCCGATCA	D07	TAGTGCCA	CGTCAAGA	D11	GAAGACTG	CACCATGA
E03	ATCGTCTC	AGAAGGAC	E07	TCGTGCAT	CTGTACCA	E11	TCTAGTCC	CGGTAATC
F03	CCAACACT	GACGAACT	F07	CTACATCC	TCACCTAG	F11	CTCGACTT	GAGTGTGT
G03	TCTAGGAG	TTGCAACG	G07	CATACGGA	AACACCAC	G11	CTAGTCTA	AACTGAGG
H03	CTCGAACA	CCAACGAA	H07	TGGTAAC	CGTCTTCA	H11	TCCAACCTG	TGTGTCAG
A04	ACGGACTT	ATCGGAGA	A08	CAGGTTCA	AACGTAGC	A12	GACATCTC	TGTCACAC
B04	CTAAGACC	CCTAACAG	B08	AGAACCAG	GCAACCAT	B12	ACTGCACT	AGATCGTC
C04	AACCGAAC	CATACTCG	C08	GAATGGCA	GATCCACT	C12	GTTCCATG	CAATGCCA
D04	CCTTAGGT	TGCCTCAA	D08	AGGCAATG	ACCTAGAC	D12	ACCAAGCA	TGCTTGCT
E04	CCTATACC	TACAGAGC	E08	TAGGAGCT	CTAGCAGT	E12	CTCTCAGA	AATGGTCG
F04	AACGCCTT	CGAGAGAA	F08	CGAACAAC	TCGATGAC	F12	ACTCTGAG	AGTTGTGC
G04	TCCATTGC	AGGTAGGA	G08	CATTGCTC	TTGGTGCA	G12	GCTCAGTT	GTATCGAG
H04	CAAGCCAA	GAACGAAG	H08	AGCCAACCT	AGTGCATC	H12	ATCTGACC	GTACGATC

## VI. Appendix

For UDI barcodes B-D and text versions of all barcode sequences visit <https://tecangenomics.github.io/>



**Note:** Manual demultiplexing of sequencing data generated on the Illumina MiniSeq, NextSeq, NovaSeq or HiSeq 4000 (PE runs only) instruments may require inputting the reverse complement of the Index 2 (i5) sequences.

### B. Library Amplification qPCR Optimization for Standard Workflow

When using the Universal Plus kit for the first time, or when working with a new sample type or input amount, we recommend performing a qPCR step prior to Library Amplification to determine the optimum number of cycles needed and ensure there is no excess amplification.

Perform a 1/10th scale qPCR reaction as follows:

**Table 20. Library Amplification qPCR Master Mix**

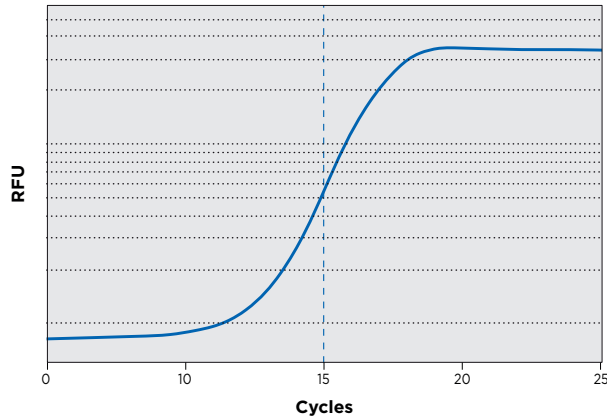
REAGENT	AMPLIFICATION REAGENT I (RED: AR1 VER 1)	AMPLIFICATION REAGENT II (RED: AR2 VER 1)	AMPLIFICATION ENZYME MIX (RED: AR3 VER 1)	STRAND SELECTION ENZYME MIX II (PURPLE: SS4)	20X EvaGreen	NUCLEASE-FREE WATER (GREEN: D1)
1X REACTION VOLUME	1 $\mu$ L	0.8 $\mu$ L	0.05 $\mu$ L	0.1 $\mu$ L	0.5 $\mu$ L	6.05 $\mu$ L

1. Prepare a PCR master mix according to the volumes shown in Table 20.
2. Aliquot 8.5  $\mu$ L of PCR master mix per sample into a 0.2 mL qPCR strip or plate.
3. Add 1.5  $\mu$ L of purified library (after **K. Strand Selection Purification** but before **L. Library Amplification A**) for a total qPCR volume of 10  $\mu$ L. Mix well by pipetting, spin and place on ice.
4. Perform real-time qPCR with the following cycling conditions:

37 °C – 10 min, 95 °C – 2 min, 2x(95 °C – 30 s, 60 °C – 90 s), 30x(95 °C – 30 s, 65 °C – 90 s), 65 °C – 5 min, hold at 4 °C

The cycle number used for subsequent library amplification should be within the exponential phase of the amplification plot (Figure 4).

**Figure 4. Stylized qPCR Amplification Plot**



To amplify the remaining 13.5  $\mu\text{L}$  of library, prepare an amplification master mix according to the volumes shown below (Table 21) for standard workflows.

**Table 21. Library Amplification Master Mix**

REAGENT	AMPLIFICATION REAGENT I (RED: AR1 VER 1)	AMPLIFICATION REAGENT II (RED: AR2 VER 1)	AMPLIFICATION ENZYME MIX (RED: AR3 VER 1)	STRAND SELECTION ENZYME MIX II (PURPLE: SS4)	NUCLEASE-FREE WATER (GREEN: D1)
<b>1X REACTION VOLUME</b>	9 $\mu\text{L}$	7.2 $\mu\text{L}$	0.45 $\mu\text{L}$	0.9 $\mu\text{L}$	58.95 $\mu\text{L}$

5. Remove Agencourt beads from storage and place on the bench top to reach room temperature for use in the next step.
6. Thaw AR1 and AR2 at room temperature. Mix by vortexing, spin and place on ice.
7. Spin down AR3 and SS4 and place on ice.
8. Make a master mix by sequentially combining D1, AR1 and AR2 in an appropriately sized capped tube according to the volumes shown in Table 21. Add AR3 and SS4 at the last moment and mix well by pipetting, taking care to avoid bubbles. Spin the tubes and place on ice.
9. On ice, add 76.5  $\mu\text{L}$  of Library Amplification Master Mix to 13.5  $\mu\text{L}$  of sample for a total of 90  $\mu\text{L}$ . Mix by pipetting, spin and place on ice.
10. Place the tubes in a pre-warmed thermal cycler programmed to run Program 9 (Library Amplification; see Table 5), using the number of cycles (N) determined by qPCR:  
37  $^{\circ}\text{C}$  - 10 min, 95  $^{\circ}\text{C}$  - 2 min, 2x(95  $^{\circ}\text{C}$  - 30 s, 60  $^{\circ}\text{C}$  - 90 s), Nx(95  $^{\circ}\text{C}$  - 30 s, 65  $^{\circ}\text{C}$  - 90 s), 65  $^{\circ}\text{C}$  - 5 min, hold at 4  $^{\circ}\text{C}$
11. Remove the tubes from the thermal cycler, spin to collect condensation and place on ice.
12. Add 10  $\mu\text{L}$  of nuclease-free water to each library. Mix well by pipetting and continue immediately to section **P. Amplified Library Purification**.

**C. Library Amplification qPCR Optimization for AnyDeplete Workflow**

When using the Universal Plus kit for the first time, or when working with a new sample type or input amount, we recommend performing a qPCR step prior to Library Amplification to determine the optimum number of cycles needed and ensure there is no excess amplification.

**Protocol**

A. Perform a 1/10th scale qPCR reaction as follows:

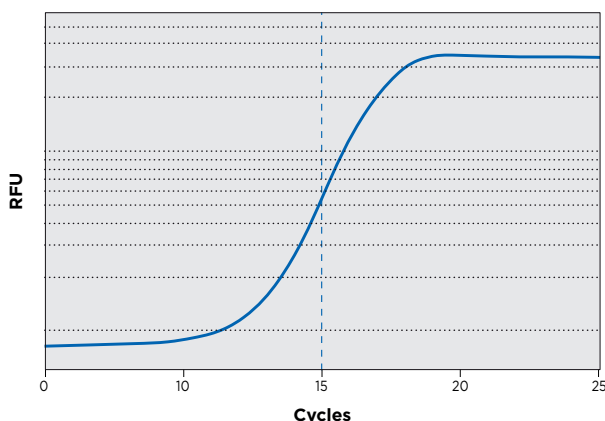
**Table 22. Library Amplification qPCR Master Mix**

REAGENT	AMPLIFICATION REAGENT I (RED: AR1 VER 1)	AMPLIFICATION REAGENT II (RED: AR2 VER 1)	AMPLIFICATION ENZYME MIX (RED: AR3 VER 1)	20X EvaGreen	NUCLEASE-FREE WATER (GREEN: D1)
1X REACTION VOLUME	1 µL	0.8 µL	0.05 µL	0.5 µL	2.65 µL

1. Prepare a PCR master mix according to the volumes shown in Table 22.
2. Aliquot 5 µL of PCR master mix per sample into a 0.2 mL qPCR strip or plate.
3. Add 5 µL of library (after **N. Targeted Depletion (optional AnyDeplete workflow)** but before **O. Library Amplification B (optional AnyDeplete workflow)**) for a total qPCR volume of 10 µL. Mix well by pipetting, spin and place on ice.
4. Perform real-time qPCR with the following cycling conditions:  
95 °C – 2 min, 2x(95 °C – 30 s, 60 °C – 90 s), 30x(95 °C – 30 s, 65 °C – 90 s), 65 °C – 5 min, hold at 4 °C

The cycle number used for subsequent library amplification should be within the exponential phase of the amplification plot (Figure 5).

**Figure 5. Stylized qPCR Amplification Plot**



- B. To amplify the remaining 45 µL of library, prepare an amplification master mix according to the volumes shown below (Table 23) for AnyDeplete workflows.

**Table 23. Library Amplification B Master Mix**

REAGENT	AMPLIFICATION REAGENT I (RED: AR1 VER 1)	AMPLIFICATION REAGENT II (RED: AR2 VER 1)	AMPLIFICATION ENZYME MIX (RED: AR3 VER 1)	NUCLEASE- FREE WATER (GREEN: D1)
<b>1X REACTION VOLUME</b>	9 $\mu$ L	7.2 $\mu$ L	0.45 $\mu$ L	28.35 $\mu$ L

1. Remove Agencourt beads from storage and place on the bench top to reach room temperature for use in the next step.
2. Thaw AR1 and AR2 at room temperature. Mix by vortexing, spin and place on ice.
3. Spin down AR3 and place on ice.
4. Make a master mix by sequentially combining D1, AR1 and AR2 in an appropriately sized capped tube according to the volumes shown in Table 23. Add AR3 at the last moment and mix well by pipetting, taking care to avoid bubbles. Spin the tubes and place on ice.
5. On ice, add 45  $\mu$ L of Library Amplification B Master Mix to 45  $\mu$ L of sample for a total of 90  $\mu$ L. Mix by pipetting, spin and place on ice.
6. Place the tubes in a pre-warmed thermal cycler programmed to run Program 12 (Library Amplification B; see Table 5), using the number of cycles (N) determined by qPCR:  
95 °C – 2 min, 2x(95 °C – 30 s, 60 °C – 90 s), Nx(95 °C – 30 s, 65 °C – 90 s), 65 °C – 5 min, hold at 4 °C
7. Remove the tubes from the thermal cycler, spin to collect condensation and place on ice.
8. Continue to section **P. Amplified Library Purification**.

## D. Frequently Asked Questions (FAQs)

### Getting Started

**Q1. What materials are provided with Universal Plus mRNA-Seq with NuQuant?**

The Universal Plus mRNA-Seq with NuQuant kit bundle provides all necessary buffers, primers and enzymes for poly(A) selection, library construction, AnyDeplete (Part No. 0521, 0521B, 0521C, 0521D, 0522, 0522B, 0522C, 0522D only) and NuQuant. SPRI purification beads and EvaGreen are not included.

**Q2. What is the concentration of the control RNA?**

The control RNA (K562 total RNA), included only with the 24 reaction kit, is provided at a concentration of 1  $\mu$ g/ $\mu$ L.

**Q3. What equipment is required or will be useful?**

A comprehensive list of required and recommended equipment can be found in Section **II.B**.

**Q4. Can this system be used with other library preparation workflows?**

Universal Plus mRNA-Seq is an end-to-end solution designed to generate libraries for Illumina sequencing starting from total RNA and has not been tested with alternative library preparation systems.

### Input Recommendations

**Q5. What methods do you recommend for RNA isolation?**

We recommend a column-based method, including:

- Norgen Biotek Total RNA Purification Kit
- Zymo Research Quick-RNA™ Kits
- Arcturus PicoPure® RNA Isolation Kit
- Ambion PureLink® RNA Mini Kit
- Qiagen RNeasy Kits

Organic methods such as TRIzol® Reagent should be subsequently followed with a column-based clean-up method.

**Q6. Can I use TRIzol or other phenol-chloroform based extractions for RNA isolation?**

We do not recommend the use of TRIzol or similar methods as any carry over of organic solvents may inhibit downstream enzyme activity. If using TRIzol extracted RNA, we recommend using a column-based purification of the RNA prior to input into the kit.

**Q7. Can I use carrier RNA during RNA isolation?**

We do not recommend the use of carriers during RNA isolation. If a carrier is required, please contact Tecan NGS Technical Support for more information.

**Q8. Can I use Universal Plus mRNA-Seq with RNA from any organism?**

Universal Plus mRNA-Seq has been designed for use with total RNA inputs containing polyadenylated transcripts. Universal Plus mRNA-Seq is compatible with polyadenylated RNA from any organism.

**Q9. Do I need to use high-quality total RNA?**

Yes, high quality RNA is required for efficient poly(A) selection.

**Q10. What input amounts can be used with the Universal Plus mRNA-Seq with NuQuant kit?**

The Universal Plus mRNA-Seq kit has been designed for inputs between 10 ng - 1 µg of total RNA as input into the mRNA Selection Module. Inputs outside of these ranges may affect reaction stoichiometry, resulting in suboptimal libraries. Lower input amounts will potentially result in insufficient yields depending on the requirements of the analytical platform.

**Q11. Can contaminating genomic DNA interfere with Universal Plus mRNA Seq?**

Contaminating genomic DNA can be incorporated into libraries and inclusion of a DNase treatment during RNA isolation is recommended.

### General Workflow

**Q12. Does this system contain a SPIA®-based amplification?**

No. The cDNA is generated with random and poly(T) primers, but no SPIA-based amplification is used.

**Q13. Is it necessary to fragment my cDNA prior to End Repair and Adaptor Ligation?**

No. Chemical fragmentation is incorporated into the workflow after poly(A) RNA selection.

**Q14. Can I combine the barcoded libraries prior to the PCR amplification step?**

This is not recommended. The stoichiometry of barcoded libraries may be adversely affected by this modification to the workflow. We suggest that the libraries be amplified and quantitated independently before being pooled for use on the sequencer.

**Q15. Where can I safely stop in the protocol?**

Samples can be placed in short-term storage at  $-20\text{ }^{\circ}\text{C}$  after second strand synthesis, end repair, strand selection or after any of the bead purification steps.

### SPRI Bead Purifications

**Q16. What is the difference between RNAClean XP and AMPure XP SPRI beads? Can both be used interchangeably?**

RNAClean XP beads are certified to be RNase and DNase free. We have tested both RNAClean XP and AMPure XP beads in our kits and observe no difference in performance between products during the purification steps in this protocol.

**Q17. What magnetic separation devices do you recommend for the SPRI bead purifications?**

Due to the large number of commercially available magnets, we do not have a comprehensive list of compatible products. However, many magnets are compatible. As long as the magnet is strong enough to clear the solution of magnetic beads, it can be applied to the system. We have the following guidelines for selecting a magnetic separation device:

- a. Use a magnet designed for 0.2 mL tubes (PCR tubes), tube strips, or plates. Compared to magnets that are designed for 1.5 mL tubes, these minimize loss that can occur when samples are transferred from one tube to another.
- b. Prior to purchasing, check the manufacturer's specifications for minimum and maximum volumes that can be effectively treated.
- c. Test the magnet with a mock purification to ensure the magnet will effectively clear the solution under the conditions in this workflow. This is also helpful to gain familiarity with the purification workflow.

**Q18. How can I ensure maximum recovery of sample from the SPRI bead purification?**

- a. Allow the SPRI beads to reach room temperature before use; cold beads result in lower yields.
- b. Ensure that the beads are fully resuspended in solution before adding to the sample.
- c. Always use fresh ethanol during the washing steps. When preparing the ethanol, measure out the ethanol and water separately to ensure the desired ethanol concentration is obtained.
- d. Mix the bead suspension and sample thoroughly to ensure maximum binding of the samples to the beads.

### Library Quantification and Qualification

**Q19. How do I measure my final library yield?**

We recommend using NuQuant to accurately quantify the final libraries for multiplex pooling using a Qubit instrument. The final library pool concentration should be determined using a qPCR-based method before loading onto an Illumina sequencer. Please refer to section **IV.Q** for guidelines on alternative library quantitative and qualitative assessments.



**Q20. How many bases do the UDI adaptors add to the library?**

The adaptors add 144 bp to the library.

### Sequencing Recommendations

**Q21. What sequencers are compatible with your libraries?**

Universal Plus mRNA-Seq libraries are compatible with Illumina sequencing platforms.

**Q22. How much material should I load into the sequencer?**

Please follow manufacturer's recommendations for library QC, quantitation, balancing and loading of the amplified library on the sequencer.

**Q23. What kind of error correction is used to minimize the impact of sequencing errors in the barcodes?**

The barcodes provided in the Universal Plus UDI 96-Plex Adaptor Plate (S02480) are a minimum edit distance of 3 from other barcodes in the adaptor plate. This means that a minimum of three edits (replacement, insertion, or deletion) must occur before one barcode becomes a different barcode. The barcodes provided in the Universal Plus UDI-B 96-Plex Adaptor Plate (S02690), Universal Plus UDI-C 96-Plex Adaptor Plate (30185200) or Universal Plus UDI-D 96-Plex Adaptor Plate (30185201) are a minimum edit distance of 2 from other barcodes in the adaptor plate. For further details on the barcode design strategy, please refer to Faircloth BC, Glenn TC (2012), Not All Sequence Tags Are Created Equal: Designing and Validating Sequence Identification Tags Robust to Indels. PLoS ONE 7(8): e42543. doi:10.1371/journal.pone.0042543.

**Q24. What kind of sequencing primers can I use with your libraries?**

The Universal Plus mRNA-Seq with NuQuant kit is designed for use with the standard Illumina sequencing primers for both single-end and paired-end sequencing applications. The final libraries contain a unique 8 base index 1 and a unique 8 base index 2 suitable for detection of "index hopping." Both indexes must be sequenced to enable detection of "index hopping."

**Q25. Can Universal Plus mRNA-Seq libraries be used with paired-end sequencing?**

Yes. The libraries produced using this kit can be used for both single-end and paired-end sequencing. Special consideration should be given to the expected insert size in the paired-end assay. Contact Tecan NGS Technical Support for additional information.

### Data Analysis

**Q26. Can I use standard alignment algorithms to analyze strand-specific sequencing data?**

Yes. Strand-specific reads can be processed and mapped to reference sequences using the same methods used for other RNA-Seq libraries. Note that in libraries generated by Universal Plus mRNA-Seq, the forward read corresponds to the sense strand.

### Custom AnyDeplete

Custom depletion designs can be tailored to any transcript, any organism. Please contact techserv-gn@tecan.com for more information.

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

**Tecan Genomics, Inc.**



**USA**

900 Chesapeake Drive  
Redwood City, CA 94063 USA

**Customer Service and**

**Technical Support:**

Toll Free Tel: 888.654.6544  
Toll Free Fax: 888.296.6544  
cservice-gn@tecan.com  
techserv-gn@tecan.com

**Europe**

P.O. Box 109,  
9350 AC Leek  
The Netherlands  
**Customer Service and**  
**Technical Support:**  
Tel: +31-13-5780215  
Fax: +31-13-5780216  
europe-gn@tecan.com

**Worldwide**

For our international  
distributors contact  
information, visit  
our website  
[www.nugen.com](http://www.nugen.com)

.....  
**Australia** +61 3 9647 4100 **Austria** +43 62 46 89 330 **Belgium** +32 15 42 13 19 **China** +86 21 220 63 206 **France** +33 4 72 76 04 80 **Germany** +49 79 51 94 170  
**Italy** +39 02 92 44 790 **Japan** +81 44 556 73 11 **Netherlands** +31 18 34 48 17 4 **Nordic** +46 8 750 39 40 **Singapore** +65 644 41 886 **Spain** +34 93 595 25 31  
**Switzerland** +41 44 922 89 22 **UK** +44 118 9300 300 **USA** +1 919 361 5200 **Other countries** +41 44 922 81 11  
.....

Tecan Group Ltd. makes every effort to include accurate and up-to-date information within this publication; however, it is possible that omissions or errors might have occurred. Tecan Group Ltd. cannot, therefore, make any representations or warranties, expressed or implied, as to the accuracy or completeness of the information provided in this publication. Changes in this publication can be made at any time without notice. For technical details and detailed procedures of the specifications provided in this document please contact your Tecan representative. This brochure may contain reference to applications and products which are not available in all markets. Please check with your local sales representative.

Tecan, DimerFree®, NuQuant®, and AnyDeplete® are registered trademarks and trademarks of Tecan Group Ltd., Männedorf, Switzerland or of Tecan Genomics, Inc., Redwood City, USA.

© 2020 Tecan Genomics, Inc., all rights reserved. For disclaimer and trademarks please visit [www.tecan.com](http://www.tecan.com)

[www.tecan.com](http://www.tecan.com)

