

NEXTFLEX[®] RNA-Seq 2.0 Unique Dual Index Barcodes

(Compatible with Illumina[®] and Element[®]
platforms)

KIT CONTAINS : 8, 24, or 96 BARCODES | 16, 48, or 192 RXNS

USER MANUAL FOR :
#NOVA-512920-eval16
#NOVA-512920-eval48
#NOVA-512920
#NOVA-512921
#NOVA-512922
#NOVA-512923

NEXTFLEX® RNA-Seq 2.0 Unique Dual Index Barcodes

GENERAL INFORMATION _____	3
Product Overview	3
Kit Overview	3
Kit Contents, Storage & Shelf Life	3
Warnings and Precautions	4
SAMPLE PREP WORKFLOW _____	5
NEXTFLEX® Rapid Directional RNA-Seq Kit 2.0 Sample Preparation Flow Chart	5
UDI Barcoded Adapter Plate Format _____	7
APPENDIX A _____	7
Oligonucleotide Sequences.....	7

This product is for research use only.

Not for use in diagnostic procedures.

This manual is proprietary to Revvity and intended only for customer use in connection with the product(s) described herein and for no other purpose. This document and its contents shall not be used or distributed for any other purpose without the prior written consent of Revvity. Follow the protocol included with the kit.

Revvity, NEXTFLEX®, NextPrep™, NextPrep-Mag™ are trademarks or registered trademarks of Revvity. All other brands and names contained herein are the property of their respective owners.

GENERAL INFORMATION

Product Overview

The NEXTFLEX® RNA-Seq 2.0 Unique Dual Index Barcodes are designed to prepare multiplexed single and paired-end cDNA libraries from total RNA, mRNA-enriched RNA or rRNA-depleted RNA for sequencing using Illumina and Element platforms. The index and flow cell binding sequences contained within the NEXTFLEX® RNA-Seq 2.0 Unique Dual Index Barcodes attach to the sample insert during adapter ligation. Sample pooling with NEXTFLEX® RNA-Seq 2.0 Unique Dual Index Barcodes allows the user to multiplex up to 384 samples when used in conjunction with other available sets of NEXTFLEX® RNA-Seq 2.0 Unique Dual Index Barcodes. NEXTFLEX® RNA-Seq 2.0 Unique Dual Index Barcodes have been performance-verified to be used in conjunction with the NEXTFLEX Rapid Directional RNA-Seq Kit 2.0.

Uniquely dual-indexed libraries are libraries prepared with adapters containing two eight base indexes:

Index 1 (P7 Index) adjacent to the P7 strand, and Index 2 (P5 Index) adjacent to the P5 strand. None of the indexes found on any given NEXTFLEX RNA-Seq 2.0 Unique Dual Index Barcode are used throughout the entire set, which prevents mis-assigned reads from appearing in final data sets.

Each lot of the NEXTFLEX® RNA-Seq 2.0 Unique Dual Index Barcodes is functionally validated and tested for index purity by sequencing.

Kit Overview

This NEXTFLEX® RNA-Seq 2.0 Unique Dual Index Barcodes Kit contains 8, 24, or 96 uniquely dual-indexed barcoded DNA adapters in tubes or plate format for a total of 16, 48, or 192 reactions.

Kit Contents, Storage & Shelf Life

Note: The 16-reaction kit contains UDI barcodes 1-8, the 48-reaction kit contains UDI barcodes 1-24, and the 192-reaction kit contains UDI barcodes 1-96, 97-192, 193-288, or 289-384.

It is recommended that UDI barcodes are stored at -20°C. The shelf life of each reagent is at least 1 year when stored properly.

Kit Contents	Cap Color	Amount (16 rxn / 48 rxn / 192 rxn)	Storage Temp.
NEXTFLEX® RNA-Seq 2.0 Unique Dual Index Barcodes* (6.25 µM)	BLUE CAP or PLATE	5 µL each	-20°C
NEXTFLEX® Primer Mix 2.0** (50 µM)	GREEN CAP	32 µL / 96 µL / 384 µL	-20°C

*These Unique Dual Index Barcodes are supplied in duplex form. Do not heat the adapters above room temperature. The UDI barcodes in the 16 and 48 rxn kits will be in vials with blue caps. The 192 rxn kits will be plated column-wise (1-8, 9-16, etc.).

** The Primer Mix 2.0 is only intended for use with the NEXTFLEX® Rapid Directional RNA-Seq Kit 2.0; if the primer mix is to be used with any other RNA-Seq Kits of choice, additional dilutions may be required.

Warnings and Precautions

- We strongly recommend that you read the following warnings and precautions. Periodically, optimizations and revisions are made to the components and manual. Therefore, it is important to follow the protocol included with the kit. If you need further assistance, you may contact your local distributor, or contact us at <https://www.revivity.com/contact-us/technical-support> and choose the “Next Gen Sequencing” category.
- Do not use the kit past the expiration date.
- Ensure pipettes are properly calibrated as library preparations are highly sensitive to pipetting error.
- Do not heat the NEXTFLEX® RNA-Seq 2.0 Unique Dual Index Barcodes above room temperature.
- Once plate has thawed, spin for one minute before use. This is to ensure all liquid settles to the bottom of the plate.
- NEXTFLEX® RNA-Seq 2.0 Unique Dual Index Barcodes have been performance-verified to be used in conjunction with the NEXTFLEX Rapid Directional RNA-Seq Kit 2.0.
- The plate seal is intended to be pierced. Do not peel the plate seal from the plate, doing so can easily lead to cross-contamination. Additional thermal heat seals may be applied upon one another to re-seal plate.
- Before use, carefully mix adapters by pipetting up and down several times using a multi-channel pipette with barrier tip. NEVER mix plates by vortexing. Placing a plate on a vortexer to mix samples or barcodes has been proven to result in cross-contamination, even if the plate appears to be securely sealed.
- Try to maintain a laboratory temperature of 20°–25°C (68°–77°F).
- The NEXTFLEX® Primer Mix 2.0 is only compatible with the NEXTFLEX® Rapid Directional RNA-Seq Kit 2.0 at the concentration it is supplied in. If the NEXTFLEX® RNA-Seq 2.0 Unique Dual Index Barcodes are used with other NEXTFLEX® kits, both the Barcodes and Primer Mix must be diluted prior to use. Make sure to verify that the Primer Mix concentration is compatible at its stock concentration or dilute with nuclease-free water prior to use. The NEXTFLEX® Primer Mix 2.0 should always be used with the NEXTFLEX® RNA-Seq 2.0 Unique Dual Index Barcodes for PCR amplification even if alternative primers are supplied with your kit of choice. Inadvertent use of an incorrect primer sequence can potentially result in elimination of the index. Primer Mix included with the NEXTFLEX® library prep kits should always be used instead of the Primer Mixes that come with the NEXTFLEX® barcodes. For additional information, please contact <https://www.revivity.com/contact-us/technical-support>.

Version	Date	Description
V20.03	March	Manual Layout
V20.11	November 2020	UDI Index Change
V20.03	March 2020	Manual Layout
V23.10	October 2023	Revvity Rebrand

SAMPLE PREP WORKFLOW

NEXTFLEX® Rapid Directional RNA-Seq Kit 2.0 Sample Preparation Flow Chart

FRAGMENTATION USING POLY (A) ENRICHED, rRNA DEPLETED, or TOTAL RNA



FIRST STRAND SYNTHESIS



SECOND STRAND SYNTHESIS



BEAD CLEANUP (Optional Stopping Point)

ADD 'A'



ADD ADAPTERS



LIGATION



BEAD CLEANUP (Optional Stopping Point)

PCR



BEAD CLEANUP

UDI Barcoded Adapter Plate Format

Representative plate layout of UDI Barcoded Adapters 1-96, 97-192, 193-288, and 289-384.

	1	2	3	4	5	6	7	8	9	10	11	12
A	1	9	17	25	33	41	49	57	65	73	81	89
B	2	10	18	26	34	42	50	58	66	74	82	90
C	3	11	19	27	35	43	51	59	67	75	83	91
D	4	12	20	28	36	44	52	60	68	76	84	92
E	5	13	21	29	37	45	53	61	69	77	85	93
F	6	14	22	30	38	46	54	62	70	78	86	94
G	7	15	23	31	39	47	55	63	71	79	87	95
H	8	16	24	32	40	48	56	64	72	80	88	96

1-96 in columns

	1	2	3	4	5	6	7	8	9	10	11	12
A	97	105	113	121	129	137	145	153	161	169	177	185
B	98	106	114	122	130	138	146	154	162	170	178	186
C	99	107	115	123	131	139	147	155	163	171	179	187
D	100	108	116	124	132	140	148	156	164	172	180	188
E	101	109	117	125	133	141	149	157	165	173	181	189
F	102	110	118	126	134	142	150	158	166	174	182	190
G	103	111	119	127	135	143	151	159	167	175	183	191
H	104	112	120	128	136	144	152	160	168	176	184	192

97-192 in columns

	1	2	3	4	5	6	7	8	9	10	11	12
A	193	201	209	217	225	233	241	249	257	265	273	281
B	194	202	210	218	226	234	242	250	258	266	274	282
C	195	203	211	219	227	235	243	251	259	267	275	283
D	196	204	212	220	228	236	244	252	260	268	276	284
E	197	205	213	221	229	237	245	253	261	269	277	285
F	198	206	214	222	230	238	246	254	262	270	278	286
G	199	207	215	223	231	239	247	255	263	271	279	287
H	200	208	216	224	232	240	248	256	264	272	280	288

193-288 in columns

	1	2	3	4	5	6	7	8	9	10	11	12
A	289	297	305	313	321	329	337	345	353	361	369	377
B	290	298	306	314	322	330	338	346	354	362	370	378
C	291	299	307	315	323	331	339	347	355	363	371	379
D	292	300	308	316	324	332	340	348	356	364	372	380
E	293	301	309	317	325	333	341	349	357	365	373	381
F	294	302	310	318	326	334	342	350	358	366	374	382
G	295	303	311	319	327	335	343	351	359	367	375	383
H	296	304	312	320	328	336	344	352	360	368	376	384

289-384 in columns

APPENDIX A

Oligonucleotide Sequences

NEXTFLEX	Sequence (5' → 3')
PCR Primer 1	AATGATACGGCGACCACCGAGATCTACAC
PCR Primer 2	CAAGCAGAAGACGGCATAACGAGAT
NEXTFLEX® UDI Barcode	AATGATACGGCGACCACCGAGATCTACACXXXXXXXX ¹ ACACTCTTCCCTA CACGACGCTCTCCGATCT GATCGGAAGAGCACACGTCTGAACTCCAGTCACXXXXXXXX ² ATCTCGTATG CCGCTCTTCTGCTTG

XXXXXXXX¹ denotes the P5 index region of adapter. The index sequences contained in each adapter are listed below.

XXXXXXXX² denotes the P7 index region of the adapter. The index sequences contained in each adapter are listed below.

When entering index sequences for the Illumina MiniSeq®, NextSeq®, HiSeq® 3000 or HiSeq® 4000 platforms, enter the P5 Index Reverse Complement. For all other Illumina platforms, enter the P5 Index in the first column.

	P5 Index	P5 Index Reverse Complement	P7 Index
UDI 1	AATAACGT	ACGTTATT	AATCGTTA
UDI 2	TTCTTGAA	TCAAGAA	GTCTACAT
UDI 3	GGCAGATC	GATCTGCC	CGCTGCTC
UDI 4	CTATGTTA	TAACATAG	GATCAACA
UDI 5	GTTGACGC	GCGTCAAC	CGAAGGAC
UDI 6	ATCTACGA	TCGTAGAT	GATGCCGG
UDI 7	CTCGACAG	CTGTCGAG	CTACGAAG
UDI 8	GAGGCTGC	GCAGCCTC	GATGCGTC
UDI 9	CCTCGTAG	CTACGAGG	CTACGGCA
UDI 10	CATAGGCA	TGCCTATG	GATTCCCTT
UDI 11	AGATGAAC	GTTTCATCT	CTACTCGA
UDI 12	CCGAGTAT	ATACTCGG	GATTCCGAG
UDI 13	AATATTGA	TCAATATT	AATCGGCG
UDI 14	GTATACCG	CGGTATAC	TTCGCCGA
UDI 15	GATCCAAC	GTTGGATC	CTGGCCTC
UDI 16	AGATACGC	GCGTATCT	GAACCTAT

UDI 17	GGTATCTT	AAGATACC	CGTATTGG
UDI 18	CCTCTGGC	GCCAGAGG	GAAGCACA
UDI 19	CCATTGTG	CACAATGG	CTTAATAC
UDI 20	ACTACGGT	ACCGTAGT	GAAGTCTT
UDI 21	AAGTGCTA	TAGCACTT	GAAGAGGC
UDI 22	GCCGAACG	CGTTCGGC	CGGATAAC
UDI 23	TGTCCACG	CGTGGACA	GAATCTGG
UDI 24	GACACACT	AGTGTGTC	CTGATTGA
UDI 25	AATATGCT	AGCATATT	AATCCGTT
UDI 26	TTCTCATA	TATGAGAA	TGCGTACA
UDI 27	TCTGTGAT	ATCACAGA	GAATCAAT
UDI 28	CCGAECTT	AAGTTCGG	TGAGTCAG
UDI 29	GTCTAACA	TGTTAGAC	GAATGCTC
UDI 30	GACGCCAT	ATGGCGTC	GAATATCC
UDI 31	GCCAATGT	ACATTGGC	CTTATGAA
UDI 32	CCAACGTC	GACGTTGG	TCGGCACC
UDI 33	GTAGATAA	TTATCTAC	AAGAAGCG
UDI 34	CTTACGGC	GCCGTAAG	CTCACGAT
UDI 35	CCAAGTGC	GCACTTGG	TCGGTCGA
UDI 36	CTAACTCA	TGAGTTAG	TCGGTAAG
UDI 37	AATATCTG	CAGATATT	AAGATACA
UDI 38	TTATATCA	TGATATAA	GTCGCTGT
UDI 39	CTGCGGAT	ATCCGCAG	TCGGATGT
UDI 40	GCGGCTTG	CAAGCCGC	CGAGCCGG
UDI 41	GAGTTGAT	ATCAACTC	CGATTATC
UDI 42	GCACTGAG	CTCAGTGC	TCGAAGCT
UDI 43	GACCACCT	AGGTGGTC	CTATCATT
UDI 44	TGGCTAGG	CCTAGCCA	CGCGCCAA
UDI 45	CCTACCGG	CCGGTAGG	CGAACGGA
UDI 46	GGAGGATG	CATCCTCC	CTACTGAC
UDI 47	CGCTGAAT	ATTCAGCG	TCTTAAGT
UDI 48	TGTGACGA	TCGTCACA	TTAGAGTC
UDI 49	AATAGATT	AATCTATT	AAGACGAA
UDI 50	TTAGCGCA	TGCGCTAA	TTATTATG
UDI 51	GCGGCCGT	ACGGCCGC	CGCTATTA
UDI 52	CAGTAACC	GGTACTG	TCTATCAG
UDI 53	GCCTAGTA	TACTAGGC	CGGTGGTA
UDI 54	CACGGCGC	GCGCCGTG	TCACCAAT

UDI 55	GGTGCAGA	TCTGCACC	CTGGAAGC
UDI 56	GTAAGTGC	CGAGTTAC	TCCTCGAT
UDI 57	CAGCCAGT	ACTGGCTG	AAGAGAGC
UDI 58	CGTCAACC	GGTTGACG	TCAACGAG
UDI 59	GCCGGCGA	TCGCCGGC	TGCGAGAC
UDI 60	GCCTCCGG	CCGGAGGC	CCTGGTGT
UDI 61	AATAGTCC	GGACTATT	AAGTAAGT
UDI 62	TTAGACGT	ACGTCTAA	TGACTGAA
UDI 63	GTGGACTA	TAGTCCAC	AAGACTGT
UDI 64	CACGGACG	CGTCCGTG	CAATGATG
UDI 65	CACTAGAG	CTCTAGTG	CACAGTAA
UDI 66	GCAGATGG	CCATCTGC	TGGTCATT
UDI 67	CTCTACG	CGTGAGAG	CAACCGTG
UDI 68	GGAATCAC	GTGATTCC	TGGTGAC
UDI 69	CGTTGACG	CGTCAACG	CCACAATG
UDI 70	CATCAGGT	ACCTGATG	TGTGTGCC
UDI 71	CGTTGTAA	TTACAACG	CACCACGG
UDI 72	GGCACGGT	ACCGTGCC	TGTGTTAA
UDI 73	AATAGCAA	TTGCTATT	AAGTTATC
UDI 74	TGATCGGT	ACCGATCA	GTACAGCT
UDI 75	AGTAGTAT	ATACTACT	CAACTGCT
UDI 76	GTTAGAGG	CCTCTAAC	CATGATGA
UDI 77	CCTTACAG	CTGTAAGG	TGACTACT
UDI 78	GTACATTG	CAATGTAC	CAGAAGAT
UDI 79	GGAGACCA	TGGTCTCC	TGAGGCGC
UDI 80	CGAACACC	GGTGTTCCG	CAGGTTCC
UDI 81	GAGAACAA	TTGTTCTC	TGAACAGG
UDI 82	TGTGAATC	GATTCACA	CAGTGTGG
UDI 83	GGTTAAGG	CCTTAACC	TTCCACCA
UDI 84	AGACCGCA	TGCGGTCT	CCGCTGTT
UDI 85	AATACAGG	CCTGTATT	AAGTTGGA
UDI 86	TGATGGCC	GGCCATCA	GGACAACG
UDI 87	TGTCACCT	AGGTGACA	TTCGAACC
UDI 88	GCTTCGGC	GCCGAAGC	CAGACCAC
UDI 89	CCAGTGGT	ACCACTGG	TTCTGGTG
UDI 90	GCACACGC	GCGTGTGC	CAATCGAA
UDI 91	GTCACGTC	GACGTGAC	AAGTACAG
UDI 92	GCAGCTCC	GGAGCTGC	CCGTGCCA

UDI 93	CATGCAGC	GCTGCATG	CATTGCAC
UDI 94	ACGATTGC	GCAATCGT	TTACCTGG
UDI 95	GACATTCG	CGAATGTC	CTGCAACG
UDI 96	GCGAATAC	GTATTTCG	TACTGTTA
UDI 97	AATTAATG	CATTAATT	AAGTGTAT
UDI 98	TTAATTCC	GGAATTAA	GTCCACTC
UDI 99	GAAGGAAC	GTTCTTC	CTTGCTAT
UDI 100	CAGGCATA	TATGCCTG	TACATAGA
UDI 101	CCAATACT	AGTATTGG	TAGCCGAT
UDI 102	TCGCGCAT	ATGCGCGA	CGATCCAC
UDI 103	GTGTGAAC	GTTACAC	TAGCGTTG
UDI 104	GGTGGCAC	GTGCCACC	CTCATCAC
UDI 105	TAGGTGCT	AGCACCTA	TATGCGGT
UDI 106	GAGCGGTT	AACCGCTC	TAACTCGC
UDI 107	AGGATTAG	CTAATCCT	CGTACGTT
UDI 108	GTGAGCCA	TGGCTCAC	TAAGTACC
UDI 109	AATTATGC	GCATAATT	AAGTCGTG
UDI 110	TTACGCCG	CGGCGTAA	TTCAGAAC
UDI 111	GTAGTGAT	ATCACTAC	GTTATATA
UDI 112	TGTTGCA	TGCAACCA	ACCGCTAT
UDI 113	GTCCGACC	GGTCGGAC	ACCGTCCT
UDI 114	GCCAGCCG	CGGCTGGC	TGTCTAAC
UDI 115	CTCGTGTC	GACACGAG	ACCGAGGT
UDI 116	GCCGTGGC	GCCACGGC	ACCGATTA
UDI 117	GTAGCCAC	GTGGCTAC	GTTCTACT
UDI 118	GACTTATA	TATAAGTC	ACCTGACT
UDI 119	GAACGTCG	CGACGTTT	GGTAATCG
UDI 120	ATGCGTAG	CTACGCAT	ACCTTAGA
UDI 121	AATTAGAT	ATCTAATT	AAGGAGTT
UDI 122	TTACGATA	TATCGTAA	TGAAGCCA
UDI 123	CTGTACAA	TTGTACAG	GTGTTGTA
UDI 124	CGTGGTTG	CAACCACG	ACCACACG
UDI 125	TGTAGCGG	CCGCTACA	ACCAGGAC
UDI 126	GTGTCATG	CATGACAC	TTGGCAGG
UDI 127	GAACCGGT	ACCGGTTT	ACCATTAA
UDI 128	CACTCTGA	TCAGAGTG	TTGTAGAT
UDI 129	TACACGTT	AACGTGTA	ACCATATC
UDI 130	GAGTGCCT	AGGCACTC	TTGGTGGC

UDI 131	GACGGTCC	GGACCGTC	ACCAACAT
UDI 132	CTGATGCT	AGCATCAG	GTCTGTGC
UDI 133	AATTACCA	TGGTAATT	AAGGCAAT
UDI 134	GTCCTGTT	AACAGGAC	TGCATTGC
UDI 135	GTTAACAG	CTGTTAAC	ACGCCACT
UDI 136	AGCCTGTA	TACAGGCT	AAGTCTCC
UDI 137	ACGTTGCT	AGCAACGT	GGATCTCT
UDI 138	GGTGTGTC	GCAACACC	ACGCGATC
UDI 139	TGGTACTA	TAGTACCA	GGATAATA
UDI 140	CATCATTG	CAATGATG	ACGCTTAT
UDI 141	GACGATTG	CAATCGTC	TTAGGTTG
UDI 142	CGGCCGAA	TTCGGCCG	ACGCACAA
UDI 143	CTGCCGCC	GGCGGCAG	TGAATATA
UDI 144	ACTAACCG	CGGTTAGT	TGCTCCGC
UDI 145	AATTGAGA	TCTCAATT	AAGCAATA
UDI 146	TTCAAGAG	CTCTTGAA	TTATGTAT
UDI 147	GTCCGCAA	TTGCGGAC	ACGGCAGA
UDI 148	CAAGGTAG	CTACCTTG	GGACTCTG
UDI 149	AGCGACTC	GAGTCGCT	GTACGTAC
UDI 150	GTAGTATC	GATACTAC	GGAAGGTA
UDI 151	GTGGCGAA	TTCGCCAC	ACGTCCAT
UDI 152	ACGTGCGC	GCGCACGT	ACGTGTTG
UDI 153	GAAGACAG	CTGTCTTC	TGAAGAAT
UDI 154	GCGCAGAT	ATCTGCGC	ACGTAGTC
UDI 155	AGAGACAT	ATGTCTCT	AAGGATAA
UDI 156	CACTTGTT	AACAAGTG	GGCGAGGA
UDI 157	AATTGGTC	GACCAATT	AACAAGGC
UDI 158	TTAACTACT	AGTGTTAA	TTGGTCCG
UDI 159	GCTCAGCG	CGCTGAGC	ACGAGCCT
UDI 160	GTTCAGAC	GTCTGAAC	GGAGATTC
UDI 161	CAGTTGCG	CGCAACTG	TGCGCGCT
UDI 162	GCAATATG	CATATTGC	ACGATCTA
UDI 163	CGCAACGT	ACGTTGCG	TGCTGAGG
UDI 164	GGTGAGAT	ATCTCACC	ACTCCTCC
UDI 165	CCAGCTGA	TCAGCTGG	ACTCTACG
UDI 166	GTTGAGCA	TGCTCAAC	TTCGTTCT
UDI 167	CTGTAGTG	CACTACAG	TGCAGTCG
UDI 168	GTGCATCC	GGATGCAC	TGCCGTAA

UDI 169	AATTCACT	AGTGAATT	AACAGTTG
UDI 170	TTCAGTGA	TCACTGAA	TGTTAACA
UDI 171	AGTCGCGA	TCGCGACT	ACTTCCTG
UDI 172	CCATATCG	CGATATGG	TGAATGCG
UDI 173	AAGCATGG	CCATGCTT	ACTTGTAC
UDI 174	CGTGTCTA	TAGACACG	ACTACTTA
UDI 175	CCGTCTAA	TTAGACGG	CTCTTCGT
UDI 176	CGAGAGTG	CACTCTCG	ACTAGCTC
UDI 177	CGAGTACG	CGTACTCG	TTATCAAC
UDI 178	GTGTCTCA	TGAGACAC	TGCTTGTC
UDI 179	CCATCGCT	AGCGATGG	AAGGTAGG
UDI 180	CTTCGCTC	GAGCGAAG	TGGTATGG
UDI 181	AATTCTAA	TTAGAATT	AACACATA
UDI 182	TTCATCTG	CAGATGAA	TGGTGTCT
UDI 183	ACAGTGAA	TTCACTGT	ACACTAAC
UDI 184	AGATGTGA	TCACATCT	TTGTGTTC
UDI 185	GGCAACTG	CAGTTGCC	ACACACCT
UDI 186	CAGCGTCT	AGACGCTG	GGTGTCCG
UDI 187	CTGTCGGT	ACCGACAG	ACACATTC
UDI 188	TCGGCGTT	AACGCCGA	TTGCTTAA
UDI 189	CACACGCG	CGCGTGTG	ACAGCCTT
UDI 190	AGCCAACA	TGTTGGCT	TTGCAGTA
UDI 191	AGCAAGTT	AACTTGCT	ACAGGCAG
UDI 192	ACGCGTCA	TGACGCGT	TTGCCATC
UDI 193	AATGAACC	GGTTCATT	ATAATGTA
UDI 194	GGAAGTGG	CCACTTCC	TCGTACCG
UDI 195	CGCTAAGC	GCTTAGCG	GTTATTAT
UDI 196	GTCATAGC	GCTATGAC	ACAGAAGC
UDI 197	GTGACCAT	ATGGTCAC	TTGATAAT
UDI 198	GAAGTTGA	TCAACTTC	ACATCGGA
UDI 199	GCGGTTAA	TTAACCGC	ACCTCTCA
UDI 200	CGCCTCTT	AAGAGGCG	GTGAGTGT
UDI 201	GTTCTGGA	TCCAGAAC	ACATCACC
UDI 202	ACTGCATC	GATGCAGT	CTCGCTTG
UDI 203	GGTTAGCC	GGCTAACC	ACATGGAT
UDI 204	TCCACAGC	GCTGTGGA	TCAGGATC
UDI 205	AATGTGGC	GCCACATT	ATAAGGCT
UDI 206	GTCACAAG	CTTGTGAC	TCTTACTC

UDI 207	CGTAGATA	TATCTACG	ACATGATA
UDI 208	TGCGGCTA	TAGCCGCA	TTGCCTCT
UDI 209	ACAACTCT	AGAGTTGT	ACATTCTC
UDI 210	GATCGCAG	CTGCGATC	ACATATGT
UDI 211	TGGAACAT	ATGTTCCA	GTTGGTAG
UDI 212	CACATACC	GGTATGTG	ACAACCGC
UDI 213	GCTAAGAA	TTCTTAGC	AAGCAGAC
UDI 214	CTTAACCA	TGGTTAAG	ACAAGGTG
UDI 215	GCCTTCAA	TTGAAGGC	ACAAGAGT
UDI 216	AACTGGAA	TTCCAGTT	TTGGAATT
UDI 217	AATGGAAG	CTTCCATT	ATAACAGA
UDI 218	TTCAATTC	GAATTGAA	TATTGTCG
UDI 219	CGTCCTTA	TAAGGACG	ACAATGCC
UDI 220	CACCGGTA	TACCGGTG	TATGGTTC
UDI 221	AGGAACCA	TGGTTCTT	AGCCACAG
UDI 222	GATCTTCC	GGAAGATC	GCTATCGA
UDI 223	CGGCTTGT	ACAAGCCG	AGCGCTCG
UDI 224	GCCGACAC	GTGTCGGC	TGCGCTTC
UDI 225	CGGCGTAC	GTACGCCG	TCTACGCC
UDI 226	GCCTCATT	AATGAGGC	TATCACAA
UDI 227	CGGCACGA	TCGTGCCG	AGCTTGGT
UDI 228	GCCTGTAG	CTACAGGC	AGCTTATG
UDI 229	AATGGCCT	AGGCCATT	ATAACGAC
UDI 230	TTCCAGGC	GCCTGGAA	GATGGAGT
UDI 231	TTAACCTC	GAGGTAA	TGATCTAA
UDI 232	CGTTAGGA	TCCTAACG	AGCTATAT
UDI 233	GCCATAAT	ATTATGGC	AGCAGAGA
UDI 233	GCCATAAT	ATTATGGC	AGCAGAGA
UDI 234	GCCAAGCC	GGCTTGGC	GACGTAAG
UDI 235	CTCGGTCA	TGACCGAG	AGGCCTGA
UDI 236	ACGTAACA	TGTTACGT	CCGCCTCG
UDI 237	GCGTCCTC	GAGGACGC	AGGCGAAT
UDI 238	CTACAAGT	ACTTGTAG	TTGTTCTT
UDI 239	GCGTGTCC	GGACACGC	TACTCCTG
UDI 240	GCGACTGA	TCAGTCGC	GACTGGCG
UDI 241	AATGCTTG	CAAGCATT	ATATATAC
UDI 242	TTAATCAA	TTGATTAA	TAGATCGG
UDI 243	CTCTGGAC	GTCCAGAG	AGGCAAGC

UDI 244	GCTCCAGT	ACTGGAGC	TCCAGGTA
UDI 245	CGCGTTAT	ATAACGCG	AGGTCAAG
UDI 246	GCTCAATC	GATTGAGC	TACACTGG
UDI 247	CGCATGTG	CACATGCG	TCCTTGCG
UDI 248	GCTGCCAA	TTGGCAGC	AGGACCTC
UDI 249	CGAACTAT	ATAGTTCG	TAAGCATT
UDI 250	GCTGTCTC	GAGACAGC	AGGATATT
UDI 251	CGACGGTT	AACCGTCG	TCATTCGG
UDI 252	GCTTCACA	TGTGAAGC	AGTCCGAC
UDI 253	AATGCGCA	TGCGCATT	ATATTACA
UDI 254	TGAATCGT	ACGATTCA	TCTCCGTG
UDI 255	CTAACCCAG	CTGGTTAG	TAATAGGA
UDI 256	GCTTGTGA	TCACAAGC	TTGTAAGA
UDI 257	CTGGCACT	AGTGCCAG	GAATGTAA
UDI 258	GCACGGTC	GACCGTGC	AGTCAATT
UDI 259	CTTGATT	AATCCAAG	GAACAATT
UDI 260	GCACATCA	TGATGTGC	AGTGGTCA
UDI 261	CTTCAATA	TATTGAAG	TACAATAT
UDI 262	TTACTGGT	ACCAGTAA	AAGCGCGT
UDI 263	CTGAGTCC	GGACTCAG	GAACCTTA
UDI 264	GCATACTT	AAGTATGC	AGTGTACT
UDI 265	AATGCCAC	GTGGCATT	TATACCAT
UDI 266	GGCCGTTG	CAACGGCC	ATCGTGGA
UDI 267	GCATAGAC	GTCTATGC	GCAAGACA
UDI 268	CATTGCGC	GGCGAATG	TACAGGCC
UDI 269	GGCCGCGT	ACGCGGCC	AGTTCGGT
UDI 270	CCTGAAGG	CCTTCAGG	TCAATTAT
UDI 271	GGCCTTCA	TGAAGGCC	AGTTCATA
UDI 272	GGCTTACT	AGTAAGCC	TCACCGGC
UDI 273	TCTGCTCG	CGAGCAGA	GACACCGA
UDI 274	GGCACTCC	GGAGTGCC	AGTTGGCT
UDI 275	CATGTGCT	ACGACATG	TACGTCTC
UDI 276	TACCTATT	AATAGGTA	CCGGTACA
UDI 277	AATCTGAA	TTCAGATT	TATACATC
UDI 278	TGCGAAGT	ACTTCGCA	ATCGTTAG
UDI 279	CAAGGAGA	TCTCCTTG	GAATTACG
UDI 280	GGTCTTAG	CTAAGACC	AGTACTGC
UDI 281	CAAGTAAT	ATTACTTG	TACTAACG

UDI 282	GGTCAGTA	TACTGACC	AGTATGTC
UDI 283	CCAGTCAC	GTGACTGG	TAAGTGTG
UDI 284	GGTCATGT	ACATGACC	ATATAAGG
UDI 285	CAACGACC	GGTCGTTG	ATAGAATA
UDI 286	GGTTCTTG	CAAGAACC	TCTAGAGA
UDI 287	CAACATAC	GTATGTTG	ATAGGCCA
UDI 288	GGTTGATT	AATCAACC	ATAGCGGT
UDI 289	AATAATAG	CTATTATT	ATATCCTA
UDI 290	GTCGGATA	TATCCGAC	GATAGGAT
UDI 291	CACATTGT	ACAATGTG	ATACTGCG
UDI 292	GGTTACAA	TTGTAACC	TCTTCCGA
UDI 293	CACGAATT	AATTCGTG	ATACCTAT
UDI 294	GGTATGCA	TGCATACC	TCGTTATA
UDI 295	CAGGACCT	AGGTCCTG	TAATTAGT
UDI 296	GGACTGGC	GCCAGTCC	ATTATTCCG
UDI 297	CAGAGCGA	TCGCTCTG	TAATAATC
UDI 298	GGAGCTAG	CTAGCTCC	ATTATCGC
UDI 299	CAGCTTAG	CTAAGCTG	TAATATAG
UDI 300	CCGCCAAC	GTTGGCGG	ATTAGACA
UDI 301	CCTATGCG	CGCATAGG	ATAGATCT
UDI 302	GGATCCGA	TCGGATCC	TAGAGCTC
UDI 303	GGATTGTT	AACAATCC	ATTACAAT
UDI 304	CAGGTGAA	TTCACCTG	ATTGAAGT
UDI 305	GGAACATT	AATGTTCC	ATTGATTC
UDI 306	CAAGACTC	GAGTCTTG	TAACTAAG
UDI 307	GTGCGGCG	CGCCGCAC	ATTGACAA
UDI 308	CAATCTTC	GAAGATTG	ATTGTGTT
UDI 309	GTGCTCAG	CTGAGCAC	ATTGCTGA
UDI 310	CAACGCAT	ATGCGTTG	TACATCCT
UDI 311	GTGTAGGC	GCCTACAC	ATTCACGG
UDI 312	CAACTATA	TATAGTTG	TAATTGAC
UDI 313	GCATTGCA	TGCAATGC	ATAGTCTG
UDI 314	CTTACATG	CATGTAAG	TATAAGAC
UDI 315	GTGTATAG	CTATACAC	ATTCTTAC
UDI 316	AACAATCA	TGATTGTT	TCGATTCA
UDI 317	CACTATCT	AGATAGTG	ATTCTCTA
UDI 318	GTGATCTC	GAGATCAC	ATTCGATG
UDI 319	CAATCCGT	ACGGATTG	TAAGCTAC

UDI 320	GTTTCGTCA	TGACGAAC	ATTCGTGT
UDI 321	CCACTTAA	TTAAGTGG	ATGAATAT
UDI 322	GTTGAATG	CATTCAAC	ATGAAGGA
UDI 323	CCAGCAAAG	CTTGCTGG	ATGAACTG
UDI 324	GTTAGTTC	GAACTAAC	ATGAGCAC
UDI 325	CCAGGCTA	TAGCCTGG	ATAGGAAT
UDI 326	GGTCCGCT	AGCGGACC	TAGACGGC
UDI 327	CACTTCCA	TGGAAGTG	GAATAGTG
UDI 328	GTTAAGTT	AACTTAAC	ATGACACC
UDI 329	AATCTAGT	ACTAGATT	ATGACGTT
UDI 330	GTTAATGA	TCATTAAC	ATGTATTA
UDI 331	CATACCTA	TAGGTATG	ATGTACCT
UDI 332	GTATGAGT	ACTCATAC	TCAATGGA
UDI 333	TAGATTAC	GTAATCTA	GAGGTCAC
UDI 334	GACCGAAT	ATTCGGTC	TAACCAGA
UDI 335	CGTACTT	AAGTAACG	ATGTGATT
UDI 336	GACCTGAC	GTCAGGTC	TAACAGCC
UDI 337	TTCGGCAT	ATGCCGAA	TATCTGTC
UDI 338	CCGAAGGC	GCCTTCGG	AGATAACT
UDI 339	GACCATGA	TCATGGTC	ATGTGGCA
UDI 340	GACGTGCA	TGCACGTC	TAACACTG
UDI 341	CGGCGACA	TGTCGCCG	ATGTCTGG
UDI 342	GACTAGGT	ACCTAGTC	TCAAGCAC
UDI 343	CGGTAGAC	GTCTACCG	ATGGTAAC
UDI 344	GACAGCTT	AAGCTGTC	ATGGTTCA
UDI 345	CGCTCGCA	TGCGAGCG	TCCAACGG
UDI 346	GAGATAGT	ACTATCTC	GAACAGAA
UDI 347	AACACTAC	GTAGTGTT	ATGGCCAG
UDI 348	GTGTGCGG	CCGCACAC	TCACGTGA
UDI 349	ACCGTG TG	CACACGGT	ATACAACC
UDI 350	TGGACACA	TGTGTCCA	TATGTTGG
UDI 351	AACTACTT	AAGTAGTT	ATGCTAGA
UDI 352	TTGAGGAA	TTCCTCAA	ATGCTGTC
UDI 353	AACTTAAT	ATTAAGTT	GATTGACC
UDI 354	TTGACTTA	TAAGTCAA	ATCATACT
UDI 355	AACTTGCC	GGCAAGTT	ATCAGCTA
UDI 356	GTGCACGT	ACGTGCAC	ATCACGCA
UDI 357	AACTTCGG	CCGAAGTT	GAGTTCTG

UDI 358	TGTACGAA	TTCGTACA	ATCACCGT
UDI 359	AACCTCAGC	GCTGAGTT	TCTTGAA
UDI 360	TGTATTCT	AGAATACA	ATCTAATC
UDI 361	GTCGTCGG	CCGACGAC	ATACATGA
UDI 362	CCGACGCA	TGCGTCGG	TATTGAAT
UDI 363	AACGAAGA	TCTTCGTT	GCTAGTCT
UDI 364	TGTCGGAT	ATCCGACA	ATCTAGCG
UDI 365	AACGACCG	CGGTCGTT	TCTCGCTA
UDI 366	TGTCGTGC	GCACGACA	ATCTGAAG
UDI 367	AACGTCAA	TTGACGTT	TCTAATGC
UDI 368	TGTAAGGT	ACCTTACA	GATAGCGC
UDI 369	ACTGAGAC	GTCTCAGT	CTCCTACA
UDI 370	TTGAATCG	CGATTCAA	TATCAAGC
UDI 371	AACGGTGT	ACACCGTT	ATCGCATT
UDI 372	AACCAAGTG	CACTGGTT	ATCCAGAA
UDI 373	AATATAAC	GTTATATT	ATACTATT
UDI 374	GGCGCGTG	CACGCGCC	GATGATAC
UDI 375	TTGTTACC	GGTAACAA	GCGGTATT
UDI 376	AACCTTGC	GCAAGGTT	ATCCGTTC
UDI 377	TGTGCAAT	ATTGCACA	AGAATTCA
UDI 378	AACCGAGG	CCTCGGTT	TATGACTT
UDI 379	TTGGATAC	GTATCCAA	GAGTCAGA
UDI 380	AACCGCAC	GTGCGGTT	AGAAGACG
UDI 381	TGTTAGTG	CACTAACA	AGAAGCTT
UDI 382	TATTTCGTA	TACGAATA	GAGTAGCA
UDI 383	ATAATTGT	ACAATTAT	AGAACCAA
UDI 384	TAGTCCAA	TTGGACTA	GCTTGGTG

Low Level Multiplexing Guidelines

Barcodes 1 and 2, 13 and 14, 25 and 26, 37 and 38, 49 and 50, 61 and 62, 73 and 74, and 85 and 86 are fully color balanced and are suitable to be used in a pool of two samples. When designing low-plexity index pools, always include two libraries barcoded with a set of two unique and fully color balanced barcodes to avoid laser color complexity issues during de-multiplexing. Additional libraries may be safely multiplexed with one set of fully color balanced barcodes in a pool.



The information provided in this document is for reference purposes only and may not be all-inclusive. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume liability for the accuracy or completeness of the information contained herein. Users should exercise caution when handling materials as they may present unknown hazards. Revvity shall not be liable for any damages or losses resulting from handling or contact with the product, as Revvity cannot control actual methods, volumes, or conditions of use. Users are responsible for ensuring the product's suitability for their specific application. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF.

www.revvity.com

revvity