

Version: 01

Update: 09/27/2022

CytoSinct™ Streptavidin Nanobeads

Contents

| | |
|--|---|
| I. <u>Product Description</u> | 1 |
| II. <u>Product Specification</u> | 1 |
| III. <u>Requirement Materials</u> | 2 |
| IV. <u>Protocol</u> | 2 |

I. Product Description

The CytoSinct™ Streptavidin Nanobeads are used for cell separation from fresh or frozen peripheral blood mononuclear cells (PBMCs), leukapheresis products or single cell suspension in combination with biotinylated antibody or ligand. The surface of Nanobeads are labeled with streptavidin protein. To begin the isolation, the cells are stained with a biotinylated primary antibody or ligand. Subsequently, Nanobeads are added to the cells. The streptavidin preconjugated to the Nanobeads can bind the biotinylated ligand on cell surface. The cells/beads mixed suspension is loaded onto a CytoSinct™ Column which is effectively magnetized by an external magnetic field from CytoSinct™ Magnet, or other compatible cell isolation columns and magnets. The Nanobeads-labeled cells are retained within the column and enriched during the wash step when Isolation Buffer is used to flush out the cells. After removing magnetic field, the target cells can be easily eluted from the column.

II. Product Specification

| Cat. No. | Name | Size | Capacity |
|------------|-----------------------------------|--------|---|
| L00898-0.5 | CytoSinct™ Streptavidin Nanobeads | 500 µL | for up to 5X10 ⁸ total cells |
| L00898-1 | CytoSinct™ Streptavidin Nanobeads | 1 mL | for up to 1X10 ⁹ total cells |

Product format Bio-degradable matrix coated nanoparticle conjugated with streptavidin supplied in water.

Application Positive selection or depletion of target cells from leukapheresis, PBMC or cell cultures in combination with biotinylated ligand.

Storage Store at 2-8 °C. Do not freeze.

III. Requirement Materials

1. **Isolation Buffer:** Prepare a solution containing phosphate-buffered saline (PBS), pH 7.2, 0.5% bovine serum albumin (BSA) and 2 mM EDTA.

- Keep Isolation Buffer cold (2–8°C).
- BSA can be replaced by human serum albumin (HSA), human serum or fetal bovine serum (FBS).
- EDTA can be replaced by sodium citrate.
- PBS containing Ca²⁺ or Mg²⁺ is not recommended.

2. **Columns and separators:**

- For samples containing less than 2×10⁸ total MNCs or less than 10⁷ labeled cells, use CytoSinct™ gM Column and CytoSinct™ M1 Magnet or CytoSinct™ M8 Magnet, or other compatible columns and magnets.
- For samples containing less than 2×10⁹ total cells or less than 10⁸ labeled cells, use CytoSinct™ gL Column and CytoSinct™ L1 Magnet or CytoSinct™ L4 Magnet or other compatible columns and magnets.

IV. Protocol

All procedures are to be performed at room temperature unless otherwise instructed in this protocol.

1. **Prepare Nanobeads**

Gently mix the Nanobeads by pipetting for several times.

2. **Prepare samples**

2.1 Prepare PBMCs.

- When working with anticoagulated peripheral blood, peripheral blood mononuclear cells (PBMCs) should be isolated by density gradient centrifugation method (for example, using Ficoll-Paque™ PLUS density gradient media) and washed by Isolation Buffer to remove interfering factors.
- When working with frozen PBMCs, resuscitate frozen PBMCs and then proceed with the protocol. When dead cells are found to be considerable, apply density gradient centrifugation method (for example, using Ficoll-Paque™ PLUS density gradient media) to remove dead cells, or culture cells in medium overnight before proceeding with this protocol.

2.2 Centrifuge PBMCs at 300 × g for 10 minutes at room temperature (15 - 25°C). Aspirate the supernatant completely. Determine cell number by using a hemocytometer or other suitable methods.

3. **Magnetic labeling**

3.1 Transfer desired number of cells into a new tube and resuspend into single cell suspension at

10^8 mononuclear cells (MNCs) per 1 mL in Isolation Buffer.

- When working with less than 10^7 MNCs, use Isolation Buffer volume of 100 μ L.

3.2 Because of different biotinylation process, proper amount of biotinylated antibody should be added at the titer recommended by the manufacture. For optimal cell labeling, it is recommended to test different antibody amount. Usually, for labeling 10^7 total MNCs, 0.15 to 5 μ g antibody can be added. Typically, labeling at 4°C for 15 minutes is sufficient. Mix the antibody and cells well by gently pipetting or tapping on the bottom of the tube.

3.3 Wash cells once by adding 1-2 mL of Isolation Buffer per 10^7 MNCs, mix well by gentle pipetting, and centrifuge at $300 \times g$ for 10 minutes. Aspirate supernatant completely.

3.4 Add 10 μ L Nanobeads for each 100 μ L cell suspension of 10^7 total MNCs.

- When working with higher number of cells, scale up the volume of Nanobeads accordingly (e.g. for 2×10^7 total MNCs, use 20 μ L Nanobeads).

- When working with less than 10^7 MNCs, use the same Nanobeads volume of 10 μ L, as that in 10^7 MNCs.

3.5 Mix the Nanobeads and cells well by gently pipetting or tapping on the bottom of the tube, and incubate for 15 min at 2–8 °C.

3.6 Wash cells once by adding 1-2 mL of Isolation Buffer per 10^7 MNCs, mix well by gentle pipetting, and centrifuge at $300 \times g$ for 10 minutes. Aspirate supernatant completely.

3.7 Resuspend up to 10^8 cells in 500 μ L of Isolation Buffer.

- Scale up the volume of Isolation Buffer accordingly when more than 10^8 MNCs are to be processed.

4. Magnetic separation

4.1 Choose an appropriate CytoSinct™ Column and CytoSinct™ Magnet or other compatible columns and magnet according to the number of total cells and the number of target cells as instructed in Section III.

4.2 Assemble the column onto the suitable CytoSinct™ Magnet or other compatible magnet (please refer to manual of CytoSinct™ Magnet or other compatible magnets for assembly instructions).

4.3 Rinse the column once with Isolation Buffer (500 μ L for CytoSinct™ gM column, 3 mL for CytoSinct™ gL Column, or other compatible columns) and let the buffer run through it but not run dry.

4.4 Transfer the cell suspension onto the prepared CytoSinct™ gM or CytoSinct™ gL Column or other compatible columns using a pipette and collect the unlabeled cells in flow-through.

4.5 Wash the column with Isolation Buffer (500 μ L \times 3 for CytoSinct™ gM Column, 3 mL \times 3 for CytoSinct™ gL Column, or other compatible columns). Collect unlabeled cells in flow-through with

a suitable tube (for example, a 2 mL or 15 mL conical tube). Repeat the washing step for another two times. Add new Isolation Buffer when the column stops dripping but not run dry.

4.6 Remove the column from the magnet and place it on a new tube with suitable size (for example, a 15 mL or 50 mL conical tube).

4.7 Pipette the Isolation Buffer onto the column (1 mL for CytoSinct™ gM Column, 5 mL for CytoSinct™ gL Column, or other compatible columns). Immediately flush out the fraction with the magnetically labeled cells by firmly applying the plunger through the column chamber supplied with the column.

4.8 The cells can then be counted, analyzed to assess the purity or used in down-stream applications. The Nanobeads do not need to be removed. To ensure cell viability, the desired cell fraction should be immediately resuspended in cell culture medium

For research use only. Not intended for human and animal therapeutic or diagnostic use.

生产商：南京金斯瑞生物科技有限公司 江苏省南京市江宁区科学园雍熙路 28 号

Manufacturer: Nanjing GenScript Biotech Co., Ltd. No. 28 Yongxi Road, Jiangning District, Nanjing, Jiangsu, China