

# LIBRARY PREPARATION

## NEBNext<sup>®</sup> Ultra<sup>™</sup> II End Repair/dA-Tailing Module

Instruction Manual

NEB #E7546S/L  
24/96 reactions  
Version 2.0 4/18



NEW ENGLAND  
**BioLabs**<sup>®</sup> Inc.

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

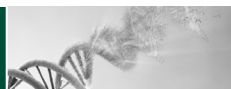
This product is intended for research purposes only. This product is not intended to be used for therapeutic or diagnostic purposes in humans or animals.



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 [gbd@neb.com](mailto:gbd@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.

ILLUMINA® is a registered trademark of Illumina, Inc.

© Copyright 2018, New England Biolabs, Inc; all rights reserved.



## Table of Contents:

---

Applications.....	2
Protocol.....	3
Kit Components.....	4
Revision History.....	5

## The NEBNext Ultra End Repair/dA-Tailing Module Includes:

---

*The volumes provided are sufficient for preparation of up to 24 reactions (NEB #E7546S) and 96 reactions (NEB #E7546L). All reagents should be stored at -20°C. Colored bullets represent the color of the cap of the tube containing the reagent.*

- (green) NEBNext Ultra II End Prep Reaction Buffer
- (green) NEBNext Ultra II End Prep Enzyme Mix

## The NEBNext Ultra II End Repair/dA-Tailing Module Is Designed For Use with the Following:

---

NEBNext Singleplex or Multiplex Oligos for Illumina®

(NEB #E7350, #E7335, #E7500, #E7600 or #E7535)

NEBNext Ultra II Ligation Module (NEB #E7595)

NEBNext Ultra II Q5® Master Mix (NEB #M0544)

## Applications:

The NEBNext Ultra II End Repair/dA-Tailing Module is optimized to convert 500 pg-1  $\mu$ g of fragmented DNA to repaired DNA having 5' phosphorylated, 3' dA-tailed ends.


**Lot Control:** The lots provided in the NEBNext Ultra II End Repair/dA-Tailing Module for Illumina are managed separately and 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 set of reagents is functionally validated together through construction and sequencing of an indexed DNA library on the 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.

# Protocol:

## 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.

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

 Stopping points in the protocol.

**Starting Material:** 500 pg–1 µg fragmented DNA. We recommend that DNA be sheared in 1X TE. If the DNA volume post shearing is less than 50 µl, add 1X TE to a final volume of 50 µl. Alternatively, 10 mM Tris-HCl, pH 8.0 or 0.1X TE can be used.

### 1.1 NEBNext End Prep

1. Add the following components to a sterile nuclease-free tube:

● (green) NEBNext Ultra II End Prep Enzyme Mix	3 µl
● (green) NEBNext Ultra II End Prep Reaction Buffer	7 µl
Fragmented DNA	50 µl
<hr/>	
Total volume	60 µl

2. Set a 100 µl or 200 µl pipette to 50 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.


**Note: It is important to mix well. The presence of a small amount of bubbles will not interfere with performance.**

3. Place in a thermocycler, with the heated lid set to  $\geq 75^{\circ}\text{C}$ , and run the following program:

30 minutes @  $20^{\circ}\text{C}$

30 minutes @  $65^{\circ}\text{C}$

Hold at  $4^{\circ}\text{C}$

 **If necessary, samples can be stored at  $-20^{\circ}\text{C}$ ; however, a slight loss in yield (~20%) may be observed. We recommend continuing with adaptor ligation before stopping.**

4. Proceed directly to NEBNext Ultra II Ligation Module (NEB #E7595) (available in September 2015).

# Kit Components

## NEB #E7546S Table of Components

NEB #	PRODUCT	VOLUME
E7646A	NEBNext Ultra II End Prep Enzyme Mix	0.072 ml
E7647A	NEBNext Ultra II End Prep Reaction Buffer	0.168 ml

## NEB #E7546L Table of Components

NEB #	PRODUCT	VOLUME
E7646AA	NEBNext Ultra II End Prep Enzyme Mix	0.288 ml
E7647AA	NEBNext Ultra II End Prep Reaction Buffer	0.672 ml

## Revision History:

REVISION #	DESCRIPTION	DATE
1.0	N/A	8/15
2.0	Create "Kit Component – Table of Components" for small and large size kits. Delete individual component information pages.	4/18

DNA CLONING  
DNA AMPLIFICATION & PCR  
EPIGENETICS  
RNA ANALYSIS  
LIBRARY PREP FOR NEXT GEN SEQUENCING  
PROTEIN EXPRESSION & ANALYSIS  
CELLULAR ANALYSIS



## USA

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

## CANADA

New England Biolabs, Ltd.  
Telephone: (905) 665-4632  
Toll Free: 1-800-387-1095  
Fax: (905) 665-4635  
Fax Toll Free: 1-800-563-3789  
e-mail: [info.ca@neb.com](mailto:info.ca@neb.com)  
[www.neb.ca](http://www.neb.ca)

## CHINA

New England Biolabs (Beijing), Ltd.  
Telephone: 010-82378265/82378266  
Fax: 010-82378262  
e-mail: [info@neb-china.com](mailto:info@neb-china.com)  
[www.neb-china.com](http://www.neb-china.com)

## FRANCE

New England Biolabs France  
Free Call: 0800-100-632  
Free Fax: 0800-100-610  
e-mail: [info.fr@neb.com](mailto:info.fr@neb.com)  
[www.neb-online.fr](http://www.neb-online.fr)

## GERMANY & AUSTRIA

New England Biolabs GmbH  
Telephone: +49/(0)69/305 23140  
Free Call: 0800/246 5227 (Germany)  
Free Call: 00800/246 52277 (Austria)  
Fax: +49/(0)69/305 23149  
Free Fax: 0800/246 5229 (Germany)  
e-mail: [info.de@neb.com](mailto:info.de@neb.com)  
[www.neb-online.de](http://www.neb-online.de)

## JAPAN

New England Biolabs Japan, Inc.  
Telephone: +81 (0)3 5669 6191  
Fax: +81 (0)3 5669 6192  
e-mail: [info.jp@neb.com](mailto:info.jp@neb.com)  
[www.nebj.jp](http://www.nebj.jp)

## SINGAPORE

New England Biolabs Pte. Ltd.  
Telephone: +65 638 59623  
Fax: +65 638 59617  
e-mail: [sales.sg@neb.com](mailto:sales.sg@neb.com)  
[www.neb.sg](http://www.neb.sg)

## UNITED KINGDOM

New England Biolabs (UK) Ltd.  
Telephone: (01462) 420616  
Call Free: 0800 318486  
Fax: (01462) 421057  
Fax Free: 0800 435682  
e-mail: [info.uk@neb.com](mailto:info.uk@neb.com)  
[www.neb.uk.com](http://www.neb.uk.com)