

Automated Dynabeads™-Based Isolation of Human Epithelial Cells using the KingFisher™ Apex Purification System

Catalog Numbers 16102, 16103D

Pub. No. MAN1000760 Rev. A



WARNING! Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. Safety Data Sheets (SDSs) are available from [thermofisher.com/support](https://www.thermofisher.com/support).

Product description

The Invitrogen™ Dynabeads™ Epithelial Enrich are superparamagnetic beads coupled to a monoclonal mouse anti-human EpCAM antibody. The beads are designed to isolate epithelial circulating tumor cells (CTCs) from liquid biopsy samples. The isolation process is simple. The sample is mixed with the beads for targeted cell capture. After a short incubation and subsequent washing, the bead-bound cells are separated by the magnetic head of a KingFisher™ instrument. The bead-bound cells are then ready for further downstream analysis, including lysis and processing for protein or nucleic acid analysis.

This guide describes automated magnetic bead-based isolation of human epithelial cells from peripheral blood mononuclear cells (PBMCs) or whole blood using the KingFisher™ Apex Purification System.

The automated protocol is completed in 45 minutes, allowing up to 96 samples to be processed simultaneously in a single run.

Contents and storage

Table 1 Dynabeads™ Epithelial Enrich

| Item | Cat. No. | Amount | Total PBMC capacity | Whole blood capacity | Storage |
|------------------------------|----------|--------|---------------------|----------------------|---------|
| Dynabeads™ Epithelial Enrich | 16103D | 2 mL | 8×10^8 | 80 mL | 2–8°C |
| | 16102 | 5 mL | 2×10^9 | 200 mL | |

Required materials not supplied

Unless otherwise indicated, all materials are available through [thermofisher.com](https://www.thermofisher.com). "MLS" indicates that the material is available from [fisherscientific.com](https://www.fisherscientific.com) or another major laboratory supplier.

Catalog numbers that appear as links open the web pages for those products.

IMPORTANT! Use only barcoded plates and tip combs with the KingFisher™ Apex instrument. Non-barcoded deep-well plates or tip combs will increase processing times, may damage the system, void the warranty, and reduce performance.

| Item | Source |
|--|-------------------------|
| Instrument | |
| KingFisher™ Apex with 96 Deep-Well Head | 5400930 |
| Equipment | |
| Laboratory mixer, vortex, or equivalent | MLS |
| Single and multichannel adjustable pipettors (1 µL–1 mL) | MLS |

| Item | Source |
|---|---------------------------|
| Plates, combs, plate seals, and pipetting consumables | |
| KingFisher™ 96 Tip Comb for Deep-Well Magnets, Barcoded | 97002534B |
| KingFisher™ 96 Deep-Well Plates, Barcoded | 95040450B |
| MicroAmp™ Adhesive Film Applicator | 4333183 |
| MicroAmp™ Clear Adhesive Film | 4306311 |
| Pipette tips (10 µL–1 mL), sterile | MLS |
| Reagents | |
| Isolation Buffer: Calcium and magnesium-free phosphate buffered saline (PBS) with 0.1% BSA and 2 mM EDTA, pH 7.4 ^[1] | MLS |

^[1] BSA can be replaced by human serum albumin (HSA) or fetal calf serum (FCS). EDTA can be replaced by sodium citrate.

Set up the instrument

Ensure that the instrument is set up for processing with the magnetic head (96 deep-well) and heat block (96 deep-well).

1. On the product web page (at thermofisher.com, search by catalog number), scroll to the **Documents & Downloads** section.
2. Locate and download the appropriate script file for your instrument and install it onto the instrument.
Note: The script serves as proof-of-concept and can be modified to meet specific user needs.
3. Refer to your KingFisher™ instrument user guide or contact Technical Support at <http://thermofisher.com/support> for detailed instructions on script installation.

Prepare magnetic beads

Vortex the magnetic beads in the vial for 30 seconds, then rotate the vial for 5 minutes.

Note: Vortex and rotate the beads immediately before use to ensure optimal performance. Ensure that no beads remain adhered to the vial.

Prepare samples

Target cells can be directly captured from PBMC suspensions.

Note: Keep anti-coagulated whole blood containing EDTA or ACD on ice.

Set up the processing plates

1. Label the side of five deep-well plates according to the Plate IDs in Table 2.
2. Set up the processing plates according to the following table.

Table 2 Processing plate set-up

| Plate ID | Plate loading position ^[1] | Reagents | Volume per well | Plate type |
|----------|---------------------------------------|--|-----------------|-----------------|
| Tip | 1 | Place a Tip Comb into an empty deep-well plate | | Deep-well plate |
| Beads | 2 | Isolation Buffer Dynabeads™ Epithelial Enrich | 175 µL 25 µL | |
| Wash | 3 | Isolation Buffer | 500 µL | |
| Sample | 4 | PBMC suspension | 1000 µL | |
| Elution | 5 | Isolation Buffer | 100 µL | |

^[1] Position on the instrument.

3. Proceed to “Process samples on the instrument”.

Process samples on the instrument

1. Select the appropriate script on the instrument.
2. Start the run, then load the prepared processing plates in their positions when prompted by the instrument.

Note: Once all five plates are loaded, the instrument will start processing.

IMPORTANT! When loading plates, ensure that the A1 position of the plate aligns with the A1 position indicated on the instrument turntable.

3. At the end of the run, remove the Elution Plate from the instrument. If the samples are not processed immediately, cover the plate with MicroAmp™ Clear Adhesive Film.

For downstream analysis, the beads can be resuspended directly in the desired lysis buffer instead of isolation buffer. Optimize the volume of lysis buffer as needed.

Related documentation

Table 3 Product-specific user manual

| Document | Publication number | Cat. No. |
|---|----------------------------|--|
| Dynabeads™ Epithelial Enrich User Guide | MAN1000855 | 16102 , 16103D |

Customer and technical support

Visit [thermofisher.com/support](https://www.thermofisher.com/support) for the latest service and support information.

- Worldwide contact telephone numbers
- Product support information
 - Product FAQs
 - Software, patches, and updates
 - Training for many applications and instruments

- Order and web support
- Product documentation
 - User guides, manuals, and protocols
 - Certificates of Analysis
 - Safety Data Sheets (SDSs; also known as MSDSs)

Note: For SDSs for reagents and chemicals from other manufacturers, contact the manufacturer.

Limited product warranty

Life Technologies Corporation and its affiliates 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. If you have questions, contact Life Technologies at 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.

Revision history: Pub. No. MAN1000759 A

| Revision | Date | Description |
|----------|----------------|--|
| A | 3 January 2025 | New document for Automated Dynabeads-Based Isolation of Human Epithelial Cells using the KingFisher™ Apex Purification System. |

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.

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.

©2025 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.