

# Whole-exosome sequencing - Purification of amplified post capture libraries for Colibri™ DNA Library Prep Kits for Illumina® systems

**Catalog Numbers** A38614196, A38612024, A38614096, A38613024, A43611024, A43612024, A43613024, A38607096, A38605024, A38607196, A38606024, A43605024, A43606024, A43607024, A38605024W, A38607096W, A38612024W, A38614096W

Pub. No. MAN0026045 Rev. A.0

**Note:** For safety and biohazard guidelines, see the “Safety” appendix in the Colibri™ DNA Library Prep Kit for Illumina® User Guides (Pub. Nos. MAN0025533, MAN0025534). Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

## About the protocol

This section describes bead-based purification of amplified post-capture DNA libraries for whole-exosome sequencing (WES) samples.

## Required materials

Components from the Colibri™ ES DNA Library Prep Kits for Illumina®
DNA Cleanup Beads
Wash Buffer (diluted with 96% ethanol)
Elution Buffer
Other consumables
Amplified post-capture DNA libraries
96% ethanol, molecular biology grade (used for diluting the Wash Buffer before first use)
Eppendorf™ 1.5-mL DNA LoBind microcentrifuge tubes
Microcentrifuge
Magnetic rack

## Before starting

- Ensure that the appropriate volume of 96% ethanol (as noted on the bottle) was added to the Wash Buffer before first use.
- Ensure that the DNA Cleanup Beads, Wash Buffer, and Elution Buffer are at room temperature.
- Gently vortex the DNA Cleanup Beads to completely resuspend the magnetic beads in the solution.

## Purify the amplified post-capture DNA library

Perform all the cleanup steps at room temperature.

### Bind the library

1. Mix amplified post-capture library (50  $\mu$ L) with 50  $\mu$ L of DNA Cleanup Beads by vortexing until you have obtained a homogeneous suspension.

**Note:** It is **not** necessary to recover the supernatant or remove the Target Enrichment and Amplification Beads from the amplified PCR product.

2. Briefly centrifuge the tube containing the amplified DNA library and bead mixture to collect all the droplets at the bottom, then incubate for 5 minutes at room temperature.

---

**IMPORTANT! Do not** extend the binding step for more than 5 minutes. Over-incubation can result in lower DNA yields.

---

3. Briefly centrifuge the tube to collect all the droplets at the bottom, then place the tube in a magnetic rack for 2 minutes or until the beads have formed a tight pellet.

**Note:** The time required for complete capture of the cleanup beads can vary depending on the reaction vessel and the magnet used. Optimize the capture time accordingly.

4. Keep the reaction tube in the magnetic rack, then carefully remove and discard the supernatant using a pipette. Ensure that all the supernatant is removed.

**Note:** If the bead pellet is disturbed, mix the sample and allow the beads to form a tight pellet on the magnet again.

### Wash the magnetic bead pellet

1. Keep the reaction tube in the magnetic rack and add 200  $\mu$ L of Wash Buffer (pre-mixed with ethanol).

---

**IMPORTANT! Do not** resuspend the magnetic beads in Wash Buffer.

---

2. Incubate for 30 seconds at room temperature, then carefully remove and discard the supernatant using a pipette.
3. Repeat steps 1–2.
4. Briefly centrifuge the reaction tube and place it in a magnetic rack. Without disturbing the pellet, carefully remove any remaining supernatant and residual ethanol.
5. Keep the reaction tube in the magnetic rack and air dry the magnetic beads for 1 minute at room temperature, or until there are no droplets of ethanol left on the walls of the tube.

---

**IMPORTANT! Do not** dry the magnetic beads for more than 5 minutes. Over-drying significantly decreases the elution efficiency.

---

### Elute the library

1. Remove the tube from the magnetic rack and add 32  $\mu$ L of Elution Buffer, then mix the suspension thoroughly by pipetting up and down or vortexing.
2. Briefly centrifuge the tube to collect all the droplets at the bottom, then incubate for 1 minute at room temperature.
3. Place the tube in a magnetic rack for 2 minutes or until the beads have formed a tight pellet. Wait for the solution to clear before proceeding to the next step..

4. Keep the tube in the magnetic rack, and transfer 30  $\mu$ L of the supernatant (i.e., the eluate) to a new tube for storage.

**Note:** If the bead pellet is disturbed, mix the sample and allow the beads to form a tight pellet on the magnet again.


**Note:** Use Nuclease-free 1.5-mL tubes, such as Eppendorf™ 1.5-mL DNA LoBind microcentrifuge tubes (or equivalent) for long term storage.

5. 14. Proceed to the assessment of the DNA library size and yield.

6. After purification, proceed immediately to sequencing, or store the enriched libraries at  $-20^{\circ}\text{C}$ .

# Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale at [www.thermofisher.com/us/en/home/global/terms-and-conditions.html](http://www.thermofisher.com/us/en/home/global/terms-and-conditions.html). If you have any questions, please contact Life Technologies at [www.thermofisher.com/support](http://www.thermofisher.com/support).

 Thermo Fisher Scientific Baltics UAB | V.A. Graiciuno 8, LT-02241 | Vilnius, Lithuania  
For descriptions of symbols on product labels or product documents, go to [thermofisher.com/symbols-definition](http://thermofisher.com/symbols-definition).

The information in this guide is subject to change without notice.

**DISCLAIMER:** TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

**Revision history:** Pub. No. MAN0026045

Revision	Date	Description
A.0	14 January 2022	Replacement post-capture purification protocol for Colibri DNA Library Prep Kits for Whole-Exome Sequencing.

**Important Licensing Information:** These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2022 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.