

# NEBNext<sup>®</sup> Ultra<sup>™</sup> II Ligation Module

## for use with NEBNext Multiplex Oligos for Illumina<sup>®</sup> (Unique Dual Index UMI Adaptors DNA Set 1, NEB #E7395)

NEB #E7595S/L

24/96 reactions

Version 1.0\_3/20

### Table of Contents

|  |   |
|--|---|
| Section 1  |   |
| Protocol for use with NEBNext Ultra II End Repair/dA-Tailing Module..... | 2 |
| Section 2  |   |
| Protocol for use with NEBNext Ultra II FS DNA Module .....               | 4 |
| Kit Components .....   | 5 |
| Revision History .....   | 5 |

### The NEBNext Ultra II Ligation Module Includes

*The volumes provided are sufficient for preparation of up to 24 reactions (NEB #E7595S) and 96 reactions (NEB #E7595L). All reagents should be stored at –20°C. Colored bullets represent the color of the cap of the tube containing the reagent.*

- (red) NEBNext Ultra II Ligation Master Mix
- (red) NEBNext Ligation Enhancer

### The NEBNext Ultra II Ligation Module is Designed for use with the Following

NEBNext Multiplex Oligos for Illumina<sup>®</sup> (Unique Dual Index UMI Adaptors DNA Set 1, #E7395)

NEBNext Ultra II End Repair/dA-Tailing Module (NEB #E7546)

NEBNext Ultra II FS DNA Module (NEB #E7810)

NEBNext Ultra II Q5<sup>®</sup> Master Mix (NEB #M0544)

### Applications

The NEBNext Ultra II Ligation Module is optimized for use with the NEBNext Ultra II End Repair/dA-Tailing Module (NEB #E7546) or the NEBNext Ultra II FS DNA Module (NEB #E7810).

Each module component must pass rigorous quality control standards, and for each new lot the entire set of reagents is functionally validated together with NEB #E7546 or NEB #E7810 along with NEB #M0544 to construct indexed libraries that are sequenced 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](mailto:OEM@neb.com) for further information.

## Section 1

### Protocol for use with NEBNext Ultra II End Repair/dA-Tailing Module (NEB #E7546)

#### Symbols



This caution sign signifies a step in the protocol that has two paths leading to the same end point but is dependent on a user variable, like the amount of input DNA.



This is a point where you can safely stop the protocol and store the samples prior to proceeding to the next step in the protocol.



Colored bullets indicate the cap color of the reagent to be added.

#### Starting Material

500 pg–1 µg fragmented DNA that has been end-repaired and dA-tailed using the NEBNext Ultra II End Repair/dA-Tailing Module (NEB #E7546).



If DNA input is  $\leq 100$  ng, dilute the NEBNext UMI Adaptors for Illumina in the UMI Dilution Buffer provided in the kit as indicated in Table 1.1.

**Note: The appropriate adaptor dilution for your sample input and type may need to be optimized experimentally. The dilutions provided here are a general starting point.**

**Table 1.1: Adaptor Dilution.**

| INPUT          | ADAPTOR DILUTION<br>(VOLUME OF ADAPTOR: TOTAL VOLUME) | WORKING ADAPTOR<br>CONCENTRATION |
|----------------|---|----------------------------------|
| 1 µg–101 ng    | No Dilution   | 20 µM                            |
| 100 ng–5 ng    | 10-Fold (1:10)  | 2 µM                             |
| less than 5 ng | 50-Fold (1:50)  | 0.4 µM                           |

1.1. Add the following components directly to the End Prep Reaction Mixture:

| COMPONENT                                      | VOLUME (µl) PER REACTION |
|--|--------------------------|
| End Prep Reaction Mixture                      | 60 µl                    |
| NEBNext UMI Adaptors for Illumina*             | 2.5 µl                   |
| • (red) NEBNext Ultra II Ligation Master Mix** | 30 µl                    |
| • (red) NEBNext Ligation Enhancer              | 1 µl                     |
| Total Volume                                   | 93.5 µl                  |

\* The NEBNext UMI adaptors are provided in NEBNext Multiplex Oligos for Illumina (Unique Dual Index UMI Adaptors DNA Set 1, NEB #E7395). Please refer to the NEB #E7395 manual for valid barcode combinations.

\*\* Mix the Ultra II Ligation Master Mix by pipetting up and down several times prior to adding to the reaction.

**Note: The Ligation Master Mix and Ligation Enhancer can be mixed ahead of time and is stable for at least 8 hours @ 4°C. We do not recommend premixing the Ligation Master Mix, Ligation Enhancer and adaptor prior to use in the Adaptor Ligation Step.**

1.2. Set a 100 µl or 200 µl pipette to 80 µl and then pipette the entire volume up and down at least 10 times to mix thoroughly. Perform a quick spin to collect all liquid from the sides of the tube.

**(Caution: The NEBNext Ultra II Ligation Master Mix is very viscous. Care should be taken to ensure adequate mixing of the ligation reaction, as incomplete mixing will result in reduced ligation efficiency. The presence of a small amount of bubbles will not interfere with performance).**

1.3. Incubate at 20°C for 15 minutes in a thermal cycler with the heated lid off.

1.4. DNA is now ready for size selection or cleanup.

**Note: Please see NEB #E7645 manual for recommended size selection/cleanup and PCR amplification protocols.**



**Samples can be stored overnight at  $-20^{\circ}\text{C}$ .**

## Section 2

### Protocol for use with NEBNext Ultra II FS DNA Module (NEB #E7810)

#### Symbols



This caution sign signifies a step in the protocol that has two paths leading to the same end point but is dependent on a user variable, like the amount of input DNA.



This is a point where you can safely stop the protocol and store the samples prior to proceeding to the next step in the protocol.



Colored bullets indicate the cap color of the reagent to be added.

#### Starting Material

100 pg–500 ng fragmented, end repaired and dA-Tailed DNA generated using the NEBNext Ultra II FS DNA Module (NEB #E7810).



If DNA input is  $\leq 100$  ng, dilute the NEBNext UMI Adaptors for Illumina in the UMI Dilution Buffer provided in the kit as indicated in Table 2.1.

**Note: The appropriate adaptor dilution for your sample input and type may need to be optimized experimentally. The dilutions provided here are a general starting point.**

**Table 2.1 Adaptor Dilution.**

| INPUT          | ADAPTOR DILUTION<br>(VOLUME OF ADAPTOR: TOTAL VOLUME) | WORKING ADAPTOR<br>CONCENTRATION |
|----------------|---|----------------------------------|
| 100 ng–500 ng  | No Dilution   | 20 $\mu$ M                       |
| 5 ng–99 ng     | 10-Fold (1:10)  | 2 $\mu$ M                        |
| 1 ng–5 ng      | 50-Fold (1:50)  | 0.4 $\mu$ M                      |
| less than 1 ng | 100-Fold (1:100)                                      | 0.2 $\mu$ M                      |

2.1. Add the following components directly to the FS Reaction Mixture

| COMPONENT                                      | VOLUME ( $\mu$ l) PER REACTION |
|--|--------------------------------|
| FS Reaction Mixture                            | 35 $\mu$ l                     |
| NEBNext UMI Adaptors for Illumina*             | 2.5 $\mu$ l                    |
| • (red) NEBNext Ultra II Ligation Master Mix** | 30 $\mu$ l                     |
| • (red) NEBNext Ligation Enhancer              | 1 $\mu$ l                      |
| Total Volume                                   | 68.5 $\mu$ l                   |

\* The NEBNext UMI adaptors are provided in NEBNext Multiplex Oligos for Illumina (Unique Dual Index UMI Adaptors DNA Set 1, NEB #E7395). Please refer to the NEB #E7395 manual for valid barcode combinations.

\*\* Mix the Ultra II Ligation Master Mix by pipetting up and down several times prior to adding to the reaction.

**Note: The Ligation Master Mix and Ligation Enhancer can be mixed ahead of time and is stable for at least 8 hours @ 4°C. We do not recommend premixing the Ligation Master Mix, Ligation Enhancer and adaptor prior to use in the Adaptor Ligation Step.**

2.2. Set a 100  $\mu$ l or 200  $\mu$ l pipette to 50  $\mu$ l and then pipette the entire volume up and down at least 10 times to mix thoroughly. Perform a quick spin to collect all liquid from the sides of the tube.

**(Caution: The NEBNext Ultra II Ligation Master Mix is very viscous. Care should be taken to ensure adequate mixing of the ligation reaction, as incomplete mixing will result in reduced ligation efficiency. The presence of a small amount of bubbles will not interfere with performance).**

- 2.3. Incubate at 20°C for 15 minutes in a thermal cycler with the heated lid off.
- 2.4. DNA is now ready for size selection or cleanup.

**Note: Please see NEB #E7805 manual for recommended size selection/cleanup and PCR amplification protocols.**



Samples can be stored overnight at –20°C.

## Kit Components

Each set of reagents is functionally validated together with NEB #E7546 or NEB #E7810 along with NEB #M0544 to construct index libraries that are sequenced on an Illumina sequencing platform.

### NEB #E7595S Table of Components

| NEB #  | PRODUCT                              | VOLUME   |
|--------|--------------------------------------|----------|
| E7648A | NEBNext Ultra II Ligation Master Mix | 0.72 ml  |
| E7374A | NEBNext Ligation Enhancer            | 0.024 ml |

### NEB #E7595L Table of Components

| NEB #   | PRODUCT                              | VOLUME       |
|---------|--------------------------------------|--------------|
| E7648AA | NEBNext Ultra II Ligation Master Mix | 3 x 0.960 ml |
| E7374AA | NEBNext Ligation Enhancer            | 0.096 ml     |

## Revision History

| REVISION # | DESCRIPTION | DATE |
|------------|-------------|------|
| 1.0        | N/A         | 3/20 |

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. 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 the buyer to obtain additional third party intellectual property rights for certain applications.

B CORPORATION™ is a trademark of B Lab, Inc.

ILLUMINA® is a registered trademark of Illumina, Inc.

© Copyright 2020, New England Biolabs, Inc.; all rights reserved



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



be INSPIRED  
drive DISCOVERY  
stay GENUINE