

NEBNext® Multiplex Oligos for Illumina® (Unique Dual Index UMI Adaptors RNA Set 1)

NEB #E7416S/L

96/384 reactions

Version 4.0_2/24

Table of Contents

Workflow	2
Library Preparation Kits for use with NEBNext Multiplex Oligos for Illumina	3
NEBNext Adaptor for Illumina Overview	3
Section 1	
Setting up the PCR Reactions	4
Section 2	
Index Pooling Guidelines	5
Kit Components	11
Revision History	11

The NEBNext Multiplex Oligos for Illumina (Unique Dual Index UMI Adaptors RNA Set 1) Includes

*The volumes provided are sufficient for preparation of up to 96 reactions (NEB #E7416S) and 384 reactions (NEB #E7416L). All reagents should be stored at -20°C.**

- NEBNext Primer Mix
- NEBNext UMI Adaptor Dilution Buffer
- NEBNext UMI RNA Adaptor Plate
 - Each well contains a unique dual index UMI adaptor (S size contains 1 plate, L size contains 4 plates)

For the list of additional materials required, please check the manual for your NEBNext Library Prep Kit.

- * If the adaptor plate is thawed upon arrival, we recommend centrifuging the 96 well plate to collect the adaptor in the bottom of the well before re-freezing. If the plate arrived frozen, we recommend to store it at -20°C right away and centrifuge the plate prior to the first use to avoid unnecessary freeze/ thaw cycles.

Overview

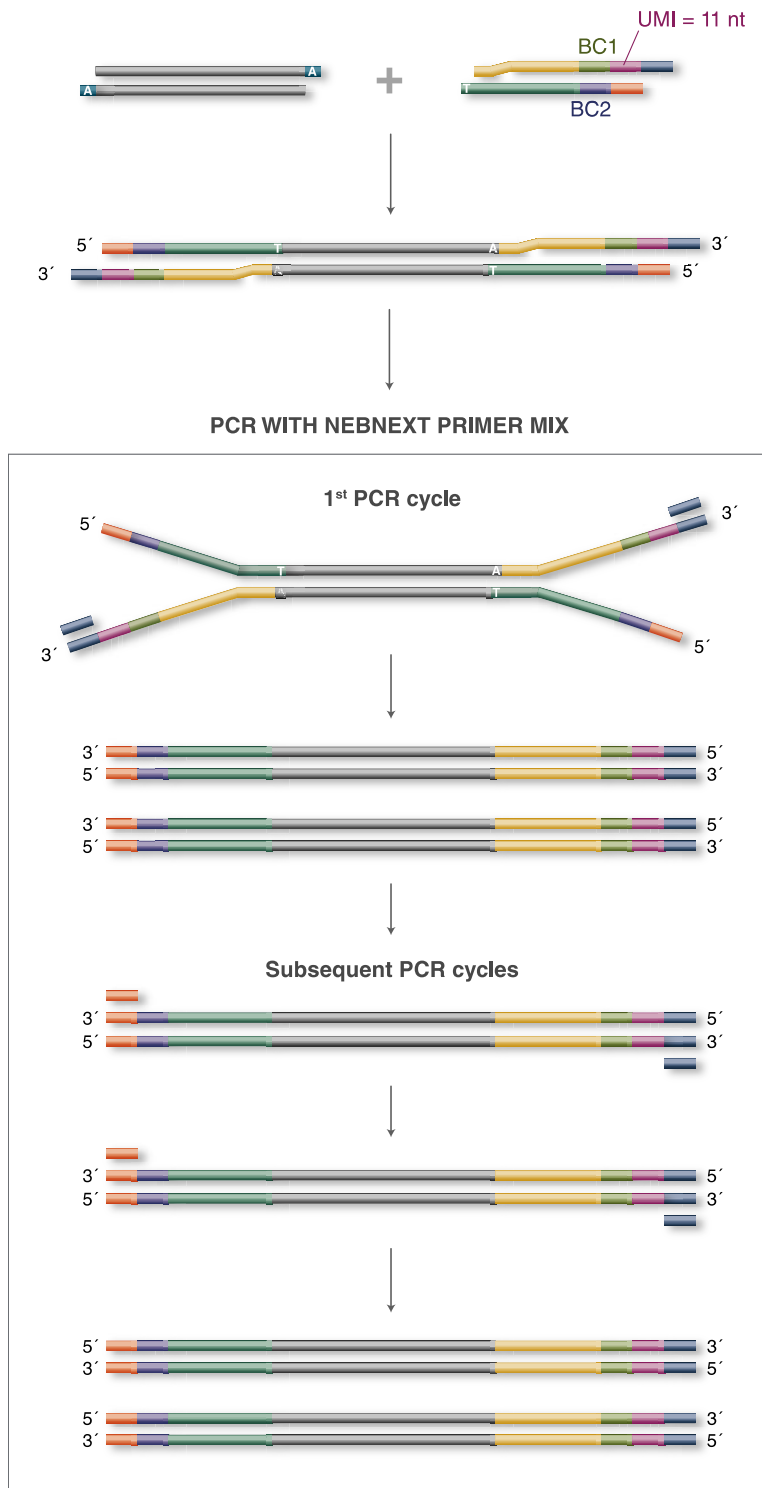
The NEBNext Multiplex Oligos for Illumina (Unique Dual Index UMI Adaptors RNA Set 1) contains adaptors and primers that are ideally suited for multiplex sample preparation for next-generation sequencing on the Illumina platform (Illumina, Inc.). Each kit component must pass rigorous quality control standards, and for each new lot the entire set of reagents is functionally validated together by construction and sequencing of indexed libraries on an Illumina sequencing platform.

Where larger volumes, customized or bulk packaging are required, we encourage consultation with the Customized Solutions team at NEB. Please complete the NEB Custom Contact Form at www.neb.com/CustomContactForm to learn more.

Workflow

Designed for use in library prep for cDNA and RNA (but not Small RNA), the NEBNext Unique Dual Index UMI Adaptors enable high-efficiency adaptor ligation and high library yields. These adaptors contain all necessary sequences for sequencing on the Illumina platform and sample pooling prior to PCR amplification. The incorporation of a 12-base unique molecular identifier (UMI) allows 1) accurate identification and removal of duplicate reads, and 2) consensus sequence building and error correction, ideally suited for accurate analysis of quantitative NGS data analysis. The 96 8-base unique dual index UMI adaptors included in this kit are packaged in a single-use 96-well plate with a pierceable foil seal. NEBNext Oligos can be used with NEBNext products, and with other standard Illumina-compatible library preparation protocols that are based on TA single base overhang ligation.

Figure 1. Workflow demonstrating the use of NEBNext Multiplex Oligos for Illumina (Unique Dual Index UMI Adaptors RNA Set 1).



Library Preparation Kits for use with NEBNext Unique Dual Index UMI Adaptors RNA

Please refer to the kit specific **library preparation kit manual** for using the NEBNext Multiplex Oligos for Illumina **for additional required materials that are not included.**

For compatibility of NEBNext Multiplex Oligos please refer to the NEBNext Multiplex Oligos Selection Chart at www.neb.com/oligos

NEBNext Adaptor for Illumina Overview

NEBNext Adaptor for Illumina sequence:

5'-/5Phos/GAT CGG AAG AGC ACA CGT CTG AAC TCC AGT CdUA CAC TCT TTC CCT ACA CGA CGC TCT TCC GAT C-s-T-3'

The following sequences are used for adaptor trimming of NEBNext adaptors for Illumina:

Read 1 AGATCGGAAGAGCACACGTCTGAACTCCAGTCA

Read 2 AGATCGGAAGAGCGTCGTGTAGGGAAAGAGTGT

Section 1

Setting up the Ligation Reactions

Symbols



This caution sign signifies a step in the protocol that has multiple paths leading to the same end point but is dependent on a user variable, like the number of samples to be processed.

1.1. Ligation



For < 96 samples, follow the protocol in Section 1.1A. For 96 samples, follow the protocol in Section 1.1B.

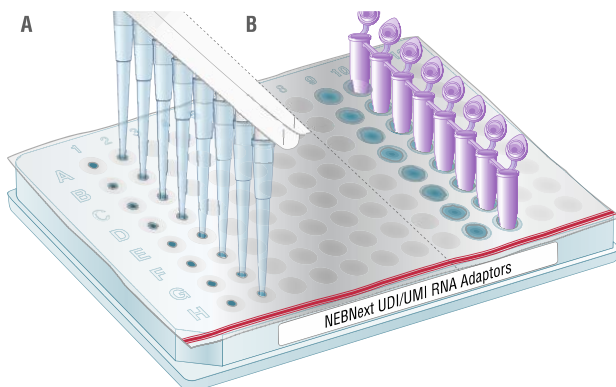
1.1A. Setting up the ligation reactions (< 96 samples)

- 1.1A.1. Determine the number of libraries that will be ligated and pooled for subsequent sequencing.
- 1.1A.2. Ensure that you choose a valid combination of barcode adaptors based on color balance guidelines in Section 2.
- 1.1A.3. Thaw the NEBNext UMI RNA Adaptor Plate for 10–15 minutes on ice.
- 1.1A.4. Remove the hard plastic plate cover. If necessary centrifuge the plate ($280 \times g$ for ~1 min) to collect all of the adaptor at the bottom of each well.
- 1.1A.5. Orient the NEBNext UMI RNA Adaptor Plate as indicated in Figure 1.1 (red stripe towards the user). With a pipette tip, pierce the desired well(s) (Figure 1.1A) and transfer the volume of adaptor mix required for the ligation reaction to the ligation plate/tubes (see specific library construction manual for protocol). It is important to change pipette tips before piercing a new well to avoid cross contamination of indexed adaptors. Alternatively, the wells can be pierced using the bottom of clean PCR strip tubes (see Figure 1.1B) prior to pipetting the adaptor mix. Use a new, clean strip tube for each new well to be pierced.

Note: Each well contains a unique pair of index adaptors. There is enough adaptor in each well for one library. Do not reuse adaptor if the seal has been previously pierced to avoid contamination with other indexed adaptors.

- 1.1A.6. Proceed with the ligation reaction according to the specific library construction manual.

Figure 1.1. NEBNext UMI RNA Adaptor Plate



1.1B. Setting up the ligation reactions (96 samples)

- 1.1B.1. Thaw the NEBNext UMI RNA Adaptor Plate for 10-15 minutes on ice.
- 1.1B.2. Remove the hard plastic plate cover. If necessary, centrifuge the plate ($280 \times g$ for ~1 min) to collect all of the adaptor at the bottom of each well.
- 1.1B.3. Orient the NEBNext UMI RNA Adaptor Plate as indicated in Figure 1.1 (red stripe towards the user). With a pipette tip, pierce the wells (Figure 1.1A) and transfer the volume of adaptor required for the ligation reaction to a 96 well plate (see specific library construction manual for protocol). It is important to change pipette tips before piercing a new well to avoid cross contamination of indexed adaptors. Alternatively, the wells can be pierced using the bottom of clean PCR strip tubes (see Figure 1.1B) prior to pipetting the adaptor mix. Use a new, clean strip tube for each new well to be pierced.

Note: Each well contains a unique pair of index adaptors. There is enough adaptor in each well for one ligation. Do not reuse adaptor if the seal has been previously pierced to avoid contamination with other indexed adaptors.

- 1.1B.4. Proceed with the ligation reaction according to the specific library construction manual.

Section 2

Index Pooling Guidelines: 96 Reaction Kit



For all HiSeq[®]/MiSeq[®] sequencers:

Illumina uses four channel chemistry with a red laser/LED to sequence bases A and C and a green laser/LED to sequence bases G and T. For each cycle, both the red and the green channel need to be read to ensure proper image registration (i.e. A or C must be in each cycle, and G or T must be in each cycle). If this color balance is not maintained, sequencing the index read could fail. The following tables list some valid combinations (up to 8-plex) for each Set that can be sequenced together. For combinations > 8 choose any column and add any plex combinations as needed.

For NovaSeq[®]6000/ NextSeq[®]/MiniSeq[®]:

Utilize red/ green or blue/ green 2 color chemistry. Valid index combinations must include some indices that do not start with GG in the first two cycles.

See Illumina document Document # 1000000041074 v12 [Chemistry and imaging on MiSeq - Illumina Knowledge](#)

For NovaSeq[®]X and X Plus:

Utilize blue/ green 2 color chemistry. Valid index combinations must include some indices that do not start with GG in the first two cycles. For additional NovaSeq X and X Plus color balancing guidelines please contact NEB technical support at info@neb.com.

Low Plex pooling options shown here are only for Illumina four channel chemistry.

Table 2.1. Index Pooling Guidelines

PLEX	WELL POSITION
< 4	Not recommended
4	A6, B6, C6, D6 A12, B12, C12, D12 B6, C6, D6, E6 B12, C12, D12, E12 C1, D1, E1, F1 C7, D7, E7, F7 E4, F4, G4, H4 E10, F10, G10, H10
5	A1, B1, C1, D1, E1 A6, B6, C6, D6, E6 A7, B7, C7, D7, E7 A12, B12, C12, D12, E12 B1, C1, D1, E1, F1 B6, C6, D6, E6, F6 B7, C7, D7, E7, F7 B12, C12, D12, E12, F12 C1, D1, E1, F1, G1 C2, D2, E2, F2, G2 C4, D4, E4, F4, G4 C7, D7, E7, F7, G7 C8, D8, E8, F8, G8 C10, D10, E10, F10, G10 D4, E4, F4, G4, H4 D10, E10, F10, G10, H10
6–7	Any 5 plex plus 1–2 adjacent wells from the same column
8	Any column

Four Channel Chemistry Color Balancing

***Forward Strand Workflow** for the following instruments: NovaSeq 6000 with v1.0 reagents kits, MiniSeq with rapid reagent kits, MiSeq®, HiSeq® 2000/2500 (pair-end flow cell), HiSeq 3000/4000 (single-read flow cell).

***Reverse Complement Workflow** for the following instruments: iSeq 100, MiniSeq with standard reagent kits, NextSeq Systems, NovaSeq 6000 with v1.5 reagent kits, HiSeq 2000/5000 (single-read flow cell), HiSeq 3000/4000 (paired-end flow cell).

See Illumina Document “Indexed Sequencing Overview” 15057455 and Illumina Guidelines for reverse complementing i5 sequences” for demultiplexing Illumina Knowledge Article #1800 [Guidelines for reverse complementing i5 sequences for demultiplexing - Illumina Knowledge](#).

Good and Bad Examples for Pooling and Color Balancing

Table 2.2. Listed below are index sequences color coded to correspond to the four color chemistry red/green channel. For combinations of valid indices, ensure that you will have signal in both the red and green channels in each cycle. See below for examples of Good and Bad index combinations based on four color chemistry guidelines:

BAD																								
WELL POSITION	EXPECTED i7 INDEX READ								EXPECTED i5 INDEX READ															
									FORWARD STRAND WORKFLOW*						REVERSE COMPLEMENT WORKFLOW*									
	E8	T	A	T	G	G	C	A	C	T	T	G	C	G	A	G	A	T	C	T	C	G	C	A
F8	G	A	A	T	C	A	C	C	G	A	A	C	G	A	A	G	C	T	T	C	G	T	T	C
G8	G	T	A	A	G	G	T	G	C	G	A	A	T	T	G	C	G	C	A	A	T	T	C	G
H8	C	G	A	G	A	G	A	A	G	G	A	A	G	A	G	A	T	C	T	C	T	T	C	C
	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	X	X	✓	✓	✓	✓	✓	✓	X	X	✓	✓	✓
A1	T	T	A	C	C	G	A	C	C	G	A	A	T	A	C	G	C	G	T	A	T	T	G	G
B1	T	C	G	T	C	T	G	A	G	T	C	C	T	T	G	A	T	C	A	A	G	G	A	C
C1	T	T	C	C	A	G	G	T	C	A	G	T	G	C	T	T	A	A	G	C	A	C	T	G
D1	T	A	C	G	G	T	C	T	T	C	C	A	T	T	G	C	G	C	A	A	T	G	G	A
	X	✓	✓	✓	✓	X	✓	✓	✓	✓	✓	✓	X	✓	✓	✓	✓	✓	✓	X	✓	✓	✓	✓
GOOD																								
WELL POSITION	EXPECTED i7 INDEX READ								EXPECTED i5 INDEX READ															
									FORWARD STRAND WORKFLOW						REVERSE COMPLEMENT WORKFLOW									
	C1	T	T	C	C	A	G	G	T	C	A	G	T	G	C	T	T	A	A	G	C	A	C	G
D1	T	A	C	G	G	T	C	T	T	C	C	A	T	T	G	C	G	C	A	A	T	G	G	A
E1	A	A	G	A	C	C	G	T	G	T	C	G	A	T	T	G	C	A	A	T	C	G	A	C
F1	C	A	G	G	T	T	C	A	A	T	A	A	C	G	C	C	G	G	C	G	T	T	A	T
	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓
A12	C	G	G	C	A	T	T	A	G	T	C	A	G	T	C	A	T	G	A	C	T	G	C	C
B12	C	A	C	G	C	A	A	T	C	C	T	T	C	C	A	T	A	T	G	G	A	A	G	G
C12	G	G	A	A	T	G	T	C	A	G	G	A	A	C	A	C	G	T	G	T	T	C	C	T
D12	T	G	G	T	G	A	A	G	C	T	T	A	C	A	G	C	G	C	T	G	T	A	A	G
	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓

The index adaptor sequences for different Illumina sequencer input sheets are indicated in Section 3.

Two Color Chemistry Color Balancing

NovaSeq 6000, NextSeq (500, 550, 1000 and 2000) and MiniSeq use red/ green or blue/ green 2 color chemistry to simplify nucleotide detection. See [Sequencing Chemistry \(illumina.com\)](http://www.illumina.com) Illumina Document # 1000000041074 v12 . For multiplexing a small number of samples, make sure the final index pool contains some indices that do not start with GG in the first two cycles. Listed here are some examples of good (signal in at least one channel for the first 2 cycles) and bad (the index read begins with GG) index combinations.

GOOD																								
WELL POSITION	EXPECTED i7 INDEX READ							EXPECTED i5 INDEX READ																
								FORWARD STRAND WORKFLOW					REVERSE COMPLEMENT WORKFLOW											
A12	C	G	G	C	A	T	T	A	G	T	C	A	G	T	C	A	T	G	A	C	T	G	C	C
B12	C	A	C	G	C	A	A	T	C	C	T	T	C	C	A	T	A	T	G	G	A	A	G	G
C12	G	G	A	A	T	G	T	C	A	G	G	A	A	C	A	C	G	T	G	T	T	C	C	T
D12	T	G	G	T	G	A	A	G	C	T	T	A	C	A	G	C	G	C	T	G	T	A	A	G
	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓

BAD																								
WELL POSITION	EXPECTED i7 INDEX READ							EXPECTED i5 INDEX READ																
								FORWARD STRAND WORKFLOW					REVERSE COMPLEMENT WORKFLOW											
C12	G	G	A	A	T	G	T	C	A	G	G	A	A	C	A	C	G	T	G	T	T	C	C	T
E12	G	G	A	C	A	T	C	A	T	A	C	C	T	G	C	A	T	G	C	A	G	G	T	A
F12	G	G	T	G	T	A	C	A	A	G	A	C	G	C	T	A	T	A	G	C	G	T	C	T
G11	G	G	T	T	G	A	A	C	T	C	C	A	C	G	T	T	A	A	C	G	T	G	G	A
	X	X	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓

Section 3 Index Sequences

Table 3.1 Index Sequences

WELL POSITION	EXPECTED i7 INDEX READ		EXPECTED i5 INDEX READ		
	i7 INDEX ID		i5 INDEX ID	FORWARD STRAND WORKFLOW	REVERSE COMPLEMENT WORKFLOW
A1	S762	TTACCGAC	S512	CGAATACG	CGTATTCG
B1	S713	TCGTCTGA	S586	GTCCTTGA	TCAAGGAC
C1	S736	TTCCAGGT	S543	CAGTGCTT	AAGCACTG
D1	S709	TACGGTCT	S575	TCCATTGC	GCAATGGA
E1	S732	AAGACCGT	S550	GTCGATTG	CAATCGAC
F1	S774	CAGGTTCA	S506	ATAACGCC	GGCGTTAT
G1	S747	TAGGAGCT	S524	GCCTTAAC	GTTAAGGC
H1	S794	TACTCCAG	S590	GGTATAGG	CCTATACC
A2	S729	AGTGACCT	S591	TCTAGGAG	CTCCTAGA
B2	S777	AGCCTATC	S526	TGCGTAAC	GTTACGCA
C2	S772	TCATCTCC	S567	CTTGCTAG	CTAGCAAG
D2	S725	CCAGTATC	S538	AGCGAGAT	ATCTCGCT
E2	S755	TTGCGAGA	S566	TATGGCAC	GTGCCATA
F2	S760	GAACGAAG	S511	GAATCACC	GGTGATTC
G2	S716	CGAATTGC	S559	GTAAGGTG	CACCTTAC
H2	S708	GGAAGAGA	S521	CGAGAGAA	TTCTCTCG
A3	S702	TCGGATTC	S523	CGCAACTA	TAGTTGCG
B3	S796	CTGTACCA	S507	CACAGACT	AGTCTGTG
C3	S757	GAGAGTAC	S545	TGGAAGCA	TGCTTCCA
D3	S783	TCTACGCA	S546	CAATAGCC	GGCTATTG
E3	S722	GCAATTCC	S578	CTCGAACA	TGTTCGAG
F3	S710	CTCAGAAG	S581	GGCAAGTT	AACTTGCC
G3	S770	GTCCTAAG	S540	AGCTACCA	TGGTAGCT
H3	S734	GCGTTAGA	S592	CAGCATAC	GTATGCTG
A4	S763	CAAGGTAC	S505	CGTATCTC	GAGATACG
B4	S797	AGACCTTG	S501	TTACGTGC	GCACGTAA
C4	S735	GTCGTTAC	S554	AGCTAAGC	GCTTAGCT
D4	S727	GTAACCGA	S598	AAGACACC	GGTGTCTT
E4	S742	GAATCCGT	S551	CAACTCCA	TGGAGTTG
F4	S795	CATGAGCA	S517	GATCTTGC	GCAAGATC
G4	S749	CTTAGGAC	S565	CTTCACTG	CAGTGAAG
H4	S773	ATCTGACC	S593	CTCGACTT	AAGTCGAG
A5	S769	TCCTCATG	S519	GTACACCT	AGGTGTAC
B5	S752	AGGATAGC	S544	CCAAGGTT	AACCTTGG
C5	S704	GGAGGAAT	S585	GAACGGTT	AACCGTTC
D5	S715	GACGTCAT	S518	CCAGTTGA	TCAACTGG
E5	S753	CCGCTTAA	S548	GTCATCGT	ACGATGAC
F5	S758	GACGAACT	S568	CAATGCGA	TCGCATTG
G5	S784	TCCACGTT	S541	GGTTGAAC	GTTCAACC
H5	S714	AACCAGAG	S520	CTTCGGTT	AACCGAAG

WELL POSITION	EXPECTED i7 INDEX READ		EXPECTED i5 INDEX READ		
	i7 INDEX ID		i5 INDEX ID	FORWARD STRAND WORKFLOW	REVERSE COMPLEMENT WORKFLOW
A6	S771	GTCAGTCA	S531	CGGCATTA	TAATGCCG
B6	S779	CCTTCCAT	S589	CACGCAAT	ATTGCGTG
C6	S788	AGGAACAC	S587	GGAATGTC	GACATTCC
D6	S739	CTTACAGC	S503	TGGTGAAG	CTTCACCA
E6	S737	TACCTGCA	S576	GGACATCA	TGATGTCC
F6	S728	AGACGCTA	S582	GGTGTACA	TGTACACC
G6	S780	CAACACAG	S530	GATAGCCA	TGGCTATC
H6	S761	GTACCACA	S533	CCACAACA	TGTTGTGG
A7	S712	CGAATACG	S562	TTACCGAC	GTCGGTAA
B7	S786	GTCCTTGA	S513	TCGTCTGA	TCAGACGA
C7	S743	CAGTGCTT	S536	TTCCAGGT	ACCTGGAA
D7	S775	TCCATTGC	S509	TACGGTCT	AGACCGTA
E7	S750	GTCGATTG	S532	AAGACCGT	ACGGTCTT
F7	S706	ATAACGCC	S574	CAGGTTCA	TGAACCTG
G7	S724	GCCTTAAC	S547	TAGGAGCT	AGCTCCTA
H7	S790	GGTATAGG	S594	TACTCCAG	CTGGAGTA
A8	S791	TCTAGGAG	S529	AGTGACCT	AGGTCACT
B8	S726	TGCGTAAC	S577	AGCCTATC	GATAGGCT
C8	S767	CTTGCTAG	S572	TCATCTCC	GGAGATGA
D8	S738	AGCGAGAT	S525	CCAGTATC	GATACTGG
E8	S766	TATGGCAC	S555	TTGCGAGA	TCTCGCAA
F8	S711	GAATCACC	S560	GAACGAAG	CTTCGTTC
G8	S759	GTAAGGTG	S516	CGAATTGC	GCAATTCC
H8	S721	CGAGAGAA	S508	GGAAGAGA	TCTCTTCC
A9	S723	CGCAACTA	S502	TCGGATTG	GAATCCGA
B9	S707	CACAGACT	S596	CTGTACCA	TGGTACAG
C9	S745	TGGAAGCA	S557	GAGAGTAC	GTA CTCTC
D9	S746	CAATAGCC	S583	TCTACGCA	TGCGTAGA
E9	S778	CTCGAACA	S522	GCAATTCC	GGAATTGC
F9	S781	GGCAAGTT	S510	CTCAGAA G	CTTCTGAG
G9	S740	AGCTACCA	S570	GTCCTAAG	CTTAGGAC
H9	S792	CAGCATA C	S534	GCGTTAGA	TCTAACGC
A10	S705	CGTATCTC	S563	CAAGGTAC	GTACCTTG
B10	S701	TTACGTGC	S597	AGACCTTG	CAAGGTCT
C10	S754	AGCTAAGC	S535	GTCGTTAC	GTAACGAC
D10	S798	AAGACACC	S527	GTAACCGA	TCGGTTAC
E10	S751	CAACTCCA	S542	GAATCCGT	ACGGATT C
F10	S717	GATCTTGC	S595	CATGAGCA	TGCTCATG
G10	S765	CTTCACTG	S549	CTTAGGAC	GTCCTAAG
H10	S793	CTCGACTT	S573	ATCTGACC	GGTCAGAT

WELL POSITION	EXPECTED i7 INDEX READ		EXPECTED i5 INDEX READ		
	i7 INDEX ID		i5 INDEX ID	FORWARD STRAND WORKFLOW	REVERSE COMPLEMENT WORKFLOW
A11	S 719	GTACACCT	S 569	TCCTCATG	CATGAGGA
B11	S 744	CCAAGGTT	S 552	AGGATAGC	GCTATCCT
C11	S 785	GAACGGTT	S 504	GGAGGAAT	ATTCCTCC
D11	S 718	CCAGTTGA	S 515	GACGTCAT	ATGACGTC
E11	S 748	GTCATCGT	S 553	CCGCTTAA	TTAAGCGG
F11	S 768	CAATGCGA	S 558	GACGAACT	AGTTCGTC
G11	S 741	GGTTGAAC	S 584	TCCACGTT	AACGTGGA
H11	S 720	CTTCGGTT	S 514	AACCAGAG	CTCTGGTT
A12	S 731	CGGCATTA	S 571	GTCAGTCA	TGACTGAC
B12	S 789	CACGCAAT	S 579	CCTTCCAT	ATGGAAGG
C12	S 787	GG AATGTC	S 588	AGGAACAC	GTGTTCTT
D12	S 703	TGGTGAAG	S 539	CTTACAGC	GCTGTAAG
E12	S 776	GGACATCA	S 537	TACCTGCA	TGCAGGTA
F12	S 782	GGTGTACA	S 528	AGACGCTA	TAGCGTCT
G12	S 730	GATAGCCA	S 580	CAACACAG	CTGTGTTG
H12	S 733	CCACAACA	S 561	GTACCACA	TGTGGTAC

Sequencing on the Illumina Platform

Pool equal molar amounts of libraries for sequencing on the Illumina platforms using the cycles settings shown in the table below.

RUN SEGMENT	CYCLE NUMBER
Read 1	X defined by users
Index 1 (i7)	8 (without UMI)
	20 (with UMI)
Index 2 (i5)	8
Read 2	X defined by users

Index Sequence File

For a link to download a sample sheet with the index sequences for use with the Illumina Experiment Manager (IEM) please go to our FAQ's tab on www.neb.com/E7416 or you can access the sample sheets by visiting the "[Usage Guidelines](#)" sub tab located under the "protocols, manuals and usage" tab on the E7416 product page.

Kit Components

The NEBNext Multiplex Oligos for Illumina (Unique Dual Index UMI Adaptors RNA Set 1) are functionally validated through library preparation using the NEBNext Library Prep Kits and sequencing on the Illumina platforms.

NEB #E7416S Table of Components

NEB #	CONCENTRATION	PRODUCT	VOLUME
E7417A	1 μ M	NEBNext UMI RNA Adaptor Plate	1 plate (5 μ l/well)
E7397A	40 μ M (Total)	NEBNext Primer Mix	0.48 ml
E7398A		NEBNext UMI Adaptor Dilution Buffer	5 ml

NEB #E7416L Table of Components

NEB #	CONCENTRATION	PRODUCT	VOLUME
E7417A	1 μ M	NEBNext UMI RNA Adaptor Plate	4 plates (5 μ l/well)
E7397AA	40 μ M (Total)	NEBNext Primer Mix	2 x 0.96 ml
E7398AA		NEBNext UMI Adaptor Dilution Buffer	20 ml

Revision History

REVISION #	DESCRIPTION	DATE
1.0	N/A	3/20
2.0	Updated tables to have the most current Illumina instrument information and removed HiSeqX.	3/21
3.0	Updated protocol.	8/22
4.0	Updated primer sequences, indexing pool guidelines, header/footer and legal footer.	2/24

This product is intended for research purposes only. This product is not intended to be used for therapeutic or diagnostic purposes in humans or animals.

Products and content are covered by one or more patents, trademarks and/or copyrights owned or controlled by New England Biolabs, Inc (NEB). The use of trademark symbols does not necessarily indicate that the name is trademarked in the country where it is being read; it indicates where the content was originally developed. See www.neb.com/trademarks. The use of these products may require you to obtain additional third-party intellectual property rights for certain applications. For more information, please email busdev@neb.com.

B CORPORATION[®] is a registered trademark of B Lab IP, LLC, Inc.

ILLUMINA[®], HISEQ[®], MISEQ[®], NEXTSEQ[®], NOVASEQ[®] and MINISEQ[®] are registered trademarks of Illumina, Inc.

© Copyright 2024, New England Biolabs, Inc.; all rights reserved

