

NEBNext® Enzymatic 5hmC-seq Conversion Module

NEB E3365S/L

24/96 reactions

Version 2.0_7/26

Table of Contents

Overview	2
Protocol for the E5hmC-seq Conversion Module.....	3
Checklist	9
Kit Components	11
Revision History	12

The Conversion Module Includes

The volumes provided are sufficient for preparation of up to 24 reactions (NEB #E3365S) or 96 reactions (NEB #E3365L).

All reagents should be stored at –20°C. Colored bullets represent the color of the cap of the tube containing the reagent.

- (lilac) Control DNA Unmethylated Lambda
- (lilac) Control DNA 5hmC T4
- (red) NEBNext Carrier DNA
- (yellow) NEBNext Glucosylation Reaction Buffer
- (yellow) UDP-Glucose
- (yellow) T4-BGT
- (yellow) Stop Reagent
- (white) Elution Buffer
- (orange) APOBEC
- (orange) Deamination Reaction Buffer
- (orange) Recombinant Albumin

Required Materials Not Included

- If DNA requires fragmentation: NEBNext UltraShear™ (NEB #M7634) or Covaris® instrument and the required tubes or other fragmentation equipment
- PCR strip tubes or 96-well plates
- Clean-up beads: SPRIselect™ Reagent Kit (Beckman Coulter®, Inc. #B23317), AMPure® XP beads (Beckman Coulter, Inc. #A63881) or preferred bead manufacturer.
- Formamide (Sigma #F9037-100 ml), Hi-Di™ Formamide (Thermo Fisher Scientific® #4401457) or 0.1 N NaOH. Formamide is preferred. If using NaOH, please see FAQ on NEB #E3365 FAQ page.
- 80% Ethanol
- 1X TE (10 mM Tris-HCl pH 8.0, 1 mM EDTA), low TE (10 mM Tris-HCl pH 8.0, 0.1 mM EDTA) or 10 mM Tris-HCl pH 7.5 or 8.0
- Nuclease-free Water
- Magnetic rack/stand, such as NEBNext Magnetic Separation Rack (NEB #S1515)
- Metal cooling block, such as Diversified Biotech® (#CHAM-1000)
- PCR machine
- Agilent® Bioanalyzer®, TapeStation® or other fragment analyzer and associated consumables

Overview

The NEBNext Enzymatic 5hmC-seq Conversion Module contains the reagents required to enzymatically modify and enable specific detection of 5-hydroxymethylcytosine (5hmC). The user-friendly workflow enables 5hmC identification from 0.1 – 200 ng of input DNA.

Each kit component must pass rigorous quality control standards, and for each new lot the entire set of reagents is functionally validated together by construction of indexed libraries that are sequenced on an Illumina® sequencing platform.

Where larger volumes, customized or bulk packaging are required, we encourage consultation with the Customized Solutions team at NEB. Please complete the NEB Custom Contact Form at www.neb.com/CustomContactForm to learn more.

Figure 1. Overview of the E5hmC-seq conversion method

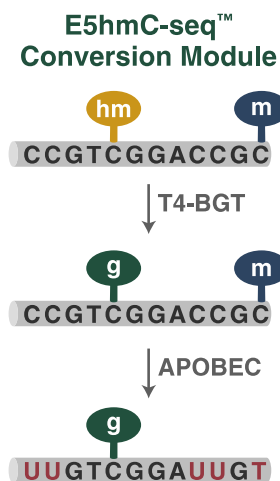


Figure 1 shows the two-step enzymatic conversion process used to detect 5hmC. The first step uses T4-BGT to glucosylate 5hmC. This protects 5hmC but not 5mC or cytosines from deamination by APOBEC. Note that the final DNA obtained from the NEBNext Enzymatic 5hmC-seq Conversion Module will contain U and glucosylated 5hmC.

Protocol for E5hmC-seq Conversion Module

Symbols



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.



This caution sign signifies a step in the protocol that has two paths leading to the same end point.



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

Important Guidelines

Reagent Handling

Remove only the kit reagents needed for a specific reaction in the EM-seq protocol. For example, for “Glucosylation of 5-Hydroxymethylcytosines (5hmC)” remove the yellow capped tubes. Frozen kit components can be thawed at room temperature. Once thawed, place immediately on ice.

Please note that some components may form a precipitate. Precipitates should disappear once the tubes are at room temperature. All components (stored at -20°C) should be stored on ice until used.

All components should be mixed before use. Typically, tubes containing enzymes can be mixed by flicking the tube and other buffers/reagents can be quickly vortexed. All tubes should be centrifuged briefly after mixing and stored on ice until used.

Master mixes can be made for most steps in E5hmC-seq unless otherwise stated in the protocol. If using a master mix, good practice is to add the component in the order specified in the protocol. Typically, enzymes are added last. Master mixes should be quickly vortexed then briefly centrifuged and stored on ice before use.

Care should be taken when handling some of the reagents to ensure the correct volumes are dispensed. For example, the Sample Purification Beads are viscous. It is difficult to accurately pipette these components. Please take care to transfer the correct volumes.

The ● (orange) Deamination Reaction Buffer can form bubbles when mixed. Centrifuging the ● (orange) Deamination Reaction Buffer can reduce their appearance making it easier to accurately dispense the buffer.

Once reactions are set up, mix either using a quick vortex or by pipetting up and down. Please note that if you choose to mix using a pipette, then we encourage you to use an appropriately sized pipette set to a volume that would ensure adequate mixing of the sample.

For example, during the Glucosylation Reaction set up, 5 µl of diluted ● (yellow) Fe(II) Solution is added, use a 200 µl pipette set to 40 µl to mix – do not mix with a pipette set to 5 µl. Briefly centrifuge the mixed reactions before incubating the reactions.

Another example where adequate mixing is required is during Sample Purification Bead clean-ups. Here, please use a 200 µl pipette set to an appropriate volume to ensure mixing. For the bead clean up after Glucosylation, a pipette set to 80 µl would be sufficient whereas for the clean-up after APOBEC deamination a 160 µl pipette volume would be more appropriate.

After each NEBNext Sample Purification Bead clean-up please ensure that the samples are centrifuged to ensure that the sample is at the bottom of the tube prior to the next reaction in the protocol. If the samples were stored at -20°C, they can be thawed at room temp or on ice prior to centrifugation.

There are safe stop points in the protocol where reactions can be stored at -20°C. If a sample has been stored at -20°C then it should be thawed and briefly centrifuged before use in the next step in the protocol.

Starting Material: 0.1–200 ng fragmented double stranded DNA

1.1. DNA Preparation

1.1.1. Sample DNA and Control DNA

DNA going into the first enzymatic reaction (protection of 5mC and 5hmC) reaction needs to be fragmented to a size appropriate for short read sequencing, regardless of downstream application.

For sequencing on an Illumina platform, refer to the Enzymatic 5hmC-seq Kit Manual (NEB #E3350) for usage recommendations. For other downstream applications and sequencing platforms, please refer to manufacturer’s guidelines.

The following table is a guide for the amount of ● (lilac) Control DNA Unmethylated Lambda and ● (lilac) Control DNA 5hmC T4 to be added to samples going into E5hmC-seq conversion. Expected read numbers along with read length should be considered to ensure that enough controls are included for the user’s individual sequencing goals or other downstream applications.

Table 1.1. Dilutions of control DNAs for a range of genomic DNA inputs.

Sample DNA Input Amount	Control DNA Dilution Recommendations
0.1 ng	1:1000
1 ng	1:250
10 ng	1:100
200 ng	1:10

For example, samples going into Illumina library preparation can be mechanically or enzymatically fragmented using NEBNext UltraShear (NEB #M7634) to an average size of ~350 bp. Control DNAs should be added prior to fragmentation to provide a means to qualify the final sequencing data as specified below. Samples can be fragmented in 1X TE (10 mM Tris-HCl pH 8.0, 1 mM EDTA), low TE (10 mM Tris-HCl pH 8.0, 0.1 mM EDTA), or 10 mM Tris-HCl pH 7.5 or 8.0. If the volume needs to be adjusted going into the Glucosylation Reaction, fragmented samples can be cleaned up using beads (~2X bead ratio) or column and eluted as specified below.

We do not recommend fragmenting input DNA in 0.1X TE (1 mM Tris-HCl pH 8.0, 0.1 mM EDTA) or water.

COMPONENT	VOLUME
Sample DNA	48 µl
• (lilac) Diluted Control DNA Unmethylated Lambda (see Table 1.1)	1 µl
• (lilac) Diluted Control DNA 5hmC T4 (see Table 1.1)	1 µl
Total Volume	50 µl

Note: While DNA and controls must be fragmented in a buffer containing 10 mM Tris-HCl (see above), the Glucosylation Reaction itself in Step 1.2. is compatible with samples in water, 1X TE, 0.1X TE, low TE, or 10 mM Tris pH 7.5 or 8.0.

Note: The volume going into conversion is 28 µl.

Regardless of sequencing depth, a minimum of 5,000 paired end reads with a read length of 76 bases for • (lilac) Control DNA Unmethylated Lambda and 15,000 paired end reads with a read length of 76 bases for • (lilac) Control DNA 5hmC T4 are needed to give enough coverage for accurate conversion estimates.

Different sequencing depths may be needed depending on the application, and therefore different strategies should be employed when deciding how much control DNA should be added. For example, some libraries may only need 2 million paired end reads whereas others may require 50 million paired end reads or even 1 billion paired end reads.

The dilutions recommended in Table 1.1. will provide sufficient coverage of controls for libraries sequenced to 10 million paired end reads and above. Dilution of controls needs to be optimized by the user if sequencing lower than 10 million paired-end reads to obtain minimum coverage for • (lilac) Control DNA Unmethylated Lambda (5,000 paired end reads) and for • (lilac) Control DNA 5hmC T4 (15,000 paired end reads). Users should be aware that deep sequencing using the dilutions recommended in Table 1.1. can result in more than the minimum required reads for the • (lilac) Control DNA Unmethylated Lambda and the • (lilac) Control DNA 5hmC T4. Ultimately, dilutions of the control DNAs should be optimized by the user.

Post fragmentation:

- **Inputs > 10 ng:** DNA must be in 28 µl of water, 1X TE, 0.1X TE, low TE, or 10 mM Tris pH 7.5 or 8.0 to go into the Glucosylation Reaction.
- **Inputs ≤ 10 ng:** For adaptor ligated DNA add 1 µl of • (red) NEBNext Carrier DNA to sample in 27 µl of water, 1X TE, 0.1X TE, low TE, or 10 mM Tris pH 7.5 or 8.0 to go into the Glucosylation Reaction.
- **Inputs ≤ 10 ng:** For DNA with no adaptors ligated, DNA must be in 28 µl of water, 1X TE, 0.1X TE, low TE, or 10 mM Tris pH 7.5 or 8.0 to go into the Glucosylation Reaction.

Note: For companion EM-seq product (NEB #E7120/E7125), there should be no EDTA in the sample going into conversion.

Note: Addition of the NEBNext Carrier DNA may not be appropriate for every application. For example, if sequencing, do not use with DNA that does not have adaptors ligated before conversion.

1.2. Glucosylation of 5-Hydroxymethylcytosines (5hmC)

- 1.2.1. On ice, add the following components directly to the fragmented DNA:

COMPONENT	VOLUME
Fragmented DNA (from Step 1.1.1.)	28 µl
Nuclease-free water	15 µl
● (yellow) NEBNext Glucosylation Reaction Buffer	5 µl
● (yellow) UDP-Glucose	1 µl
● (yellow) T4-BGT	1 µl
Total Volume	50 µl

Mix thoroughly by vortexing 1–2 seconds or by pipetting up and down at least 10 times and centrifuge briefly. For multiple reactions, a master mix of the reaction components can be prepared before addition to the sample DNA.

- 1.2.2. Place in a thermal cycler, and run the following program with the heated lid set to $\geq 45^{\circ}\text{C}$ or on:
1 hour at 37°C
Hold at 4°C

- 1.2.3. Transfer the samples to ice and add 1 µl of ● (yellow) Stop Reagent.

COMPONENT	VOLUME
Glucosylated DNA (from Step 1.2.1.)	50 µl
● (yellow) Stop Reagent	1 µl
Total Volume	51 µl

Mix thoroughly by vortexing 1–2 seconds or by pipetting up and down at least 10 times and centrifuge briefly.

- 1.2.4. Place in a thermal cycler, and run the following program with the heated lid set to $\geq 45^{\circ}\text{C}$ or on:
30 minutes at 37°C
Hold at 4°C .



Safe Stopping Point: Samples can be stored overnight at either 4°C in the thermal cycler or at -20°C in the freezer.

1.3. Clean-up of Glucosylated DNA

The ratios recommended for SPRIselect beads in this manual have been experimentally optimized for every step; this is critical since buffer compositions differ between steps and across protocols e.g., post ligation recommendations will not apply to samples post PCR. Please adhere to these guidelines and not those recommended by other sources or for other kits.

- 1.3.1. Vortex SPRIselect beads to resuspend.
- 1.3.2. Add 50 µl (1X ratio) of resuspended SPRIselect beads to each sample. Mix well by pipetting up and down at least 10 times. Be careful to expel all of the liquid out of the tip during the last mix.
- 1.3.3. Incubate samples on bench top for at least 5 minutes at room temperature.
- 1.3.4. Place the tubes against an appropriate magnetic stand to separate the beads from the supernatant.
- 1.3.5. After 5 minutes (or when the solution is clear), carefully remove and discard the supernatant. Be careful not to disturb the beads that contain DNA targets (**Caution: do not discard the beads**).
- 1.3.6. Add 200 µl of freshly prepared 80% ethanol to the tubes while on the magnetic stand. Incubate at room temperature for 30 seconds, and then carefully remove and discard the supernatant. Be careful not to disturb the beads that contain DNA targets.
- 1.3.7. Repeat the ethanol wash once for a total of two washes. Be sure to remove all visible liquid after the second wash using a p10 pipette tip.
- 1.3.8. Air-dry the beads for 30 seconds – 1 minute while the tubes are on the magnetic stand with the lid open.

Caution: Do not over-dry the beads. This may result in lower recovery of DNA targets. 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.

- 1.3.9. Remove the tubes from the magnetic stand. Elute the DNA targets from the beads by adding 17 μ l of $\circ$ (white) Elution Buffer.
- 1.3.10. Mix well by pipetting up and down 10 times. Incubate for at least 1 minute at room temperature. If necessary, quickly spin the sample to collect the liquid from the sides of the tube before placing back on the magnetic stand.
- 1.3.11. Place the tube on the magnetic stand. After 3 minutes (or whenever the solution is clear), transfer 16 μ l of the supernatant to a new PCR tube.

Caution: Carrying even a small amount of beads forward can lead to inefficient deamination.



Safe Stopping Point: Samples can be stored overnight at -20°C .

1.4. Denaturation of DNA



Denaturation Options A or B

The DNA can be denatured using either Formamide or 0.1 N Sodium Hydroxide. Use Option A for denaturing with Formamide and Option B for denaturing with 0.1 N Sodium Hydroxide.

Option 1.4A: Formamide (Recommended)

- 1.4A.1. Pre-heat thermal cycler to 85°C with the heated lid set to $\geq 105^{\circ}\text{C}$ or on.
- 1.4A.2. Add 4 μ l Formamide to the 16 μ l of glucosylated DNA (from Step 1.3.11.). Mix thoroughly by vortexing 1–2 seconds or by pipetting up and down at least 10 times and centrifuge briefly.
- 1.4A.3. Incubate at 85°C for 10 minutes in the pre-heated thermal cycler.
- 1.4A.4. Immediately place in a pre-chilled metal cooling block on ice and allow the sample to fully cool (~ 2 minutes) before proceeding to Section 1.5. If a metal cooling block is not available, place on ice and ensure proper contact with the tube walls.

Option 1.4B: Sodium Hydroxide

Optional, See FAQ about preparing NaOH.

- 1.4B.1. Prepare freshly diluted 0.1 N NaOH.
- 1.4B.2. Pre-heat thermal cycler to 85°C with the heated lid set to $\geq 105^{\circ}\text{C}$ or on.
- 1.4B.3. Add 4 μ l 0.1 N NaOH to the 16 μ l of glucosylated DNA (from Step 1.3.11.). Mix thoroughly by vortexing 1–2 seconds or by pipetting up and down at least 10 times and centrifuge briefly.
- 1.4B.4. Incubate at 85°C for 10 minutes in the pre-heated thermal cycler.
- 1.4B.5. Immediately place in a pre-chilled metal cooling block on ice and allow the sample to fully cool (~ 2 minutes) before proceeding to Section 1.5. If a metal cooling block is not available, place on ice and ensure proper contact with the tube walls.

1.5. Deamination of 5mC and Cytosines



Deamination Options A or B

Use Option A (1.5.3A.) for DNA inputs ≤ 100 ng

Option B (1.5.3B.) is required for > 100 ng DNA inputs and can be used for all inputs (0.1 ng to 200 ng).

- 1.5.1. On ice, add the following components to the denatured DNA:

COMPONENT	VOLUME
Denatured DNA (Step 1.4A.4. or 1.4B.5.)	20 µl
Nuclease-free water	14 µl
• (orange) Deamination Reaction Buffer	4 µl
• (orange) Recombinant Albumin	1 µl
• (orange) APOBEC	1 µl
Total Volume	40 µl

For multiple reactions, a master mix of the reaction components can be prepared before addition to the denatured DNA.

- 1.5.2. Mix thoroughly by vortexing 1–2 seconds or by pipetting up and down at least 10 times and centrifuge briefly.

Option 1.5.3A: (Inputs ≤ 100ng)

- 1.5.3A. Place in a thermal cycler, and run the following program with the heated lid set to ≥ 85°C or on:

10 min at 37°C
10 min at 42°C
10 min at 50°C
2 min at 75°C
Hold at 4°C

Option 1.5.3B: (Required for Inputs > 100 ng and can be used for all inputs)

- 1.5.3B. Place in a thermal cycler, and run the following program with the heated lid set to ≥ 45°C or on:

3 hours at 37°C
Hold at 4°C



Safe Stopping Point: Samples can be stored overnight at either 4°C in the thermal cycler or at –20°C in the freezer.

1.6. Clean-up of Deaminated DNA

The ratios recommended for SPRIselect beads in this manual have been experimentally optimized for every step; this is critical since buffer compositions differ between steps and across protocols e.g., post ligation recommendations will not apply to samples post PCR. Please adhere to these guidelines and not those recommended by other sources or for other kits.

Caution: The Sample Purification Beads behave differently during the APOBEC clean-up. After the bead washes, do not overdry the beads as they become very difficult to resuspend.

- 1.6.1. Vortex SPRIselect beads to resuspend.
- 1.6.2. Add 40 µl (1X ratio) of resuspended SPRIselect beads to each sample. Mix well by pipetting up and down at least 10 times. Be careful to expel all of the liquid out of the tip during the last mix.
- 1.6.3. Incubate samples on bench top for at least 5 minutes at room temperature.
- 1.6.4. Place the tubes against an appropriate magnetic stand to separate the beads from the supernatant.
- 1.6.5. After 5 minutes (or when the solution is clear), carefully remove and discard the supernatant. Be careful not to disturb the beads that contain DNA targets (**Caution: do not discard the beads**).
- 1.6.6. Add 200 µl of freshly prepared 80% ethanol to the tubes while on the magnetic stand. Incubate at room temperature for 30 seconds, and then carefully remove and discard the supernatant. Be careful not to disturb the beads that contain DNA targets.
- 1.6.7. Repeat the ethanol wash once for a total of two washes. Be sure to remove all visible liquid after the second wash using a p10 pipette tip.
- 1.6.8. Air-dry the beads for 30 seconds – 1 minute while the tubes are on the magnetic stand with the lid open.

Caution: Do not over-dry the beads. This may result in lower recovery of DNA targets. 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.

- 1.6.9. Remove the tubes from the magnetic stand. Elute the DNA targets from the beads by adding 21 μ l of $\circ$ (white) Elution Buffer or an appropriate buffer for downstream application.
- 1.6.10. Mix well by pipetting up and down 10 times. Incubate for at least 1 minute at room temperature. If necessary, quickly spin the sample to collect the liquid from the sides of the tube before placing back on the magnetic stand.
- 1.6.11. Place the tube on the magnetic stand. After 3 minutes (or whenever the solution is clear), transfer 20 μ l of the supernatant to a new PCR tube.

Checklist

1.1 DNA Preparation

1.1.1. Combine DNA and control DNA

See Table 1.1 for recommended control dilutions

- ☐ Fragment 0.1 ng – 200 ng sample DNA with ● (lilac) Control DNA Unmethylated Lambda and ● (lilac) Control DNA 5hmC T4
- ☐ 1 µl of ● (red) NEBNext Carrier DNA for inputs ≤ 10 ng (**application dependent; do not include if adaptors have not been ligated**)

1.2. Glucosylation of 5-Hydroxymethylcytosines

Add Glucosylation Reagents to 28 µl DNA

- ☐ 15 µl Nuclease-free water
- ☐ 5 µl ● (yellow) NEBNext Glucosylation Reaction Buffer
- ☐ 1 µl ● (yellow) UDP-Glucose
- ☐ 1 µl ● (yellow) T4-BGT
- ☐ Vortex 1–2 seconds or pipette mix 10 times, centrifuge briefly
- ☐ Incubate in a thermal cycler (heated lid ≥ 45°C or on)
60 minutes at 37°C
Hold at 4°C
- ☐ Add 1 µl ● (yellow) Stop Reagent
- ☐ Vortex 1–2 seconds or pipette mix 10 times, centrifuge briefly
- ☐ Incubate in a thermal cycler (heated lid ≥ 45°C or on)
30 minutes at 37°C
Hold at 4°C

1.3. Clean-up of Glucosylated DNA

- ☐ Vortex SPRIselect beads
- ☐ Add 50 µl of resuspended beads to each sample and mix by pipetting 10 times
- ☐ Incubate 5 minutes
- ☐ Place tubes on magnet for 5 minutes
- ☐ Remove and discard the supernatant, while keeping the sample on the magnet
- ☐ On magnet add 200 µl 80% ethanol, wait 30 seconds and then remove and discard the ethanol wash
- ☐ Repeat the ethanol wash
- ☐ Air-dry the beads for 30 seconds – 1 minute while on magnet
- ☐ Remove the samples from the magnet and resuspend in 17 µl of ○ (white) Elution Buffer
- ☐ Place back on the magnet, wait until the supernatant clears and transfer 16 µl of sample to fresh PCR tube

1.4. Denaturation of DNA

Use either Formamide (A) or Sodium Hydroxide (B)

A. Formamide Denaturation (Recommended)

- ☐ Pre-heat thermal cycler to 85°C (heated lid $\geq$ 105°C or on)
- ☐ Add 4 μ l Formamide to the 16 μ l glucosylated DNA
- ☐ Vortex 1 – 2 seconds or pipette mix 10 times, centrifuge briefly
- ☐ Incubate in pre-heated thermal cycler
85°C for 10 minutes
- Immediately, place in cooling block on ice until fully cooled (~2 minutes)
- ☐ Proceed immediately into Section 1.5.

B. Sodium Hydroxide Denaturation (Formamide is preferred)

- ☐ Prepare freshly diluted 0.1 N NaOH
- ☐ Pre-heat thermal cycler to 85°C (heated lid $\geq$ 105°C or on)
- ☐ Add 4 μ l 0.1 N NaOH to the 16 μ l glucosylated DNA
- ☐ Vortex 1–2 seconds or pipette mix 10 times, centrifuge briefly
- ☐ Incubate in pre-heated thermal cycler
85°C for 10 minutes
- Immediately, place in cooling block on ice until fully cooled (~2 minutes)
- ☐ Proceed immediately into Section 1.5.

1.5. Deamination of 5mC and Cytosines

Add Deamination Reagents to 20 μ l denatured DNA on ice

- ☐ 14 μ l Nuclease-free water
- ☐ 4 μ l ● (orange) Deamination Reaction Buffer
- ☐ 1 μ l ● (orange) Recombinant Albumin
- ☐ 1 μ l ● (orange) APOBEC
- ☐ Vortex 1–2 seconds or pipette mix 10 times, centrifuge briefly
- ☐ Incubate in a thermal cycler (heated lid $\geq$ 45°C or on)
3 hours at 37°C
Hold at 4°C

1.6. Clean-up of Deaminated DNA

- ☐ Vortex SPRIselect beads
- ☐ Add 40 μ l of resuspended beads to each sample and mix by pipetting 10 times
- ☐ Incubate 5 min
- ☐ Place tubes on magnet for 5 minutes
- ☐ Remove and discard the supernatant, while keeping the sample on the magnet
- ☐ On magnet add 200 μ l 80% ethanol, wait 30 seconds and then remove and discard the ethanol wash
- ☐ Repeat the ethanol wash
- ☐ Air-dry the beads for 30 seconds – 1 minute while on magnet
- ☐ Remove the samples from the magnet and resuspend in 21 μ l of ○ (white) Elution Buffer or appropriate buffer for downstream application.
- ☐ Place back on the magnet, wait until the supernatant clears and transfer 20 μ l of sample to fresh PCR tube

Kit Components

NEB #E3365S Table of Components

NEB #	PRODUCT	VOLUME
E7123A	Control DNA Unmethylated Lambda	0.024 ml
E3349A	Control DNA 5hmC T4	0.024 ml
E3351A	NEBNext Carrier DNA	0.024 ml
E3352A	NEBNext Glucosylation Reaction Buffer	0.12 ml
E3353A	UDP-Glucose	0.024 ml
E3354A	T4-BGT	0.024 ml
E7132A	Stop Reagent	0.024 ml
E7124A	Elution Buffer	2.1 ml
E7133A	APOBEC	0.024 ml
E3356A	Deamination Reaction Buffer	0.096 ml
E3357A	Recombinant Albumin	0.024 ml

NEB #E3365L Table of Components

NEB #	PRODUCT	VOLUME
E7123AA	Control DNA Unmethylated Lambda	0.096 ml
E3349AA	Control DNA 5hmC T4	0.096 ml
E3351AA	NEBNext Carrier DNA	0.096 ml
E3352AA	NEBNext Glucosylation Reaction Buffer	0.480 ml
E3353AA	UDP-Glucose	0.096 ml
E3354AA	T4-BGT	0.096 ml
E7132AA	Stop Reagent	0.096 ml
E7124AA	Elution Buffer	8.6 ml
E7133AA	APOBEC	0.096 ml
E3356AA	Deamination Reaction buffer	0.384 ml
E3357AA	Recombinant Albumin	0.096 ml

Revision History

REVISION #	DESCRIPTION	DATE
1.0	N/A	12/23
1.1	Clarified post fragmentation verbiage in Section 1.1. Updated header and footer.	2/24
2.0	Update protocol and add important guidelines.	7/26

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.

Products and content are covered by one or more patents, trademarks and/or copyrights owned or controlled by New England Biolabs, Inc (NEB). The use of trademark symbols does not necessarily indicate that the name is trademarked in the country where it is being read; it indicates where the content was originally developed. See www.neb.com/trademarks. The use of these products may require you to obtain additional third-party intellectual property rights for certain applications. For more information, please email busdev@neb.com.

B CORPORATION® is a registered trademark of B Lab IP, LLC, Inc.

AGILENT®, BIOANALYZER® and TAPESTATION® are registered trademarks of Agilent Technologies, Inc.

BECKMAN COULTER® and AMPURE® are registered trademarks of Beckman Coulter, Inc. SPRISELECT™ is a trademark of Beckman Coulter, Inc.

COVARIS® is a registered trademark of Covaris, LLC

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

DIVERSIFIED BIOTECH® is a registered trademark of Diversified Biotech, Inc.

HI-DI™ is a trademark of Thermo Fisher Scientific.

THERMO FISHER SCIENTIFIC® is a registered trademark of Thermo Fisher Scientific, Inc.

© Copyright 2026, New England Biolabs, Inc.; all rights reserved



New England Biolabs, Inc., 240 County Road, Ipswich, MA 01938-2723 | T: (978) 927-5054 F: (978) 921-1350 US Orders (Toll-free): 1-800-632-5227 US Tech Support: 1-800-632-7799 | info@neb.com