

NEBNext® Multiplex Oligos for Illumina® (96 Unique Dual Index Primer Pairs Set 3)

NEB #E6444S/L

96/384 reactions

Version 4.0_7/22

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	9
Revision History	10

The NEBNext Multiplex Oligos for Illumina (96 Unique Dual Index Primer Pairs Set 3) Includes

The volumes provided are sufficient for preparation of up to 96 reactions (NEB #E6444S) and 384 reactions (NEB #E6444L). All reagents should be stored at –20°C.

NEBNext Adaptor for Illumina

USER® Enzyme

NEBNext 96 Unique Dual Index Primer Pairs Plate (Set 3)

Each well contains a unique pair of Index Primers (S size contains 1 plate, L size contains 4 plates)

Overview

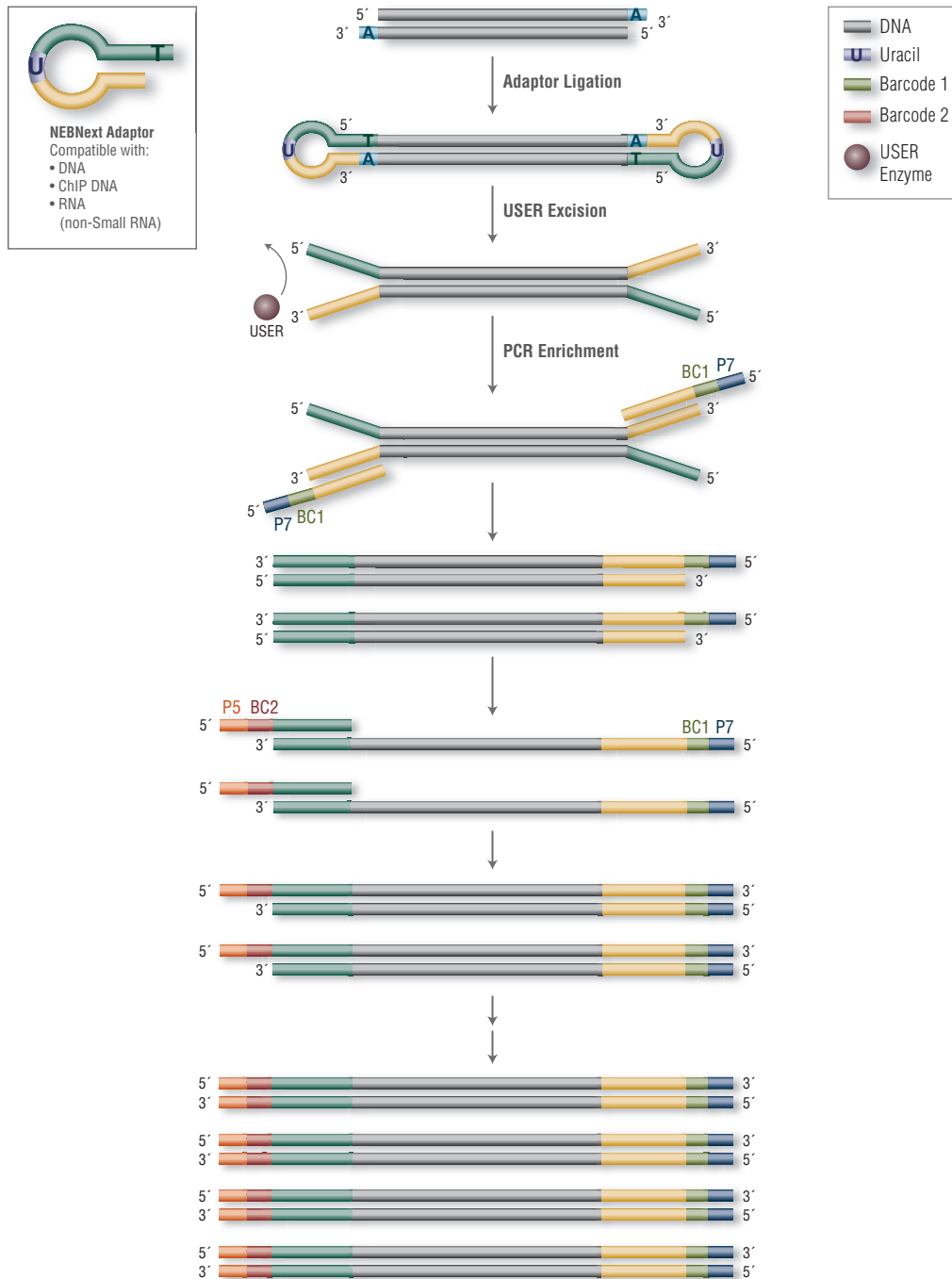
The NEBNext Multiplex Oligos for Illumina (96 Unique Dual Index Primer Pairs Set 3) contains adaptors and primers that are ideally suited for multiplex sample preparation for next-generation sequencing on the Illumina platform (Illumina, Inc.). Each kit component must pass rigorous quality control standards, 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.

For larger volume requirements, customized and bulk packaging is available by purchasing through the OEM/Bulks department at NEB. Please contact OEM@neb.com for further information.

Workflow

Designed for use in library prep for DNA, ChIP DNA and RNA (but not Small RNA), the NEBNext Adaptors enable high-efficiency adaptor ligation and high library yields, with minimized adaptor-dimer formation. Incorporating a novel hairpin loop structure, the NEBNext Adaptor ligates with increased efficiency to end-repaired, dA-tailed DNA. The loop contains a U, which is removed by treatment with USER Enzyme (a combination of UDG and Endo VIII), to open up the loop and make it available as a substrate for PCR. During PCR, barcodes can be incorporated by use of the NEBNext index primers, thereby enabling multiplexing. The 96 8-base index primer pairs included in this kit are pre-mixed and 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.

Figure 1. Workflow demonstrating the use of NEBNext Multiplex Oligos for Illumina (96 Unique Dual Index Primer Pairs).



Please Refer to the Kit Specific Protocol for using the NEBNext Multiplex Oligos for Illumina

For compatibility of NEBNext Multiplex Oligos please refer to the NEBNext Multiplex Oligos Selection Chart at 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 PCR 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 amount of input DNA.

1.1. PCR Amplification

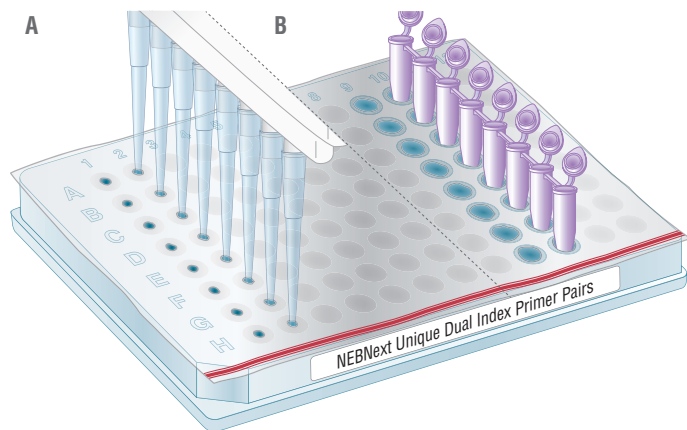


For < 96 samples, follow the protocol in Section 1.1A. For 96 samples, follow the protocol in Section 1.1B.

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

- 1.1A.1. Determine the number of libraries that will be amplified and pooled for subsequent sequencing.
- 1.1A.2. Ensure that you choose a valid combination of barcode primers based on color balance guidelines in Section 2.
- 1.1A.3. Thaw the 96 Unique Dual Index Primers Plate for 10-15 minutes at room temperature.
- 1.1A.4. Remove the hard plastic plate cover. Mix briefly by vortexing and then centrifuge the plate ($280 \times g$ for ~1 min) to collect all of the primer at the bottom of each well.
- 1.1A.5. Orient the 96 Unique Dual Index Primers Plate Set 3 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 primer mix required for the PCR reaction to the PCR 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 primers. Alternatively, the wells can be pierced using the bottom of clean PCR strip tubes (see Figure 1.1B) prior to pipetting the primer mix. Use a new, clean strip tube for each new well to be pierced.
Note: Each well contains a unique pair of index primers. There is enough primer in each well for one PCR reaction. Do not reuse primer if the seal has been previously pierced to avoid contamination with other indexed primers.
- 1.1A.6. Proceed with the PCR reaction according to the specific library construction manual.

Figure 1.1. NEBNext Unique Dual Index Pairs Plate Set 3



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

- 1.1B.1. Thaw the 96 Unique Dual Index Primer Pairs plate for 10-15 minutes at room temperature.
- 1.1B.2. Remove the hard plastic plate cover. Mix briefly by vortexing and then centrifuge the plate ($280 \times g$ for ~1 min) to collect all of the primer at the bottom of each well.
- 1.1B.3. Orient the 96 Unique Dual Index Primer Pairs 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 primer mix required for the PCR reaction to the PCR 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 primers. Alternatively, the wells can be pierced using the bottom of clean PCR strip tubes (see Figure 1.1B) prior to pipetting the primer mix. Use a new, clean strip tube for each new well to be pierced.
Note: Each well contains a unique pair of index primers. There is enough primer in each well for one PCR reaction. Do not reuse primer if the seal has been previously pierced to avoid contamination with other indexed primers.
- 1.1B.4. Proceed with the PCR reaction according to the specific library construction manual.

Section 2

Index Pooling Guidelines: 96 Reaction Kit



For a link to download a sample sheet with the index sequences for use with the Illumina Experiment Manager (IEM) please visit the "[Usage Guidelines](#)" sub tab located under the "protocols, manuals and usage" tab on the E6444 product page.

For all HiSeq[®]/MiSeq[®] sequencers, Illumina uses 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. See Table 2.1 for examples of Good and Bad Index combinations.

For the NovaSeq[®]/NextSeq[®]/MiniSeq[®] which utilize 2 color chemistry, valid index combinations must include some indices that do not start with GG in the first two cycles. See Table 2.2 for examples of Good and Bad Index combinations.

The barcoded primers are organized on the plate such that including the primers in rows A and B from any column will produce a color balanced pool. For example, if preparing 2 libraries, choose primer wells A1 and B1. For pools containing 3-8 libraries, add any other primers from that column. For pools containing more than 8 libraries, choose any column and add any other primers as needed.

***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 Strand 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).

Table 2.1. lists each index sequence color coded to correspond to the 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 HiSeq/MiSeq guidelines:

GOOD																								
WELL POSITION	EXPECTED i7 INDEX READ								EXPECTED i5 INDEX READ															
									FORWARD STRAND WORKFLOW*						REVERSE STRAND WORKFLOW*									
A1	T	G	T	C	G	T	A	G	A	A	A	G	C	T	A	A	T	T	A	G	C	T	T	T
B1	C	A	A	T	C	A	T	A	T	G	G	A	G	A	T	T	A	A	T	C	T	C	C	A
C1	G	T	T	C	T	T	A	T	A	A	T	T	A	G	A	C	G	T	C	T	A	A	T	T
D1	G	A	T	G	C	G	A	C	A	C	T	T	T	G	G	G	C	C	C	A	A	A	G	T
	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓

BAD																								
WELL POSITION	EXPECTED i7 INDEX READ								EXPECTED i5 INDEX READ															
									FORWARD STRAND WORKFLOW						REVERSE STRAND WORKFLOW									
E1	G	A	A	G	A	G	G	G	C	G	G	A	C	G	G	A	T	C	C	G	T	C	C	G
F1	T	A	G	T	A	A	T	C	G	C	A	G	A	G	C	C	G	G	C	T	C	T	G	C
G1	G	T	G	T	G	G	A	G	G	C	A	T	G	A	T	C	G	A	T	C	A	T	G	C
H1	A	C	G	T	T	G	T	A	T	C	G	A	C	C	T	A	T	A	G	G	T	C	G	A
	✓	✓	✓	X	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	X	X	✓	✓	✓	✓	✓	✓	✓

The index primer sequences, for different Illumina sequencer input sheets are indicated in Table 2.3.

Table 2.2. NovaSeq, NextSeq and MiniSeq use 2 color channel sequencing to simplify nucleotide detection. Clusters only in red or green are interpreted as C or T, respectively. Clusters in both red and green are read as A, while unlabeled clusters are G bases. 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 STRAND WORKFLOW										
A1	T	G	T	C	G	T	A	G	A	A	A	G	C	T	A	A	T	T	A	G	C	T	T	T
B1	C	A	A	T	C	A	T	A	T	G	G	A	G	A	T	T	A	A	T	C	T	C	C	A
C1	G	T	T	C	T	T	A	T	A	A	T	T	A	G	A	C	G	T	C	T	A	A	T	T
	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓

BAD																								
WELL POSITION	EXPECTED i7 INDEX READ								EXPECTED i5 INDEX READ															
									FORWARD STRAND WORKFLOW					REVERSE STRAND WORKFLOW										
D9	C	A	G	C	A	G	G	G	A	T	G	C	T	C	C	C	G	G	G	A	G	C	A	T
G9	C	G	A	T	A	C	A	T	C	G	A	G	A	A	C	C	G	G	T	T	C	T	C	G
D11	C	T	C	A	C	G	T	C	T	A	T	T	T	A	C	C	G	G	T	A	A	A	T	A
	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	X	X	✓	✓	✓	✓	✓	✓

Table 2.3. Lists each index sequence color coded to correspond to the red/green channel. For combinations of valid indices, ensure that you will have signal in both the red and green channels in each cycle.

WELL POSITION	EXPECTED i7 INDEX READ		EXPECTED i5 INDEX READ		
	i7 INDEX ID		i5 INDEX ID	FORWARD STRAND WORKFLOW	REVERSE STRAND WORKFLOW
A1	7-197	TGTCGTAG	5-245	AAAGCTAA	TTAGCTTT
B1	7-198	CAATCATA	5-246	TGGAGATT	AATCTCCA
C1	7-199	GTTCTTAT	5-247	AATTAGAC	GTCTAATT
D1	7-200	GATGCGAC	5-248	ACTTTGGG	CCCAAAGT
E1	7-201	GAAGAGGG	5-249	CGGACGGA	TCCGTCCG
F1	7-202	TAGTAATC	5-250	GCAGAGCC	GGCTCTGC
G1	7-203	GTGTGGAG	5-251	GCATGATC	GATCATGC
H1	7-204	ACGTTGTA	5-252	TCGACCTA	TAGGTCGA
A2	7-205	GCGCTAAT	5-253	ACCCTGAC	GTCAGGGT
B2	7-206	AGAGCTGC	5-254	GTTGAAGG	CCTTCAAC
C2	7-207	CATACTTA	5-255	GCACGGGA	TCCCGTGC
D2	7-208	TTGCACCG	5-256	CGTATAAA	TTTATACG
E2	7-209	GCGGGATA	5-257	AGAGACGG	CCGTCTCT
F2	7-210	GAAGTGAA	5-258	TACAAGTC	GACTTGTA
G2	7-211	CTGTTTAC	5-259	TGAATCTT	AAGATTCA
H2	7-212	GAGCACTC	5-260	GCAACTTG	CAAGTTGC
A3	7-213	TTGTTGCA	5-261	ACGACGTC	GACGTCGT
B3	7-214	CCACACTT	5-262	GTATGACG	CGTCATAC
C3	7-215	CCCGTTTG	5-263	TACAGCAA	TTGCTGTA
D3	7-216	ATGCTCCC	5-264	CAGCAGGG	CCCTGCTG
E3	7-217	GCTCAATA	5-265	GATAAATG	CATTTATC
F3	7-218	GTAGTTCG	5-266	GCATCAAG	CTTGATGC
G3	7-219	CGAGAACC	5-267	CGATACAT	ATGTATCG
H3	7-220	GCCATGTA	5-268	AACCCTAT	ATAGGGTT
A4	7-221	TTTCTCTA	5-269	ACGTCGAG	CTCGACGT
B4	7-222	CCAGCGAT	5-270	TGACTAGA	TCTAGTCA
C4	7-223	TGGGAGTG	5-271	TAGACGGG	CCCGTCTA
D4	7-224	CCCTCGTA	5-272	CTCTTCTA	TAGAAGAG
E4	7-225	CGATATGG	5-273	TACGTCCC	GGGACGTA
F4	7-226	TTGTGCCC	5-274	GATGGAAA	TTTCCATC
G4	7-227	TGTCCTCT	5-275	GTTGTCGC	CGACGAAC
H4	7-228	GTATAGTC	5-276	GAGACCAA	TTGGTCTC
A5	7-229	TTTGGGAT	5-277	ACGTGAAC	GTTCACGT
B5	7-230	CACCAAGC	5-278	TTCCCTTT	AAAGGGAA
C5	7-231	CGGAGAGG	5-279	GACGCTCG	CGAGCGTC
D5	7-232	TATTTACC	5-280	CTCACGTC	GACGTGAG
E5	7-233	TATATGGA	5-281	CTGCCAAG	CTTGGCAG
F5	7-234	GTTAACAT	5-282	ACGCCGCA	TGCGGCGT
G5	7-235	CGTCTTGG	5-283	CGCCAGTC	GACTGGCG
H5	7-236	CGTAGCGA	5-284	CTAAACAA	TTGTTTAG

WELL POSITION	EXPECTED i7 INDEX READ		EXPECTED i5 INDEX READ		
	i7 INDEX ID		i5 INDEX ID	FORWARD STRAND WORKFLOW	REVERSE STRAND WORKFLOW
A6	7-237	TAGTCACA	5-285	TATACCTC	GAGGTATA
B6	7-238	AGAA GTGG	5-286	CTCTTGAT	ATCAAGAG
C6	7-239	CGTGGATT	5-287	ACTCTTAG	CTAAGAGT
D6	7-240	GTAGATGC	5-288	GAGCAACA	TGTTGCTC
E6	7-241	TACCGCTC	5-289	CAGTGACG	CGTCACTG
F6	7-242	CGAACCAC	5-290	AAGATTGA	TCAATCTT
G6	7-243	TATTGTTT	5-291	GTGTGTTT	AAACACAC
H6	7-244	GTTGTGTG	5-292	CGTCCGAC	GTCGGACG
A7	7-245	AAAGCTAA	5-197	TGTCGTAG	CTACGACA
B7	7-246	TGGAGATT	5-198	CAATCATA	TATGATTG
C7	7-247	AATTAGAC	5-199	GTTCTTAT	ATAAGAAC
D7	7-248	ACTTTGGG	5-200	GATGCGAC	GTCGCATC
E7	7-249	CGGACGGA	5-201	GAAGAGGG	CCCTCTTC
F7	7-250	GCAGAGCC	5-202	TAGTAATC	GATTACTA
G7	7-251	GCATGATC	5-203	GTGTGGAG	CTCCACAC
H7	7-252	TCGACCTA	5-204	ACGTTGTA	TACAACGT
A8	7-253	ACCCTGAC	5-205	GCGCTAAT	ATTAGCGC
B8	7-254	GTTGAAGG	5-206	AGAGCTGC	GCAGCTCT
C8	7-255	GCACGGGA	5-207	CATACTTA	TAAGTATG
D8	7-256	CGTATAAA	5-208	TTGCACCG	CGGTGCAA
E8	7-257	AGAGACGG	5-209	GCGGGATA	TATCCCGC
F8	7-258	TACAAGTC	5-210	GAAGTGAA	TTCACTTC
G8	7-259	TGAATCTT	5-211	CTGTTTAC	GTA AACAG
H8	7-260	GCAACTTG	5-212	GAGCACTC	GAGTGCTC
A9	7-261	ACGACGTC	5-213	TTGTTGCA	TGCAACAA
B9	7-262	GTATGACG	5-214	CCACACTT	AAGTGTGG
C9	7-263	TACAGCAA	5-215	CCCGTTTG	CAAACGGG
D9	7-264	CAGCAGGG	5-216	ATGCTCCC	GGGAGCAT
E9	7-265	GATAAATG	5-217	GCTCAATA	TATTGAGC
F9	7-266	GCATCAAG	5-218	GTAGTTCG	CGAACTAC
G9	7-267	CGATACAT	5-219	CGAGAACC	GGTTCTCG
H9	7-268	AACCCTAT	5-220	GCCATGTA	TACATGGC
A10	7-269	ACGTCGAG	5-221	TTTCTCTA	TAGAGAAA
B10	7-270	TGACTAGA	5-222	CCAGCGAT	ATCGCTGG
C10	7-271	TAGACGGG	5-223	TGGGAGTG	CACTCCCA
D10	7-272	CTCTTCTA	5-224	CCCTCGTA	TACGAGGG
E10	7-273	TACGTCCC	5-225	CGATATGG	CCATATCG
F10	7-274	GATGGAAA	5-226	TTGTGCCC	GGGCACAA
G10	7-275	GTTTCGTCG	5-227	TGTCCTCT	AGAGGACA
H10	7-276	GAGACCAA	5-228	GTATAGTC	GACTATAC

WELL POSITION	EXPECTED i7 INDEX READ		EXPECTED i5 INDEX READ		
	i7 INDEX ID		i5 INDEX ID	FORWARD STRAND WORKFLOW	REVERSE STRAND WORKFLOW
A11	7-277	ACGTGAAC	5-229	TTTGGGAT	ATCCCAA
B11	7-278	TTCCCTTT	5-230	CACCAAGC	GCTTGGTG
C11	7-279	GACGCTCG	5-231	CGGAGAGG	CCTCTCCG
D11	7-280	CTCACGTC	5-232	TATTTACC	GGTAAATA
E11	7-281	CTGCCAAG	5-233	TATATGGA	TCCATATA
F11	7-282	ACGCCGCA	5-234	GTTAACAT	ATGTTAAC
G11	7-283	CGCCAGTC	5-235	CGTCTTGG	CCAAGACG
H11	7-284	CTAAACAA	5-236	CGTAGCGA	TCGCTACG
A12	7-285	TATACCTC	5-237	TAGTCACA	TGTGACTA
B12	7-286	CTCTTGAT	5-238	AGAAGTGG	CCACTTCT
C12	7-287	ACTCTTAG	5-239	CGTGGATT	AATCCACG
D12	7-288	GAGCAACA	5-240	GTAGATGC	GCATCTAC
E12	7-289	CAGTGACG	5-241	TACCGCTC	GAGCGGTA
F12	7-290	AAGATTGA	5-242	CGAACCCAC	GTGGTTCG
G12	7-291	GTGTGTTT	5-243	TATTGTTC	GAACAATA
H12	7-292	CGTCCGAC	5-244	GTTGTGTG	CACACAAC

Kit Components

The NEBNext Multiplex Oligos for Illumina (96 Unique Dual Index Primer Pairs Set 3) are functionally validated through library preparation using the NEBNext Library Prep Kits and sequencing on the Illumina platforms.

NEB #E6444S Table of Components

NEB #	CONCENTRATION	PRODUCT	VOLUME
E6612A	15 µM	NEBNext Adaptor for Illumina	0.96 ml
E6610A		USER Enzyme	0.288 ml
E6445A	5 µM each	NEBNext 96 Unique Dual Index Primer Pairs Plate (Set 3)	1 plate (10 µl/well)

NEB #E6444L Table of Components

NEB #	CONCENTRATION	PRODUCT	VOLUME
E6612A	15 µM	NEBNext Adaptor for Illumina	4 x 0.96 ml
E6610AA		USER Enzyme	2 x 0.576 ml
E6445A	5 µM each	NEBNext 96 Unique Dual Index Primer Pairs Plate (Set 3)	4 plates (10 µl/well)

Note :

For the NEBNext Adaptor for Illumina sequence, please see NEBNext Multiplex Oligos for Illumina (Index Primers Set 1), NEB #E7335, Manual.

Revision History

REVISION #	DESCRIPTION	DATE
1.0	N/A	5/20
2.0	Update concentration of E6445A in Table of Components	7/20
3.0	Updating tables to have the most current Illumina instrument information and removed HiSeqX.	2/21
4.0	Update Protocol and Tables	7/22

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.

This product is covered by one or more patents, trademarks and/or copyrights owned or controlled by New England Biolabs, Inc. The use of trademark symbols does not necessarily indicate that the name is trademarked in the country where it is being read; rather, it indicates where the document was originally developed. For more information about commercial rights, please email us at busdev@neb.com. While NEB develops and validates its products for various applications, the use of this product may require you to obtain additional third party intellectual property rights for certain applications.

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

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

© Copyright 2022, New England Biolabs, Inc.; all rights reserved



be INSPIRED
drive DISCOVERY
stay GENUINE