

## NEBNext® Multiplex Oligos for Illumina® (Dual Index Primers Set 2)

NEB #E7780S

96 reactions

Version 2.0\_8/22

### Table of Contents

Applications.....	2
Section 1	
Setting up the PCR Reaction .....	3
Section 2	
Appendix A: Principle for Use and Pooling Guide .....	5
Section 3	
Appendix B: PCR Setup Template .....	8
NEBNext i509 Primer–i516 Primer .....	9
NEBNext i713 Primer–i724 Primer .....	9
Kit Components.....	10
Revision History .....	11

### The NEBNext Multiplex Oligos for Illumina (Dual Index Primers Set 2) Includes

The volumes provided are sufficient for preparation of up to 96 reactions (NEB #E7780S). Primers are supplied at 10  $\mu$ M. All reagents should be stored at  $-20^{\circ}\text{C}$ . Colored bullets represent the color of the cap of the tube containing the reagent.

- (red) NEBNext Adaptor for Illumina
- (red) USER® Enzyme
- (white) NEBNext i509 Primer
- (white) NEBNext i510 Primer
- (white) NEBNext i511 Primer
- (white) NEBNext i512 Primer
- (white) NEBNext i513 Primer
- (white) NEBNext i514 Primer
- (white) NEBNext i515 Primer
- (white) NEBNext i516 Primer
- (green) NEBNext i713 Primer
- (green) NEBNext i714 Primer
- (green) NEBNext i715 Primer
- (green) NEBNext i716 Primer
- (green) NEBNext i717 Primer
- (green) NEBNext i718 Primer
- (green) NEBNext i719 Primer
- (green) NEBNext i720 Primer
- (green) NEBNext i721 Primer
- (green) NEBNext i722 Primer
- (green) NEBNext i723 Primer
- (green) NEBNext i724 Primer

## Applications

The NEBNext Multiplex Oligos for Illumina (Dual Index Primers Set 2) contains the adaptor and index primers that are ideally suited for multiplex sample preparation for next-generation sequencing on the Illumina platform. Each of these components must pass rigorous quality control standards and are lot controlled, both individually and as a set of reagents.

**Lot Control:** The lots provided in the NEBNext Multiplex Oligos for Illumina (Dual Index Primers Set 2) are managed separately and are qualified by additional functional validation. Individual reagents undergo standard enzyme activity and quality control assays, and also meet stringent criteria in the additional quality controls specific for each individual component.

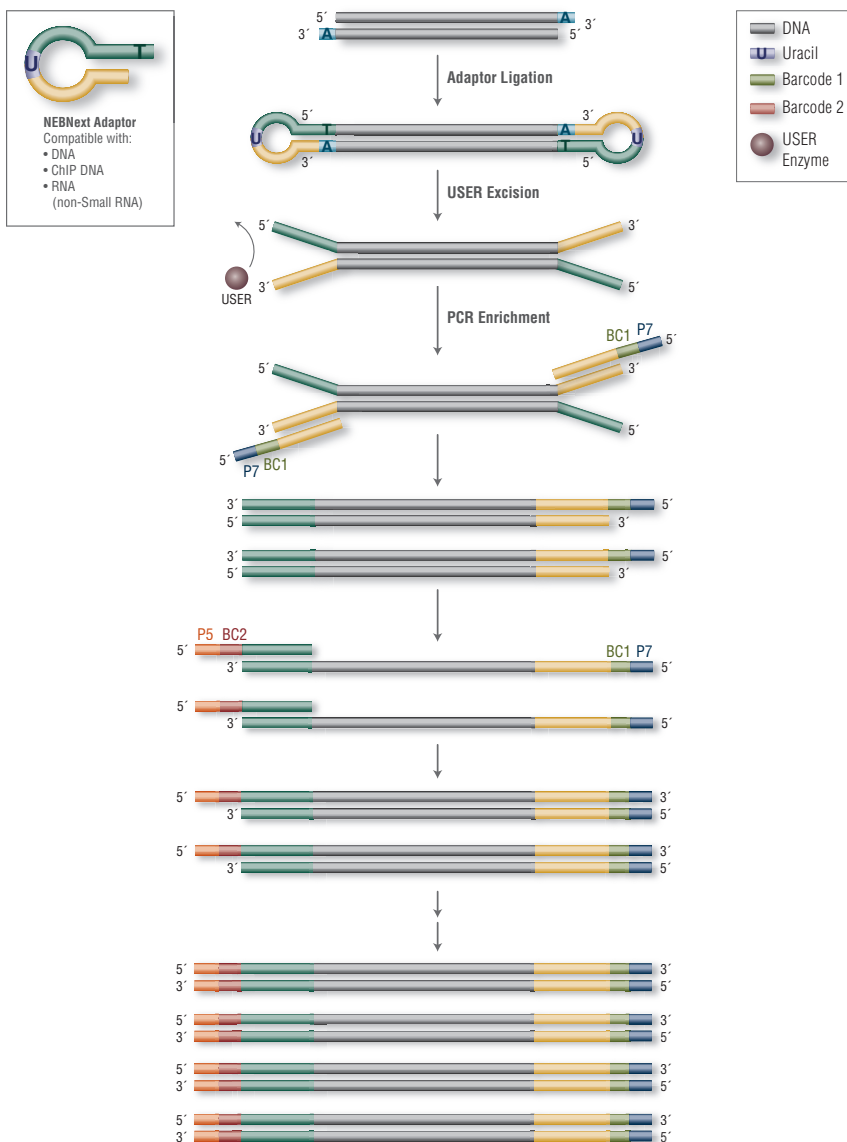
**Functionally Validated:** Each set of reagents is functionally validated together through construction and sequencing of genomic DNA libraries on the Illumina platform.

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 8-base index primers included in this kit are supplied in tubes with spare caps. 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 (Dual Index Primers).**



## 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](http://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 Reaction

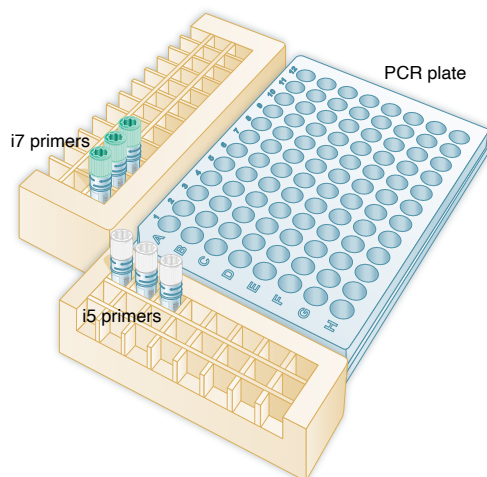
#### 1.1 PCR Amplification



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)

**Note:** We recommend using a PCR work-up rack such as the TruSeq® Index Plate Fixture (Illumina #FC-130-1005) to assist in properly combining the index primers during the PCR amplification step. Alternatively, 96-well deep well plates can be used and aligned against a PCR plate as in the diagram below.

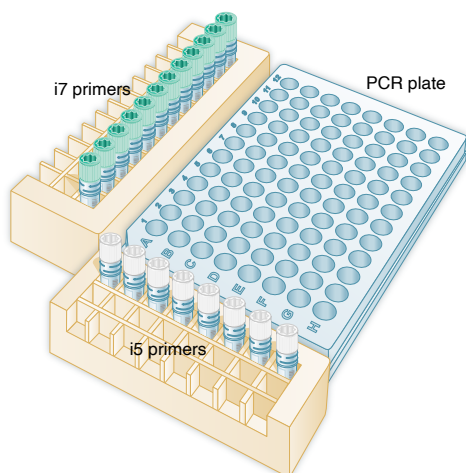


- 1.1A.1. Ensure that a valid combination of i7 and i5 primers is used. See Appendix A to verify that correct primer combinations have been selected.
- 1.1A.2. Arrange the index primers in the Index Plate Fixture as follows:
  - a. Arrange the ● (green) i7 primers in increasing order horizontally, so that the lowest number i7 index primer is in column 1, second lowest number i7 index primer is in column 2, etc.
  - b. Arrange the ○ (white) i5 primers in increasing order vertically, so that the lowest number i5 index primer is in row A, second lowest number i5 index primer is in row B, etc.
  - c. Record their positions on the PCR setup template (see Appendix B).
- 1.1A.3. Using a multichannel pipette, add desired volume of ○ (white) i5 primers to every column (as needed) of the PCR plate. **It is critical to change tips between columns to avoid cross-contamination.**
- 1.1A.4. Discard the original i5 white caps and apply new caps to avoid index cross-contamination.
- 1.1A.5. Using a multichannel pipette, add desired volume of ● (green) i7 primers to every row (as needed) of the PCR plate. **It is critical to change tips between rows to avoid cross-contamination.**
- 1.1A.6. Discard the original i7 green caps and apply new caps to avoid index cross-contamination.
- 1.1A.7. Add 25 µl ● (blue) NEBNext PCR Master Mix to each well that contains primers.

- 1.1A.8. Add desired volume of adaptor ligated DNA for a final volume of 50  $\mu$ l to the corresponding well. Gently pipette up and down 5–10 times to mix. **It is critical to change tips between samples to avoid cross-contamination.** Record each sample position on the PCR setup template (see Appendix B).
- 1.1A.9. Cover the plate with Bio-Rad<sup>®</sup> Microseal<sup>®</sup> "A" Film, and seal with a rubber roller. Quickly centrifuge.
- 1.1A.10. Perform PCR according to recommended cycling conditions.

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

**Note: We recommend using a PCR work-up rack such as the TruSeq Index Plate Fixture (Illumina #FC-130-1005) to assist in properly combining the index primers during the PCR amplification step. Alternatively, 96-well deep well plates can be used and aligned against a PCR plate as in the diagram below.**



- 1.1B.1. Arrange the index primers in the Index Plate Fixture as follows:
  - a. Arrange ● (green) i7 primers in increasing order horizontally, so that i713 is in column 1, i714 is in column 2, i715 is in column 3, etc.
  - b. Arrange the ○ (white) i5 primers in increasing order vertically, so that i509 is in row A, i510 is in row B, i511 is in row C, etc.
  - c. Record their positions on the PCR setup template (see Appendix B).
- 1.1B.2. Using a multichannel pipette, add desired volume of ○ (white) i5 primers to every column of the PCR plate. **It is critical to change tips between columns to avoid cross-contamination.**
- 1.1B.3. Discard the original i5 white caps and apply new caps to avoid index cross-contamination.
- 1.1B.4. Using a multichannel pipette, add desired volume of ● (green) i7 primers to every row of the PCR plate. **It is critical to change tips between rows to avoid cross-contamination.**
- 1.1B.5. Discard the original i7 green caps and apply new caps to avoid index cross-contamination.
- 1.1B.6. Add 25  $\mu$ l ● (blue) NEBNext PCR Master Mix to each well.
- 1.1B.7. Add desired volume of adaptor ligated DNA for a final volume of 50  $\mu$ l to the corresponding well. Gently pipette up and down 5–10 times to mix. **It is critical to change tips between samples to avoid cross-contamination.** Record each sample position on the PCR setup template (see Appendix B).
- 1.1B.8. Cover the plate with Bio-Rad Microseal "A" Film, and seal with a rubber roller. Quickly centrifuge.
- 1.1B.9. Perform PCR according to recommended cycling conditions.

## Section 2

### Appendix A: Principle for Use and Pooling Guide

#### The Principle of Dual Index Primers

The dual index primer strategy utilizes two 8 base indices within each primer; i7 primers contain indices that are adjacent to the P7 sequence; i5 primers contain indices that are adjacent to the P5 sequence. Dual indexing is enabled by adding a unique index to both ends of a sample to be sequenced. Up to 96 different samples can be uniquely indexed by combining each of the 12 i7 primers with each of the 8 i5 primers. Similarly, < 96 samples can be uniquely indexed by combining i7 primers with i5 primers as follows:

$$N = \text{Number of samples} = X(i7) * Y(i5) + \text{other primers as needed}$$

#### Examples:

##### 1. For N = 12 samples

Option 1: 4 (i7) \* 3 (i5)

From the i7 primers, choose a valid set of 4. From the i5 primers choose a valid set of 3. Use each i7 primer with each i5 primer to form 12 primer pairs for PCR amplification of 12 libraries. When setting up the sequencing run, select “Dual Index” and choose the indices used for each sample.

Option 2: 3 (i7) \* 4 (i5)

From the i7 primers, choose a valid set of 3. From the i5 primers choose a valid set of 4. Use each i7 primer with each i5 primer to form 12 primer pairs for PCR amplification of 12 libraries. When setting up the sequencing run, select “Dual Index” and choose the indices used for each sample.

Option 3: 12 (i7) \* 1 (i5)

Use all 12 i7 primers. Use any i5 primer. Use each i7 primer with the i5 primer to form 12 primer pairs for PCR amplification of 12 libraries. When setting up the sequencing run, select “Single Index”, and choose the i7 index used for each sample.

##### 1. For N = 26 samples

Option 1: 6 (i7) \* 4 (i5) + 2 (i5)

From the i7 primers, choose a valid set of 4 and add any other two i7 primers, for a total of 6 primers. From the i5 primers choose a valid set of 4 and add any other two i5 primers, for a total of 6 primers. Use each i7 primer with four of the i5 primers to form 24 primer pairs. Use any of the six i7 primers with the remaining two i5 primers to form 2 primer pairs. This will give you a total of 26 primer pairs for PCR amplification of 26 libraries. When setting up the sequencing run, select “Dual Index” and choose the indices used for each sample.

Option 2: 6 (i7) \* 5 (i5)

From the i7 primers, choose a valid set of 4 and add any other two i7 primers, for a total of 6 primers. From the i5 primers choose a valid set of 4 and add any other one i5 primer. Use each i7 primer with each i5 primer to form 30 primer pairs for PCR amplification. Use 26 of the 30 primer pairs to amplify 26 libraries. When setting up the sequencing run, select “Dual Index” and choose the indices used for each sample.

#### Low Plexity Pooling Guidelines

Illumina uses a red laser/LED to sequence A/C and a green laser/LED to sequence G/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. Table 2.1 and 2.2 list some valid combinations that can be sequenced together. Note: for 1-plex (no pooling), use any i7 primer with any i5 primer. In this case it is important to select “0” index reads in the Illumina Experiment Manager.

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.

**CAUTION: Sufficient primers are provided to generate 96 different samples if each i5 primer is used only once with each i7 primer.**

**If using subsets of i5 and i7 primers multiple times, you may have to readjust primer pairs to be able to generate 96 samples.**

**Table 2.1 Pooling: 2–12 libraries; Sequencing Workflow: Single Index  
(Select "1" Index Reads in the Illumina Experiment Manager).**

PLEX	i7 PRIMERS	i5 PRIMERS
2	i713 and i721 i714 and i724 i715 and i716 i717 and i719 i718 and i720 i722 and i723	Any i5 Primer
3	i714, i715 and i716 i715, i716 and i717 i717, i718 and i719 i718, i719 and i720 i721, i722 and i723 i722, i723 and i724	Any i5 Primer
4	i713, i714, i715 and i716 i715, i716, i717 and i718 i717, i718, i719 and i720 i718, i719, i720 and i721 i720, i721, i722 and i723 i721, i722, i723 and i724	Any i5 Primer
5–12	Any valid i7 4-plex with any other i7 Primers	Any i5 Primer

**Table 2.2 Pooling: 6+ libraries; Sequencing Workflow: Dual Index  
(Select "2" Index Reads in the Illumina Experiment Manager).**

PLEX	i7 PRIMERS	i5 PRIMERS
6–12	Any 2 plex combination from Table 2.1 with any other i7 primer(s) (as needed)	i509, i511 and i513 i510, i511 and i513 i511, i513 and i514 i511, i513 and i515 i511, i515 and i516
Greater than 12	Any 3 plex combination from the Table 2.1 with any other i7 primer(s) (as needed)	i509, i510, i511, i513 and any other i5 primer (as needed) i509, i511, i512, i513 and any other i5 primer (as needed) i509, i511, i514, i515 and any other i5 primer (as needed) i510, i511, i515, i516 and any other i5 primer (as needed) i511, i513, i515, i516 and any other i5 primer (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 (paired and 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).

**Tables 2.1 and 2.2 do not include an extensive list of all valid index combinations. Please check the sequences of each index to be used to ensure that you will have signal in both the red and green channels for every cycle. See example below:**

**Table 3.1**

GOOD																									
i7 PRIMERS				i5 PRIMERS																					
				FORWARD STRAND WORKFLOW*		REVERSE STRAND WORKFLOW*																			
i7-PCR Index 13	T	T	C	C	T	C	C	T	i5-PCR Index 9	T	T	G	C	T	T	G	C	G	C	A	A	G	C	A	A
i7-PCR Index 14	T	G	C	T	T	G	C	T	i5-PCR Index 10	G	A	G	A	G	G	T	T	A	A	C	C	T	C		C
i7-PCR Index 15	G	G	T	G	A	T	G	A	i5-PCR Index 11	A	C	C	T	G	G	T	T	A	A	C	C	A	G	G	T
i7-PCR Index 16	A	A	C	C	T	A	C	G	i5-PCR Index 13	C	G	G	A	A	C	A	A	T	T	G	T	T	C	C	G
	✓	✓	✓	✓	✓	✓	✓	✓		✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓	✓

**Table 3.2**

BAD																									
i7 PRIMERS				i5 PRIMERS																					
				FORWARD STRAND WORKFLOW *		REVERSE STRAND WORKFLOW *																			
i7-PCR Index 13	T	T	C	C	T	C	C	T	i5-PCR Index 9	T	T	G	C	T	T	G	C	G	C	A	A	G	C	A	A
i7-PCR Index 14	T	G	C	T	T	G	C	T	i5-PCR Index 10	G	A	G	A	G	G	T	T	A	A	C	C	T	C	T	C
i7-PCR Index 15	G	G	T	G	A	T	G	A	i5-PCR Index 12	A	A	G	C	G	G	A	A	T	T	C	C	G	C	T	T
i7-PCR Index 16	A	A	C	C	T	A	C	G	i5-PCR Index 14	G	G	T	A	A	G	C	T	A	G	C	T	T	A	C	C
	✓	✓	✓	✓	✓	✓	✓	✓		✓	✓	X	X	✓	X	✓	✓	✓	✓	X	✓	X	X	✓	✓

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. Within the NEB #E7780 and NEB #E7600 NEBNext Multiplex Oligos for Illumina kits there are no such combinations, but we wanted to include this information here in case of combining these kits with oligos from other index kits.

### Section 3

### Appendix B: PCR Setup Template

For each well, record: 1. DNA Sample Name \_\_\_\_\_

2. Index Primer Pairs \_\_\_\_\_

	1	2	3	4	5	6	7	8	9	10	11	12
A	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /
B	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /
C	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /
D	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /
E	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /
F	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /
G	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /
H	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /	1. _____ 2. /



## NEBNext Adaptors and Primers for Illumina

For sample sheets please see NEB.com, E7780 Product Page, "Protocols, Manuals and Usage Guidelines" Tab, [Usage Guidelines](#)

### NEBNext i509 Primer–NEBNext i516 Primer

Description: 8 Index Primers (10 µM) are included for producing barcoded libraries.

NEB #	PRODUCT	INDEX PRIMER SEQUENCE	EXPECTED INDEX	
			Forward Strand Workflow*	Reverse Strand Workflow*
#E7781A	NEBNext i509 Primer	5'-AATGATACGGCGACCACCGAGATCTACACT <b>TTGCTTGC</b> ACACTCTTCCCTACACGACGCTCTCCGATC*T-3'	TTGCTTGC	GCAAGCAA
#E7782A	NEBNext i510 Primer	5'-AATGATACGGCGACCACCGAGATCTACAC <b>GAGAGGTT</b> ACACTCTTCCCTACACGACGCTCTCCGATC*T-3'	GAGAGGTT	AACCTCTC
#E7783A	NEBNext i511 Primer	5'-AATGATACGGCGACCACCGAGATCTACAC <b>ACCTGGTT</b> ACACTCTTCCCTACACGACGCTCTCCGATC*T-3'	ACCTGGTT	AACCAGGT
#E7784A	NEBNext i512 Primer	5'-AATGATACGGCGACCACCGAGATCTACAC <b>AAGCGGAA</b> ACACTCTTCCCTACACGACGCTCTCCGATC*T-3'	AAGCGGAA	TTCCGCTT
#E7785A	NEBNext i513 Primer	5'-AATGATACGGCGACCACCGAGATCTACAC <b>CGGAACAA</b> ACACTCTTCCCTACACGACGCTCTCCGATC*T-3'	CGGAACAA	TTGTTCCG
#E7786A	NEBNext i514 Primer	5'-AATGATACGGCGACCACCGAGATCTACAC <b>GGTAAGCT</b> ACACTCTTCCCTACACGACGCTCTCCGATC*T-3'	GGTAAGCT	AGCTTACC
#E7787A	NEBNext i515 Primer	5'-AATGATACGGCGACCACCGAGATCTACAC <b>TGTGGCAT</b> ACACTCTTCCCTACACGACGCTCTCCGATC*T-3'	TGTGGCAT	ATGCCACA
#E7788A	NEBNext i516 Primer	5'-AATGATACGGCGACCACCGAGATCTACAC <b>ACTACGGA</b> ACACTCTTCCCTACACGACGCTCTCCGATC*T-3'	ACTACGGA	TCCGTAGT

\*If you are sequencing on a NovaSeq 6000 with v1.0 reagent kits, MiniSeq with Rapid reagent kits, MiSeq, HiSeq 2000/2500 (paired-end flow cell), HiSeq 3000/4000 (single-read flow cell), please follow sequence guidelines/sample sheet selection for **Forward Strand Workflow**. If you are sequencing on a iSeq 100, MiniSeq with Standard reagent kits, NextSeq Systems, NovaSeq 6000 with v1.5 reagent kits, HiSeq 2000/2500 (single-read flow cell), HiSeq 3000/4000 (paired-end flow cell), please follow sequence guidelines/ sample sheet selection for **Reverse Strand Workflow**.

### NEBNext i713 Primer–NEBNext i724 Primer

Description: 12 Index Primers (10 µM) are included for producing barcoded libraries.

NEB #	PRODUCT	INDEX PRIMER SEQUENCE	EXPECTED INDEX READ
#E7789A	NEBNext i713 Primer	5'-CAAGCAGAAGACGGCATAACGAGAT <b>AGGAGGA</b> AGTGACTGGAGTTCAGACGTGTGCTCTCCGATC*T-3'	TTCCTCCT
#E7790A	NEBNext i714 Primer	5'-CAAGCAGAAGACGGCATAACGAGAT <b>AGCAAGCA</b> GTGACTGGAGTTCAGACGTGTGCTCTCCGATC*T-3'	TGCTTGCT
#E7791A	NEBNext i715 Primer	5'-CAAGCAGAAGACGGCATAACGAGAT <b>TCATCAC</b> CGTGACTGGAGTTCAGACGTGTGCTCTCCGATC*T-3'	GGTGATGA
#E7792A	NEBNext i716 Primer	5'-CAAGCAGAAGACGGCATAACGAGAT <b>CGTAGGTT</b> GTGACTGGAGTTCAGACGTGTGCTCTCCGATC*T-3'	AACCTACG
#E7793A	NEBNext i717 Primer	5'-CAAGCAGAAGACGGCATAACGAGAT <b>TCAGATCC</b> GTGACTGGAGTTCAGACGTGTGCTCTCCGATC*T-3'	GGATCTGA
#E7794A	NEBNext i718 Primer	5'-CAAGCAGAAGACGGCATAACGAGAT <b>CGTGATCA</b> GTGACTGGAGTTCAGACGTGTGCTCTCCGATC*T-3'	TGATCACG
#E7795A	NEBNext i719 Primer	5'-CAAGCAGAAGACGGCATAACGAGAT <b>AGTCGCTT</b> GTGACTGGAGTTCAGACGTGTGCTCTCCGATC*T-3'	AAGCGACT
#E7796A	NEBNext i720 Primer	5'-CAAGCAGAAGACGGCATAACGAGAT <b>GAACGCTT</b> GTGACTGGAGTTCAGACGTGTGCTCTCCGATC*T-3'	AAGCGTTC
#E7797A	NEBNext i721 Primer	5'-CAAGCAGAAGACGGCATAACGAGAT <b>TACGCCTT</b> GTGACTGGAGTTCAGACGTGTGCTCTCCGATC*T-3'	AAGGCGTA
#E7798A	NEBNext i722 Primer	5'-CAAGCAGAAGACGGCATAACGAGAT <b>CTCATCAG</b> GTGACTGGAGTTCAGACGTGTGCTCTCCGATC*T-3'	CTGATGAG

#E7799A	NEBNext i723 Primer	5'-CAAGCAGAAGACGGCATACGAGATTCTTCTGCGTGACTGGAGTTCAGACGTGTGCTCTCCGATC*T-3'	GCAGAAGA
#E7800A	NEBNext i724 Primer	5'-CAAGCAGAAGACGGCATACGAGATGCTGGATTGTGACTGGAGTTCAGACGTGTGCTCTCCGATC*T-3'	AATCCAGC

## Kit Components

The NEBNext Dual Index Primers Set 2 is functionally validated through library preparation using the NEBNext Library Prep Kits and sequencing on the Illumina platform.

### NEB #E7780S Table of Components

NEB #	PRODUCT NAME	VOLUME
E7601A	NEBNext Adaptor for Illumina*	0.96 ml
E7602A	USER Enzyme	0.288 ml
E7781A	NEBNext i509 Primer	0.060 ml
E7782A	NEBNext i510 Primer	0.060 ml
E7783A	NEBNext i511 Primer	0.060 ml
E7784A	NEBNext i512 Primer	0.060 ml
E7785A	NEBNext i513 Primer	0.060 ml
E7786A	NEBNext i514 Primer	0.060 ml
E7787A	NEBNext i515 Primer	0.060 ml
E7788A	NEBNext i516 Primer	0.060 ml
E7789A	NEBNext i713 Primer	0.040 ml
E7790A	NEBNext i714 Primer	0.040 ml
E7791A	NEBNext i715 Primer	0.040 ml
E7792A	NEBNext i716 Primer	0.040 ml
E7793A	NEBNext i717 Primer	0.040 ml
E7794A	NEBNext i718 Primer	0.040 ml
E7795A	NEBNext i719 Primer	0.040 ml
E7796A	NEBNext i720 Primer	0.040 ml
E7797A	NEBNext i721 Primer	0.040 ml
E7798A	NEBNext i722 Primer	0.040 ml
E7799A	NEBNext i723 Primer	0.040 ml
E7800A	NEBNext i724 Primer	0.040 ml

\*15 µM Concentration

## Revision History

REVISION #	DESCRIPTION	DATE
1.0	N/A	11/17
2.0	Updated manual format	8/22

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](mailto: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.

AGENCOURT® and AMPURE® are registered trademarks of Beckman Coulter, Inc.

BIO-RAD® and MICROSEAL® are registered trademarks of Bio-Rad Laboratories, Inc.

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

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



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



*be* INSPIRED  
*drive* DISCOVERY  
*stay* GENUINE

New England Biolabs, Inc., 240 County Road, Ipswich, MA 01938-2723 Telephone: (978) 927-5054 Toll Free: (USA Orders) 1-800-632-5227 (USA Tech) 1-800-632-7799 Fax: (978) 921-1350 e-mail: [info@neb.com](mailto:info@neb.com)