

NEBNext[®] Ultra[™] II Directional RNA Second Strand Synthesis Module

NEB #E7550S/L

24/96 reactions

Version 5.0_1/20

Table of Contents

Protocol	2
Kit Components	3
Revision History	3

The NEBNext Ultra II Directional RNA Second Strand Synthesis Module Includes

The volumes provided are sufficient for preparation of up to 24 reactions (NEB #E7550S) and 96 reactions (NEB #E7550L). All reagents should be stored at -20°C . Colored bullets represent the color of the cap of the tube containing the reagent.

- (orange) NEBNext Second Strand Synthesis Enzyme Mix
- (orange) NEBNext Second Strand Synthesis Reaction Buffer with dUTP Mix (10X)

The NEBNext Ultra II Directional RNA Second Strand Synthesis Module is Designed for use with the Following:

NEBNext Ultra II RNA First Strand Synthesis Module (NEB #E7771) or
 NEBNext RNA First Strand Synthesis Module (NEB #E7525)

NEBNext Singleplex (NEB #E7350) or
 NEBNext Multiplex (NEB #E7335, #E7500, #E7710, #E7730, #E6609 or #E7600) Oligos for Illumina

NEBNext Ultra II End Repair/dA-Tailing Module (NEB #E7546) or
 NEBNext Ultra End Repair/dA-Tailing Module (NEB #E7442)

NEBNext Ultra II Ligation Module (NEB #E7595) or
 NEBNext Ultra Ligation Module (NEB #E7445)

NEBNext Ultra II Q5 Master Mix (NEB #M0544) or
 NEBNext High Fidelity 2X PCR Master Mix (NEB #M0541)

USER[®] Enzyme (NEB #M5505)

Note: This is included in the NEBNext Oligo sets; however, if you are using alternative oligos, it is necessary to purchase separately for the directional method.

Overview

The NEBNext Ultra II Directional RNA Second Strand Synthesis Module has been optimized to generate double-stranded cDNA from first-strand cDNA using the NEBNext Ultra II RNA First Strand Synthesis Module (NEB #E7771) or NEBNext RNA First Strand Synthesis Module (NEB #E7525). The dsDNA generated by the NEBNext Ultra II Directional RNA Second Strand Module can be subsequently converted to blunt-end DNA fragments using the NEBNext Ultra II End Repair/dA-Tailing Module (NEB #E7546).

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 #E7111, #E7546, #E7595 and #M0544 to construct an indexed transcriptome library that is 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 for further information.

Protocol

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 type of RNA input.



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.

Note: This protocol has changed to be compatible with the NEBNext Ultra II Directional RNA Workflow. If you need access to the previous version of the manual, please contact info@neb.com.

Starting Material: 20 μ l of first strand cDNA synthesized with the NEBNext Ultra II RNA First Strand RNA Synthesis Module (#E7771, Chapter 1).

1. Second Strand cDNA Synthesis

- 1.1. Assemble the second strand cDNA synthesis reaction **on ice** by adding the following components into the first strand synthesis reaction product.

COMPONENT	VOLUME
First-strand Synthesis Product	20 μ l
• (orange) NEBNext Second Strand Synthesis Reaction Buffer with dUTP Mix	8 μ l
• (orange) NEBNext Second Strand Synthesis Enzyme Mix	4 μ l
Nuclease-free Water	48 μ l
Total Volume	80 μ l

- 1.2. Keeping the tube on ice, mix thoroughly by pipetting the reaction up and down at least 10 times.
1.3. Incubate in a thermocycler for **1 hour at 16°C** with the heated lid set at $\leq 40^\circ\text{C}$ (or off).

2. Purification of Double-stranded cDNA using SPRIselect Beads or NEBNext Sample Purification Beads

- 2.1. Vortex SPRIselect Beads or NEBNext Sample Purification Beads to resuspend.
2.2. Add 144 μ l (1.8X) of resuspended beads to the second strand synthesis reaction (~80 μ l). Mix well on a vortex mixer or by pipetting up and down at least 10 times.
2.3. Incubate for 5 minutes at room temperature.
2.4. Briefly spin the tube in a microcentrifuge to collect any sample from the sides of the tube. Place the tube on a magnetic rack to separate beads from the supernatant. After the solution is clear, carefully remove and discard the supernatant. Be careful not to disturb the beads, which contain DNA. **(Caution: Do not discard beads)**
2.5. Add 200 μ l of freshly prepared 80% ethanol to the tube while in the magnetic stand. Incubate at room temperature for 30 seconds, and then carefully remove and discard the supernatant.
2.6. Repeat Step 2.5 once for a total of 2 washing steps.
2.7. Air dry the beads for up to 5 minutes while the tube is on the magnetic rack with the lid open.

Caution: Do not overdry the beads. This may result in lower recovery of DNA target. Elute the samples when the beads are still dark brown and glossy looking, but when all visible liquid has evaporated. When the beads turn lighter brown and start to crack they are too dry.

- 2.8. Remove the tube from the magnetic rack. Elute the DNA from the beads by adding 53 μ l 0.1X TE Buffer (provided) to the beads. Mix well on a vortex mixer or by pipetting up and down at least 10 times. Quickly spin the tube and incubate for 2 minutes at room temperature. Place the tube on the magnetic rack until the solution is clear.
2.9. Remove 50 μ l of the supernatant and transfer to a clean nuclease-free PCR tube.
2.10. Proceed to the NEBNext Ultra II End Repair/dA-Tailing Module (NEB #E7546)

Kit Components

Each set of reagents is functionally validated together with NEB #E7111, NEB #E7546, NEB #E7595, and NEB #M0544, and compared to the previous lot through construction of libraries using the minimum and maximum amount of Universal Human Reference Total RNA. The previous and current lots are sequenced together on the same Illumina flow cell and compared across various sequence metrics including individual transcript abundances, 5'→3' transcript coverage, and fraction of reads mapping to the reference

NEB #E7550S Table of Components

NEB #	PRODUCT	VOLUME
E7425A	NEBNext Second Strand Synthesis Enzyme Mix	0.096 ml
E7426A	NEBNext Second Strand Synthesis Reaction Buffer with dUTP Mix	0.192 ml

NEB #E7550L Table of Components

NEB #	PRODUCT	VOLUME
E7425AA	NEBNext Second Strand Synthesis Enzyme Mix	0.384 ml
E7426AA	NEBNext Second Strand Synthesis Reaction Buffer with dUTP Mix	0.768 ml

Revision History

REVISION #	DESCRIPTION	DATE
2.0	Change name of kit from NEBNext Ultra Directional RNA Second Strand Synthesis Module to NEBNext Ultra II Directional RNA Second Strand Synthesis Module. Protocol changed to be compatible with the NEBNext Ultra II Directional RNA Workflow.	4/17
3.0	Update "NEBNext Ultra II Directional RNA Second Strand Synthesis Module is Designed for Use with the Following:" list of products. Update protocol .	4/18
4.0	Update protocol.	6/18
5.0	New format applied.	1/20

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 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.

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

ILLUMINA® is a registered trademark of Illumina, Inc.

QIAQUICK® is a registered trademark of Qiagen.

© Copyright 2019, New England Biolabs, Inc.; all rights reserved



www.neb.com



be INSPIRED
drive DISCOVERY
stay GENUINE