



# NGS library prep automated workstation solutions

Over 120 tested and vendor qualified NGS methods



## Table of contents

### Instruments..... 3

Automated library prep throughput and walkaway-time .....	3
Fontus NGS workstation .....	4
Sciclone G3 NGSx iQ workstation .....	4
Sciclone G3 NGSx workstation. ....	5
Zephyr G3 NGS workstation. ....	5
Zephyr G3 NGS iQ workstation .....	5
BioQule NGS System .....	5

### Accessories..... 6

Workstation table.....	6
Deck Verification. ....	6

### Consumables and kits..... 7

Pipette tips .....	7
Microplates and lids .....	7
Complete NGS consumable kits .....	8
High resolution library sizing with the LabChip GX	
Touch HT nucleic acid analyzer .....	9
NEXTFLEX NGS library preparation kits.....	9
Service plans and applications development .....	9

### Verified, automated NGS methods..... 10

DNA Library Preparation Kits .....	10
RNA Library Preparation Kits .....	12
Small RNA Library Preparation Kits.....	13
Exome/Target Capture Library Preparation Kits .....	14
Metagenomics Library Preparation Kits .....	15
Supplementary Applications & Protocols.....	15

# Instruments

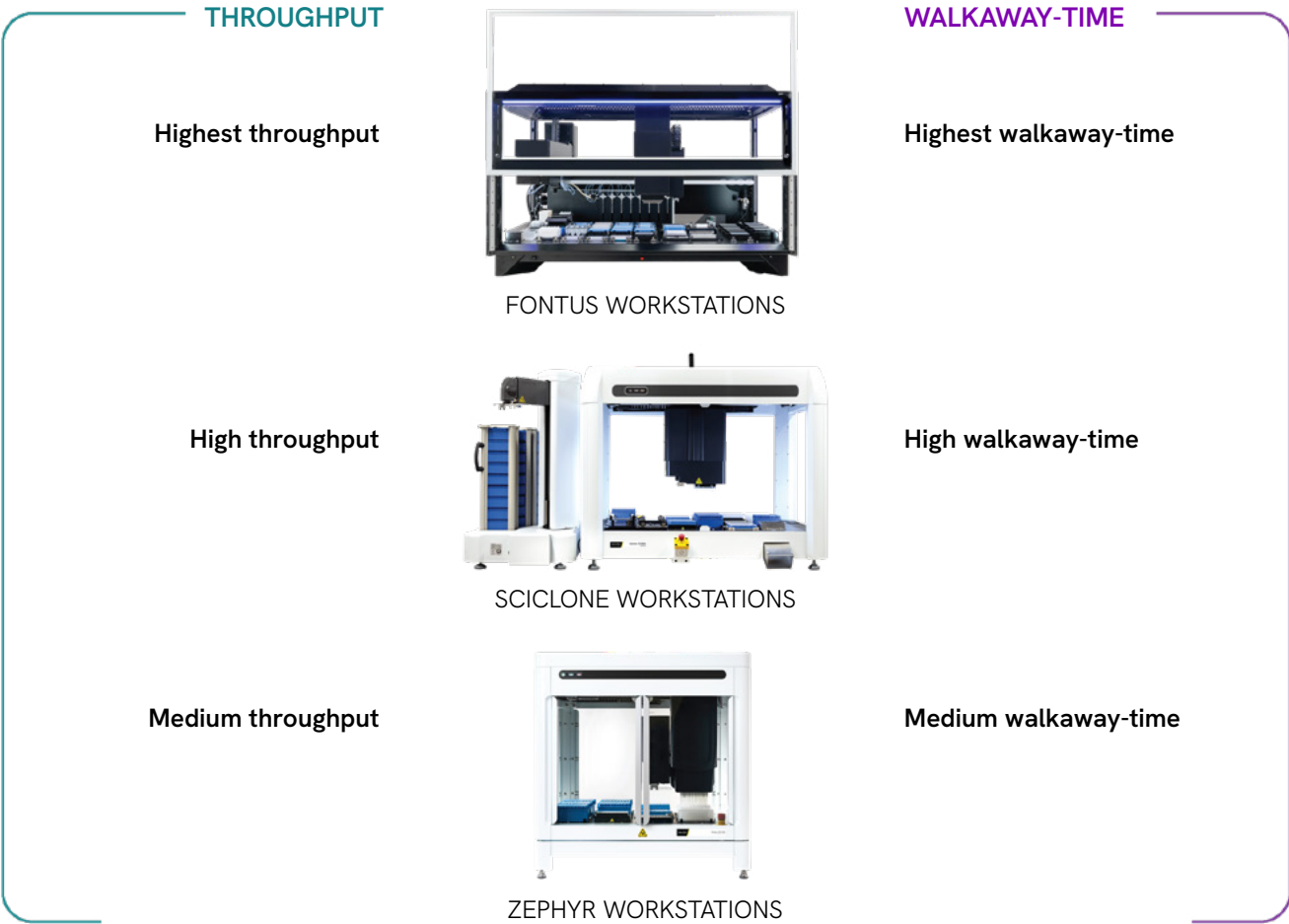
Revvity’s automated library prep solutions are designed and programmed for next generation sequencing (NGS) applications. They feature a library of over 120 pre-developed and verified or vendor qualified automated NGS preparation kits such as:

- DNA-Seq
- RNA-Seq
- Exome
- Small RNA-Seq
- Oncology
- Target Capture
- Metagenomics
- Supplementary Protocols

Pre-developed and verified NGS methods allow labs to get up and running quickly and gives confidence that their methods will perform to specifications.

Automating NGS protocols ensures precise step execution while minimizing variability and risk as well as providing extended walk-away time for researchers, freeing them up for higher value tasks. Revvity has a complete line of up NGS automation solutions from low to high throughput and partial to full walk away.

## Automated library prep throughput and walkaway-time



## Fontus NGS workstation

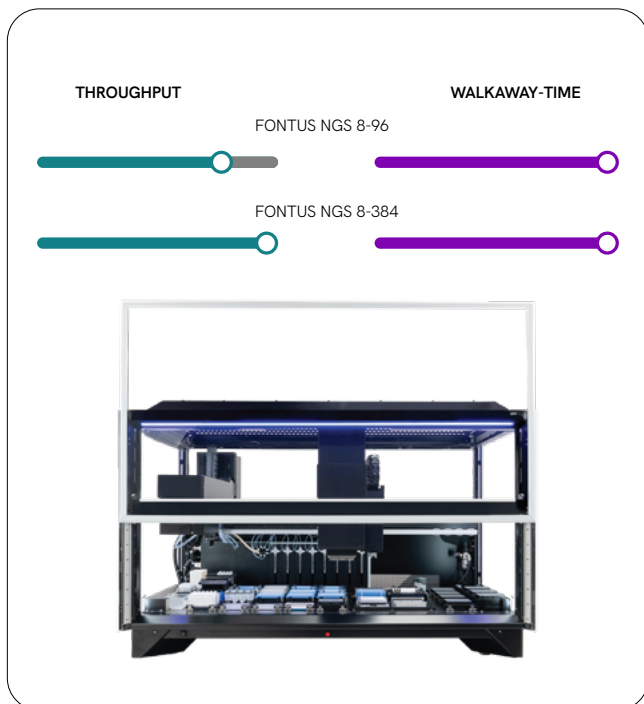
FONTUS NGS 8-96 - YFL8H02

FONTUS NGS 8-384 - YFL8L02

For the most complete automation and minimum user intervention (longest walkaway), the Fontus™ NGS workstation includes either a 96 or 384 pipetting head for processing up to 96 or 384 samples simultaneously. The Fontus system also includes a highly flexible variable 8-span pipetting arm (Varispan™), which can perform all the initial vials/tubes-to-plates reformatting as well as fast normalization and post assay sample pooling.

The Fontus NGS 8-384 enables miniaturization, which can drastically reduce reagent and per-sample costs while quadrupling throughput.

With full walkaway and the highest throughput, the Fontus NGS workstations are the ideal choice for automation of NGS library prep.

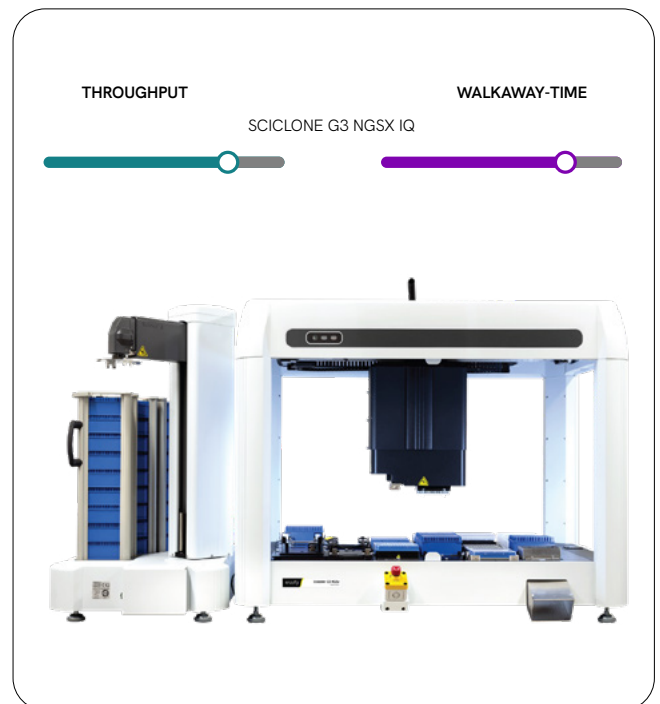


## Sciclone G3 NGSx workstation

SCICLONE G3 NGSx iQ (96) - CLS152236

Designed for longer walkaway times than the Sciclone NGSx workstation, this system shares all the same features and also includes an integrated on-deck thermal cycler and a Twister® III robotic arm. The ODTc enables full automation of PCR amplification and heated incubations and the Twister III robot stores and loads up to 27 tip boxes of various labware. This provides increased walkaway time for users and lowers variability and risks related to off-deck incubation and PCR.

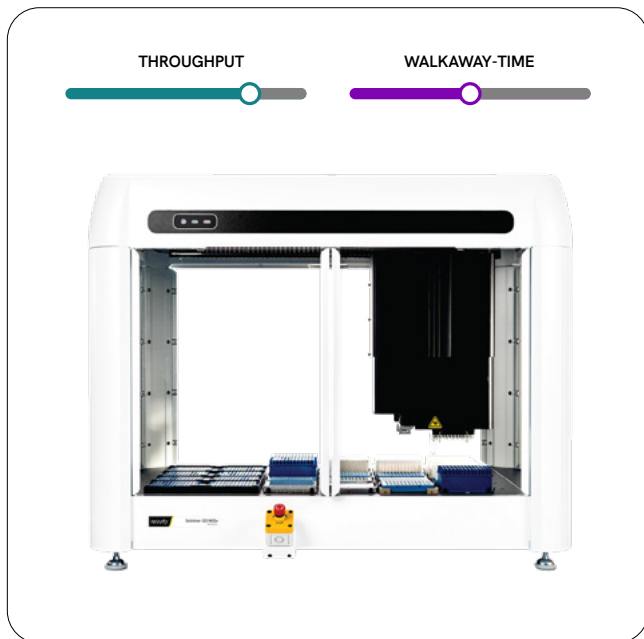
For very high throughput NGS protocols, the Sciclone® G3 NGSx HT workstation includes all the walkaway benefits of the iQ workstation but with a 384 pipetting head enabling assay miniaturization, which can drastically reduce reagent and per-sample costs.



## Sciclone G3 NGSx workstation

CLS152932

Equipped with a high precision, 96-channel pipetting head, this enables batch processing of up to 96 samples simultaneously. It includes an integrated gripper which facilitates moving labware around the deck. Below deck storage allows enough tip storage to accommodate large NGS protocols and a tip chute is included to seamlessly remove consumables from the deck. This fully enclosed workstation also includes temperature, magnetic and shaking options for precise, standardized NGS library preparation.



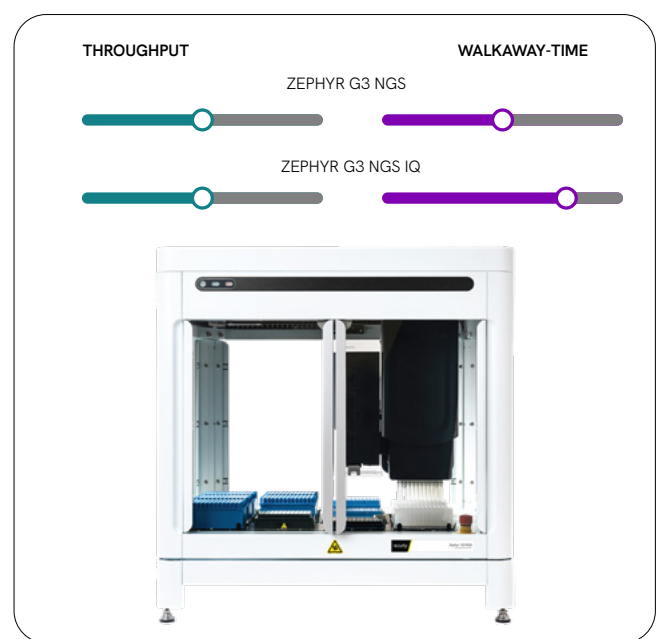
## Zephyr G3 NGS workstation

ZEPHYR G3 NGS - CLS145300

ZEPHYR G3 NGS iQ - CLS158259

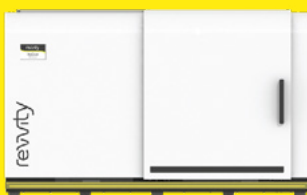
For low to medium-throughput NGS needs, the Zephyr® G3 NGS workstation is a compact, semiautomated solution designed to construct up to 96 NGS libraries per day. This compact workstation is equipped with a highly precise, 96-channel pipetting head and includes an integrated gripper. Also included in this compact workstation is on-deck temperature and shaking integrated peripherals and a waste tip chute.

The Zephyr G3 NGS iQ workstation includes all the same benefits of the Zephyr G3 NGS, but also includes a addition of an integrated thermocycler and a Twister robot to enable a higher level of automation and longer unattended NGS protocol runs.



## BioQule NGS system

[CLS155700](#)



The BioQule™ NGS system is designed for low throughput NGS library prep and is ideal for labs moving into automation from manual processing. It seamlessly processes up to 8 samples per run and features pre-plated kits, plates and consumables. Just add your samples, reagents, and run!



## Accessories

These optional add-ons and accessories can be used to further integrate or provide additional functionality to the liquid handlers for NGS workflows.

### Workstation table

The specifically designed tables hold the liquid handler workstation and also has compartments for organizing the under-table hardware in a clean, organized format. The table includes specific compartments for:

- INHECO® ODT controller
- Compressor
- Tip Waste Bucket
- Fuse box
- Computer
- Wheel Lock
- Network power and air distribution

[Buy online](#)

#### Fontus NGS workstation table (HH20230001)

Width: 158 cm | Height: 92 cm | Depth: 73 cm

#### Sciclone G3 NGSx workstation table (119877)

Width: 120 cm | Height: 91 cm | Depth: 66 cm

#### Sciclone G3 NGSx iQ and HT workstation table (CLS152481)

Width: 171 cm | Height: 91 cm | Depth: 66 cm

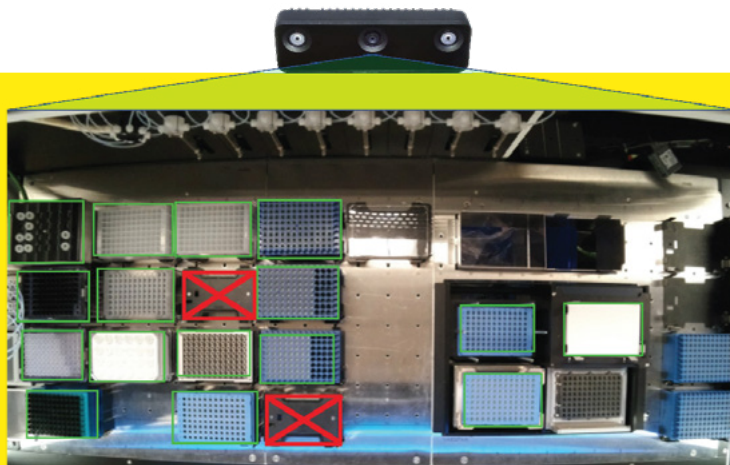
#### Zephyr G3 NGS iQ workstation table (CLS158557)

Width: 171 cm | Height: 91 cm | Depth: 66 cm

[More info](#)

### Fontus™ Deck Verification

**Identify potential errors before they happen.**



Deck Verification is an optional vision system enabled through the WinPREP™ software. A camera integrated above the deck automatically verifies the placement of labware relative to the selected protocol prior to a run.

This ensures correct deck configuration and labware setup is correct prior to starting the run. This can eliminate mid-run errors due to empty tip boxes, misplaced labware and reagents and avoid potential collisions. The vision system has the ability to detect the correct consumables, the presence or absence of disposable tips and determine the number of tips racks present in the recessed deck.

## Consumables and kits

Our consumables are manufactured and designed to provide a dependable standardized NGS preparation performance. Our NGS consumable kits combine consumable tips, microplates and lids to perform a single run of 96 samples for a variety of standardized protocols.

### Pipette tips

[Buy tips online](#)

Revvity manufactures and provides high quality pipette tips for the Fontus, Sciclone, and Zephyr workstations. These are made to the highest standards and are tested and verified to ensure peak performance of NGS protocols.

All of our sterile tips are sterilized using irradiation to the current regulatory maximum of  $10^{-6}$  SAL which is equivalent to the sterility assurance level (SAL) required by the medical device industry. This sterilization ensures the product is free from the following contaminants: human DNA, DNase and RNase and bioburden. Additionally, our filter tips include a filter membrane which functions as a barrier between the pipetting head and the liquid as it is aspirating and dispensing. The filters ensure no-cross contamination is possible from aerosols and reduces overall aerosol intake into the pipetting head.



### Microplates and lids

[Buy online](#)

The Fontus, Sciclone, and Zephyr workstations are designed to work with SBS format labware. Widely utilized plates on these workstations include:

- Deepwell microplates: Storing large volumes and for washing steps
- U-bottom microplates: General use
- V-bottom microplates: Reduces dead volume in pipetting steps
- PCR plates: Preparing samples for NGS libraries  
Revvity makes a complete line of labware optimized for NGS assays and automation. Please visit [www.revvity.com](http://www.revvity.com) to view our offering.



## Complete NGS consumable kits

Revvity offers consumable kits specific to NGS library prep methods. These packs can help user ensure high quality libraries and no missing labware items.



### KIT A (6060008)

- NEBNext® Ultra™ II DNA Library Prep Kit for Illumina® (E7645L)
- QIAGEN QIAseq FX DNA Library Kit (180479180482, 180484)

### KIT B (6060009)

- Illumina® TruSeq™ RNA Sample Prep (RS-122-2001/ RS-122-2002)
- Illumina® TruSeq™ Stranded Total RNA LT Sample Prep (RS-122-2201/RS-122-2202/RS-122-2301/RS-122-2302)
- Illumina® TruSeq™ Stranded Total mRNA LT Sample Prep (RS-122-2101/RS-122-2102/RS-122-2103)
- Illumina® Stranded mRNA Prep, Ligation (20040534)
- Illumina® Stranded Total RNA Prep, Ligation with Ribo-Zero Plus (20040529)

### KIT C (6060010)

- Agilent® SureSelectXT Library Prep and Target Enrichment for Illumina Sequencing (G9641B/G9641C/G9642B/G9642C)
- Agilent® SureSelectXT2 Library Prep and Target Enrichment for Illumina Sequencing (G9621A/G9621B/G9621C)
- Agilent® SureSelect XT HS Target Enrichment System for Illumina Sequencing (G9702C)
- Agilent® SureSelect XT HS2 DNA Library Preparation and Target Enrichment for Illumina Sequencing (G9985A-D or G9956A-D, G9957B)
- Illumina® RNA Prep with Enrichment, (L) Tagmentation (20040537)

### KIT D (6060017)

- NEXTFLEX Variant-Seq™ SARS-CoV-2 V1 (NOVA5349-096)
- NEXTFLEX Variant-Seq™ SARS-CoV-2 V2 (NOVA5350-096)
- NEXTFLEX Rapid XP Kit v2 (NOVA-5149-21, NOVA5149-22, NOVA-5149-23)
- NEXTFLEX Rapid DNA-Seq Kit 2.0 (NOVA-5188-01/ NOVA-5188-02/NOVA-5188-03/ NOVA-5188-11/ NOVA-5188-12/NOVA-5188-13)

### KIT E (6060018)

- NEXTFLEX Rapid Directional RNA-Seq Kit 2.0 (NOVA-5189-01/NOVA-5198-02/NOVA-5198-03)
- WITH: NEXTFLEX RiboNaut rRNA Depletion Kit (NOVA-512961/NOVA-512962/NOVA-512963)
- OR: NEXTFLEX Poly(A) Beads 2.0 (NOVA-512991/NOVA-512992/NOVA-512993)
- QIAGEN QIAseq miRNA UDI Library Kit (331505)



## High resolution library sizing with the LabChip GX Touch nucleic acid analyzer

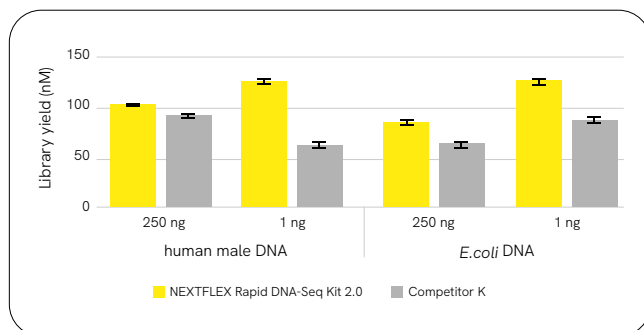
The LabChip® GX Touch™ Nucleic Acid Analyzer provides precise and reliable sizing distribution of NGS libraries to determine your library quality. This advanced system utilizes assays optimized for NGS library quality control and can run up to 192 NGS libraries per run with a per sample analysis time of 68 seconds.



## NEXTFLEX NGS library preparation kits

High-quality library preparation is a critical step in Next Generation Sequencing (NGS). Revvity provides a complete portfolio of DNA and RNA NGS library preparation kits and barcodes designed to increase the sensitivity, flexibility, and speed of library prep for the Illumina® and Ion Torrent™ sequencing platforms. We have focused our attention on optimizing our products to cater to the customer's need for a robust and cost-effective solution to achieve reproducible, high-quality sequencing results every time.

Pre-plated automation ready versions of our library prep kits are available for high throughput needs while manual versions cater to the needs of lower throughput users.



Rapid DNA library prep for sequencing on Illumina® Platforms.

Learn more about rapid DNA library prep



## Service plans and applications development

Revvity provides dependable coverage for our workstations through service plans and applications development.

Our service plans are designed to enhance your lab's productivity by ensuring critical maintenance requirements aren't compromised.

Our service engineers have an average of 16 years of experience and are cross trained to provide increased support and efficient resolution of workstation issues. Our service plans guarantee a rapid response from our expert engineers to ensure your workstation is in top shape for your important work.

Our applications team provides development services for NGS kit automation. These scientists specialize in automating NGS kits on our Fontus, Sciclone, and Zephyr workstations. Application services can also support development of protocols to automate any new kits for your laboratory processes.

## Verified, automated NGS methods

A variety of NGS reagent kits are automated on Revvity liquid handlers. Each automated protocol is qualified with sequencing data prior to deploying to customers. After we develop the method and thorough wet testing, reagents are tested on the liquid handler with control sample inputs. The final NGS libraries are quality controlled and sequenced to confirm proper library prep by the automation. Our application scientists and subject matter experts ensure the automated liquid handler produces acceptable results for a range of sample inputs, with no cross contamination between wells or column bias.

## DNA library preparation kits

List is current as of Nov. 2024, inquire for updated list.

Vendor	Library Prep Kit	Sciclone™ G3 NGSx	Sciclone™ G3 NGSx iQ™	Zephyr™ G3 NGS	Fontus™ NGS 96	BioQule™
10x Genomics®	Genomic Preparation	X				
Agilent®	SureSelect XT HS2 DNA	X*	X*			
ArcherDx®	Universal DNA Reagent Kit v2	X				
CareDx®	AlloSeq TZ			X		
Celeomics®	Whole Exome Sequencing (WES), Library Preparation, Target Enrichment	X				
Claret Bioscience	ClaretBio® SRSly® PicoPlus Kit					X
IDT®	xGen DNA Library Prep Kit EZ	X	X			
IDT®	xGen cfDNA & FFPE DNA Library Prep MC v2 Kit	X				
IDT®	xGen™ Pre-Hybridization Capture Normalase™ Module with xGen DNA Library Prep Kit EZ	In Dev				
IDT®	Accel-NGS® 2S Plus DNA Library Kit	X				
IDT®	xGen Amplicon Panels	X	X			
Illumina®	Illumina® Complete Long Reads with Enrichment	X				
Illumina®	Illumina® DNA Prep with Exome 2.0 Plus Enrichment	X	X	X**		
Illumina®	Illumina® DNA PCR-Free Library Prep Tagmentation	X	X		X*	X
Illumina®	Illumina® DNA Prep (previously called Nextera DNA Flex)	<a href="#">[X]</a>	X	X**		X
Illumina®	Illumina® DNA Prep with Enrichment	X	X			
Illumina®	Illumina® DNA Prep with Exome 2.5 Enrichment	<a href="#">[X]</a>	<a href="#">[X]</a>			
Illumina®	Nextera® XT DNA Kit	X	X	X		
Illumina®	TruSeq® DNA PCR-Free Kit	X	X			
Illumina®	TruSeq® Nano DNA Kit	X				
New England Biolabs® / NEB®	NEBNext® dsDNA Fragmentase® Kit	X				
NEB®	NEBNext® Ultra™ DNA Kit	X		X		
NEB®	NEBNext® Ultra™ II DNA Library Prep for Illumina® Instruments	X	X			
NuGen® / Tecan®	Celero™ DNA-Seq Library Prep Kit		X			
Oxford Nanopore Technologies (ONT)	Ligation Sequencing Kit V14 Protocol	In Dev	In Dev			X

## DNA library preparation kits (continued)

Vendor	Library Prep Kit	Sciclone™ G3 NGSx	Sciclone™ G3 NGSx iQ™	Zephyr™ G3 NGS	Fontus™ NGS 96	BioQule™
PacBio®	HiFi SMRTbell® Libraries with SMRTbell Express Template Prep Kit 2.0	X	X			
PacBio®	HiFi SMRTbell® Libraries with SMRTbell Prep Kit 3.0	<a href="#">X</a>				
PacBio®	Multiplexed Microbial Library Prep with SMRTbell® Express Template Prep Kit 2.0	X	X			
QIAGEN®	QIAGEN® Targeted DNA v3 Panel	X	X			
QIAGEN®	QIAseq® FX DNA Library Kit	X				
QIAGEN®	QIAseq® Normalizer Kit	In Dev				
Revvity	NEXTFLEX™ Cell Free DNA-Seq Kit 2.0	X	X	X		
Revvity	NEXTFLEX™ Rapid DNA-Seq Kit 2.0	X	X	X		
Revvity	NEXTFLEX™ Rapid XP V2 DNA-seq kit for Metagenomics	<a href="#">X</a>	X	X	X**	
Revvity	PG-Seq™ Rapid v2 Kit	<a href="#">X</a>				
Roche®	Kapa EvoPlus Kit	X				
Roche®	DNA HyperPlus™ Kit	X	X	X	X**	
Roche®	DNA HyperPrep™ Kit	X	X	X	X**	X
seqWell®	plexWell™ 96	In Dev				
Thermo Fisher®	Ion Express® Plus gDNA Fragment Library Prep Kit	X		X		
Thermo Fisher®	Ion Torrent™ Ion Express® Fragment Library Prep Kit	X		X		
TwinStrand® Bioscience	TwinStrand® Duplex Sequencing: Pre-PCR (Mechanical Shearing)	X				
TwinStrand® Bioscience	TwinStrand® Duplex Sequencing: Pre-PCR (Enzymatic Shearing)	X				
TwinStrand® Bioscience	TwinStrand® Duplex Sequencing: Post-PCR			X		
Twist®	Library Preparation Kit (Mechanical Fragmentation)	X	X			
Twist®	Library Preparation Enzymatic Fragmentation (EF) Kit 2.0	X	X		X**	
Watchmaker Genomics	DNA Library Prep	X				
Watchmaker Genomics	DNA Library Prep with Fragmentation	X				

\* Chemistry tested at customer site

\*\* Water tested

In Dev = in development

## RNA library preparation kits

Vendor	Automated qualified methods	Sciclone G3 NGSx	Sciclone G3 NGSx iQ	Zephyr G3 NGS	Fontus NGS 96
10X Genomics®	Chromium® Single Cell 3' Library Kit v3			X	
Agilent®	SureSelect® Strand-Specific RNA Library Prep Kit		X		
Agilent®	SureSelect® XT HS2 RNA		In Dev		
Illumina®/ Epicentre	Ribozero™/Scriptseq™ Stranded RNA Kit	X			
Illumina®	16S Metagenomic Sequencing Library Preparation				X**
Illumina®	RNA Prep with Enrichment (L) Tagmentation	X	X		
Illumina®	Stranded mRNA Prep, Ligation Kit	<a href="#">[X]</a>	X		
Illumina®	Stranded Total RNA Prep, Ligation with Ribo-Zero Plus	<a href="#">[X]</a>	X*		
Illumina®	TruSeq® RNA Access Kit	X			
Illumina®	TruSeq® RNA Library Prep Kit v2	X	X	X	
Illumina®	TruSeq® Stranded mRNA (poly-A selection) Kit	X	X	X	
Illumina®	TruSeq® Stranded Total RNA™ Kit	X	X	X	
Illumina®	TruSeq® RNA Library Prep Kit v2	X	X	X	
Illumina®	TruSight Oncology 500	X*			
Illumina®	TruSight Oncology 500 HT				In Dev
Lexogen®	autoSENSE™ mRNA-Seq Library Prep Kit for Illumina® Sequencing	X			
New England Biolabs® / NEB®	NEBNext® Enzymatic Methyl-seq Library Prep	X	X		
NEB®	NEBNext® Ultra™ RNA Library Prep Kit for Illumina® Sequencing	X		X	
NEB®	NEBNext® Ultra™ RNA II Library Prep Kit for Illumina® Sequencing	X			
NuGen® / Tecan®	Ovation® Human Blood RNA-Seq Library Systems	X			
NuGen® / Tecan®	Ovation® Human FFPE RNA-Seq Multiplex System	X			
NuGen® / Tecan®	Ovation® RNA-Seq System v2	X			
NuGen® / Tecan®	Ovation® Universal RNA-Seq System (strand-specific)	X			
Oxford Nanopore Technologies (ONT)	PCR Tiling of SARS-CoV-2 Virus with rapid barcoding and Midnight RT PCR Expansion	X			
PacBio®	HiFIViral SARS-CoV-2	<a href="#">[X]</a>			
QIAGEN®	QIAseq® DIRECT SARS-CoV-2	X	X		
QIAGEN®	QIAseq® miRNA	X	X*		
Revvity	NEXTFLEX™ Rapid Directional RNA-Seq Kit with Poly(A) Beads 2.0	X	X	X	
Revvity	NEXTFLEX™ Rapid Directional RNA-Seq Kit 2.0 with RiboNaut™ rRNA Depletion Kit	X	X	X	
Roche®	KAPA mRNA HyperPrep™ Kit	X		X	
Roche®	KAPA RNA HyperPrep™ Kit	X	X	X	
Roche®	KAPA RNA HyperPrep™ Kit with RiboErase™ (HMR)	X		X	
Roche®	Stranded mRNA-Seq Kit	X		X	
Roche®	Stranded RNA-Seq Kit	X		X	
Roche®	Stranded RNA-Seq Kit with RiboErase™	X		X	

## RNA library preparation kits (continued)

Vendor	Automated qualified methods	Sciclone G3 NGSx	Sciclone G3 NGSx iQ	Zephyr G3 NGS	Fontus NGS 96
ScaleBio	ScaleBio Single Cell RNA Sequencing Kit v1.1				Coming
Takara® formerly Clontech	SMART-Seq® HT Kit		NGSx HT		
Takara®	SMART-Seq® HT PLUS Kit			X	
Takara®	SMART-Seq® v4 PLUS Kit	In Dev			
Takara®	SMART-Seq® v4 Ultra® Low Input RNA Kit for Sequencing	X			
Takara®	SMARTer® Stranded Total RNA-Seq Kit v2 Pico Input Mammalian		X		
Takara®	SMARTer® Stranded Total RNA-Seq Kit v3 Pico Input Mammalian			X	
Twist®	cDNA Library Preparation Kit for ssRNA Virus Detection	X	X		
Watchmaker Genomics	Watchmaker mRNA Library Prep Kit	<a href="#">[X]</a>			
Watchmaker Genomics	Watchmaker RNA Library Prep Kit with Polaris Depletion	In Dev			

\* Chemistry tested at customer site

\*\* Water tested

In Dev = in development

## Small RNA library prep kits automated on Revvity liquid handlers

Vendor	Library Prep Kit	Sciclone G3 NGSx	Sciclone G3 NGSx iQ	Zephyr G3 NGS	Fontus NGS 96
Revvity	NEXTFLEX™ Small RNA-Seq Kit v4	<a href="#">[X]</a>	<a href="#">[X]</a>		



Revvity NGS library prep kits

[Buy online](#)



## Exome/Targeted Capture library preparation kits

Vendor	Library Prep Kit	Sciclone G3 NGSx	Sciclone G3 NGSx iQ	Zephyr G3 NGS	Fontus NGS 96
10X Genomics®	Exome Preparation	X			
Agilent®	HaloPlex® Kit	X			
Agilent®	SureSelect® XT HS Target Enrichment Kit	X	X		
Agilent®	SureSelect® XT Reagent Kit	X	X		X**
Agilent®	SureSelect® XT2 Reagent Kit	X	X	X	
Agilent®	SureSelect® QXT Reagent Kit	X		X	
Cergentis® / Solvias®	TLA Targeted Locus Amplification Kit	X			
Integrated DNA Technologies / IDT™	XGen™ Hybridization and Wash Kit	X	X		X**
IDT™	xGen™ Exome Hyb Panel v2	X	X		
Illumina®	COVIDSeq™ Test	X	X	X	
Illumina®	Illumina® DNA Prep with Exome 2.5 Enrichment	X	X		
Illumina®	Nextera® Rapid Capture Exome Kit	X		X	
Illumina®	TruSeq® Cancer HotSpot Panel v2	X			
Illumina®	TruSight® Cancer Kit	X			
Illumina®	TruSeq® Comprehensive Panel v3	X			
Illumina®	TruSeq® Custom Amplicon v1.5 Kit			X	
Illumina®	TruSeq® Custom Amplicon Low Input Kit	X		X	
Illumina®	TruSeq® Exome Kit	X			
Illumina®	TruSeq® Focus Panel	X			
Illumina®	TruSight® Rapid Capture Kit	X			
Illumina®	TruSeq® Rapid Exome Kit	X			
New England Biolabs® / NEB®	NEBNext® Direct® Cancer HotSpot Panel	X			
NuGen® / Tecan®	Ovation® Target Enrichment System	X			
Revvity	NEXTFLEX™ Neo NGS RUO Panel 1				X**
Roche®	KAPA HyperCap Workflow v3	<a href="#">X</a>	<a href="#">X</a>		
Roche®	SeqCap® EZ Kit	X			
Twist®	Fast Hybridization Target Enrichment	X	X		X**
Twist®	Human Core Exome Kit	X	X		X**
Twist®	TWIST® Custom Panels	X	X		

\* Chemistry tested at customer site

\*\* Water tested

## Metagenomics library preparation kits

Vendor	Library Prep Kit	Sciclone G3 NGSx	Sciclone G3 NGSx iQ	Zephyr G3 NGS	Fontus NGS 96
NuGen® / Tecan®	Ovation® Ultralow Methyl-Seq Library System	X			
Revvity	NEXTFLEX™ 16S V1-V3 Amplicon-Seq Kit	X			
Revvity	NEXTFLEX™ 16S V3-V4 Amplicon-Seq Kit 2.0	X	X		In Dev
Revvity	NEXTFLEX™ 16S V4 Amplicon-Seq Kit 2.0	X			
Revvity	NEXTFLEX™ Rapid XP DNA-Seq Kit v2	<a href="#">X</a>	X	X	X*

\* Water tested

In Dev = in development

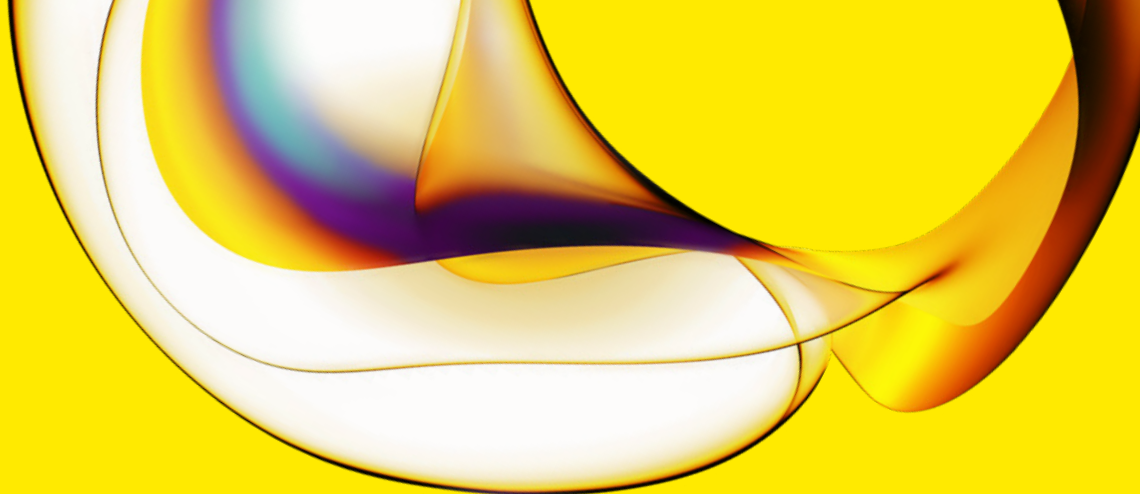
## Supplementary applications & protocols

Vendor	Library Prep Kit	Sciclone G3 NGSx	Zephyr G3 NGS	Fontus NGS 96
Beckman Coulter®	SPRI® Purification/Bead Size Selection Protocol	X	X	X
NuGen® / Tecan®	Ovation® FFPE WTA System	X		
NuGen® / Tecan®	Ovation® Pico WTA System v2	X		
Revvity	NEXTFLEX™ Poly(A) Beads 2.0	X	X	
Revvity	Normalization & Pooling Protocol	X	X	
Roche®	KAPA Library Quantification Kit	X	X	
Roche®	KAPA SYBR® FAST qPCR Kit	X		
Thermo Fisher®	PicoGreen® Quantitation	X	X	
Thermo Fisher®	Qubit® Quantitation		X	
Thermo Fisher®	Quant-iT® Quantitation		X	



Learn more about our automated liquid handling solutions.

[Learn more](#)



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

revvity