

NEBNext[®] Multiplex Oligos for Illumina[®] (Index Primers Set 3)

NEB #E7710S/L

24/96 reactions

Version 5.0_6/22

Table of Contents

NEBNext Adaptors and Primers for Illumina	4
Kit Components	5
Revision History	6

The NEBNext Multiplex Oligos for Illumina (Index Primers Set 3) Includes:

The volumes provided are sufficient for preparation of up to 24 reactions (NEB #E7710S) and 96 reactions (NEB #E7710L). All reagents should be stored at -20°C.

NEBNext Adaptor for Illumina

USER[®] Enzyme

NEBNext Universal PCR Primer for Illumina

NEBNext Index 17 Primer for Illumina

NEBNext Index 24 Primer for Illumina

NEBNext Index 26 Primer for Illumina

NEBNext Index 28 Primer for Illumina

NEBNext Index 29 Primer for Illumina

NEBNext Index 31 Primer for Illumina

NEBNext Index 32 Primer for Illumina

NEBNext Index 33 Primer for Illumina

NEBNext Index 34 Primer for Illumina

NEBNext Index 35 Primer for Illumina

NEBNext Index 36 Primer for Illumina

NEBNext Index 42 Primer for Illumina

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

Applications

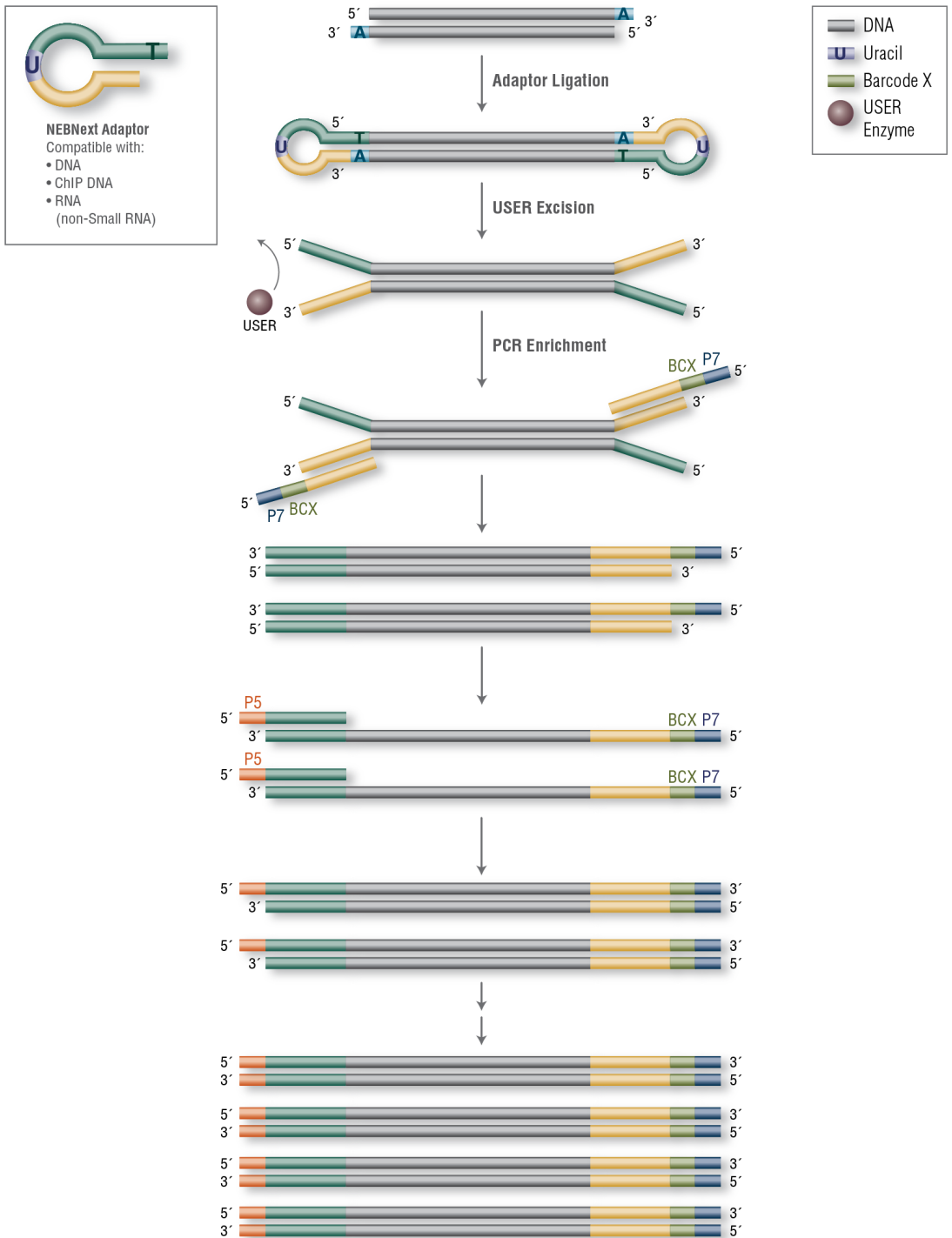
The NEBNext Multiplex Oligos for Illumina (Index Primers Set 3) contains adaptors and primers that are ideally suited for multiplex sample preparation for next generation sequencing on the Illumina platform (Illumina, Inc.). Each of these components must pass rigorous quality control standards and is lot controlled, both individually and as a set of reagents.

Lot Control: The lots provided in the NEBNext Multiplex Oligos for Illumina (Index Primers Set 3) 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 listed on each individual component page.

Functionally Validated: Each of the components is functionally validated through construction and sequencing of a 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.

Figure 1. Workflow demonstrating the use of NEBNext Multiplex Oligos for Illumina (Index Primers Set 3)



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

NEBNext Adaptors and Primers for Illumina

For sample sheets please see NEB.com, E7710 Product Page, "Protocols, Manuals and Usage Guidelines" Tab, [Usage Guidelines](#)

PRODUCT	INDEX PRIMER SEQUENCE	EXPECTED INDEX PRIMER SEQUENCE READ
NEBNext Index 17 Primer for Illumina (10 μM)	5'-CAAGCAGAAGACGGCATAACGAGAT <u>CTCTAC</u> GTGACTGGAGTTCAGACGTGTGCTCTTCCGATC-s-T-3'	GTAGAG
NEBNext Index 24 Primer for Illumina (10 μM)	5'-CAAGCAGAAGACGGCATAACGAGAT <u>GCTACC</u> GTGACTGGAGTTCAGACGTGTGCTCTTCCGATC-s-T-3'	GGTAGC
NEBNext Index 26 Primer for Illumina (10 μM)	5'-CAAGCAGAAGACGGCATAACGAGAT <u>GCTCAT</u> GTGACTGGAGTTCAGACGTGTGCTCTTCCGATC-s-T-3'	ATGAGC
NEBNext Index 28 Primer for Illumina (10 μM)	5'-CAAGCAGAAGACGGCATAACGAGAT <u>CTTTTG</u> GTGACTGGAGTTCAGACGTGTGCTCTTCCGATC-s-T-3'	CAAAAG
NEBNext Index 29 Primer for Illumina (10 μM)	5'-CAAGCAGAAGACGGCATAACGAGAT <u>TAGTTG</u> GTGACTGGAGTTCAGACGTGTGCTCTTCCGATC-s-T-3'	CAACTA
NEBNext Index 31 Primer for Illumina (10 μM)	5'-CAAGCAGAAGACGGCATAACGAGAT <u>ATCGTG</u> GTGACTGGAGTTCAGACGTGTGCTCTTCCGATC-s-T-3'	CACGAT
NEBNext Index 32 Primer for Illumina (10 μM)	5'-CAAGCAGAAGACGGCATAACGAGAT <u>TGAGTG</u> GTGACTGGAGTTCAGACGTGTGCTCTTCCGATC-s-T-3'	CACTCA
NEBNext Index 33 Primer for Illumina (10 μM)	5'-CAAGCAGAAGACGGCATAACGAGAT <u>CGCCTG</u> GTGACTGGAGTTCAGACGTGTGCTCTTCCGATC-s-T-3'	CAGGCG
NEBNext Index 34 Primer for Illumina (10 μM)	5'-CAAGCAGAAGACGGCATAACGAGAT <u>GCCATG</u> GTGACTGGAGTTCAGACGTGTGCTCTTCCGATC-s-T-3'	CATGGC
NEBNext Index 35 Primer for Illumina (10 μM)	5'-CAAGCAGAAGACGGCATAACGAGAT <u>AAAATG</u> GTGACTGGAGTTCAGACGTGTGCTCTTCCGATC-s-T-3'	CATTTT
NEBNext Index 36 Primer for Illumina (10 μM)	5'-CAAGCAGAAGACGGCATAACGAGAT <u>TGTTGG</u> GTGACTGGAGTTCAGACGTGTGCTCTTCCGATC-s-T-3'	CCAACA
NEBNext Index 42 Primer for Illumina (10 μM)	5'-CAAGCAGAAGACGGCATAACGAGAT <u>CGATTAG</u> GTGACTGGAGTTCAGACGTGTGCTCTTCCGATC-s-T-3'	TAATCG
NEBNext Adaptor for Illumina (15 μM)	5'-/5Phos/GAT CGG AAG AGC ACA CGT CTG AAC TCC AGT CUA CAC TCT TTC CCT ACA CGA CGC TCT TCC GAT C-s-T-3'	N/A
NEBNext Universal PCR Primer for Illumina (10 μM)	5'-AAT GAT ACG GCG ACC ACC GAG ATC TAC ACT CTT TCC CTA CAC GAC GCT CTT CCG ATC-s-T-3'	N/A

Where -s- indicates phosphorothioate bond.

Note: If fewer than 12 indexes are used in a lane for sequencing, it is recommended to use the following indexes:

Pool of 2 samples: Index #24 and 31; Index #26 and 42

Pool of 3 samples: Index #24, 32 and 33; Index #24, 34 and 42

Pool of 4 samples: Index #24, 32, 35 and 36; Index #17, 26, 36 and 42

To determine possible combinations for low plex pooling, please also see <https://indexoligo.neb.com/>

Kit Components

NEB #E7710S Table of Components

NEB #	PRODUCT	VOLUME
E7337A	NEBNext Adaptor for Illumina	0.24 ml
E7338A	USER Enzyme	0.072 ml
E6861A	NEBNext Universal PCR Primer for Illumina	0.120 ml
E7585A	NEBNext Index 17 Primer for Illumina	0.010 ml
E7592A	NEBNext Index 24 Primer for Illumina	0.010 ml
E7713A	NEBNext Index 26 Primer for Illumina	0.010 ml
E7714A	NEBNext Index 28 Primer for Illumina	0.010 ml
E7715A	NEBNext Index 29 Primer for Illumina	0.010 ml
E7717A	NEBNext Index 31 Primer for Illumina	0.010 ml
E7718A	NEBNext Index 32 Primer for Illumina	0.010 ml
E7719A	NEBNext Index 33 Primer for Illumina	0.010 ml
E7720A	NEBNext Index 34 Primer for Illumina	0.010 ml
E7721A	NEBNext Index 35 Primer for Illumina	0.010 ml
E7722A	NEBNext Index 36 Primer for Illumina	0.010 ml
E7736A	NEBNext Index 42 Primer for Illumina	0.010 ml

NEB #E7710L Table of Components

NEB #	PRODUCT	VOLUME
E7337AA	NEBNext Adaptor for Illumina	0.96 ml
E7338AA	USER Enzyme	0.288 ml
E6861AA	NEBNext Universal PCR Primer for Illumina	0.480 ml
E7585AA	NEBNext Index 17 Primer for Illumina	0.040 ml
E7592AA	NEBNext Index 24 Primer for Illumina	0.040 ml
E7713AA	NEBNext Index 26 Primer for Illumina	0.040 ml
E7714AA	NEBNext Index 28 Primer for Illumina	0.040 ml
E7715AA	NEBNext Index 29 Primer for Illumina	0.040 ml
E7717AA	NEBNext Index 31 Primer for Illumina	0.040 ml
E7718AA	NEBNext Index 32 Primer for Illumina	0.040 ml
E7719AA	NEBNext Index 33 Primer for Illumina	0.040 ml
E7720AA	NEBNext Index 34 Primer for Illumina	0.040 ml
E7721AA	NEBNext Index 35 Primer for Illumina	0.040 ml
E7722AA	NEBNext Index 36 Primer for Illumina	0.040 ml
E7736AA	NEBNext Index 42 Primer for Illumina	0.040 ml

Revision History

REVISION #	DESCRIPTION	DATE
1.0	N/A	6/16
1.1	Updated sequences for NEBNext Index Primers 17 and 24.	12/16
2.0	Delete "Required Materials Not Included for DNA or ChIP Libraries" and "Required Materials Not Included for RNA Libraries" and replace with guidance note. Update the list of kits using the NEBNext Multiplex Oligos for Illumina. Create "Kit Component – Table of Components" for small and large size kits. Change Index Primer sequences table to include adaptor, and NEBNext Universal PCR Primer and rename it as NEBNext Adaptors and Primers for Illumina. Delete individual component information pages. Updated the functionally validated statement on page 2.	1/18
3.0	Updated to new manual format.	1/20
4.0	Updated adaptors and primers table for Illumina	9/21
5.0	Updated adaptors and primers table for Illumina	6/22

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® is a registered trademark of Illumina, Inc.

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



www.neb.com



be INSPIRED
drive DISCOVERY
stay GENUINE