

MojoSort™ Human NK Cell Isolation Kit

Catalog# / Size	480053 / 20 tests 480054 / 200 tests
Regulatory Status	RUO
Description	Non-natural killer (NK) cells are depleted by incubating the sample with the biotin antibody cocktail followed by incubation with magnetic Streptavidin Nanobeads. The magnetically labeled fraction is retained by the use of a magnetic separator. The untouched NK cells (CD56 ⁺ , CD3 ⁻) are collected by decanting the liquid in a clean tube. These are the cells of interest; do not discard the liquid. Some of the downstream applications include functional assays, gene expression, phenotypic characterization, etc.

Kit Contents

Kit Contents	For Cat# 480053: <ul style="list-style-type: none">• 200 µl Biotin-Antibody Cocktail• 200 µl Streptavidin Nanobeads For Cat# 480054: <ul style="list-style-type: none">• 2 vials of 1 ml Biotin-Antibody Cocktail each• 2 vials of 1 ml Streptavidin Nanobeads each
---------------------	---

Product Details

Verified Reactivity	Human
Formulation	Cocktail: Phosphate buffer solution containing 0.09% sodium azide, pH 7.2. Streptavidin Nanobeads: Aqueous solution containing BSA and 0.05% sodium azide.
Preparation	The antibodies were purified by affinity chromatography, and conjugated with biotin under optimal conditions. Streptavidin Nanobeads: Streptavidin-coated magnetic beads.
Storage & Handling	Antibody cocktail and Streptavidin Nanobeads should be stored undiluted between 2°C and 8°C.
Application	Cell Separation (MojoSort™) - Quality tested
Recommended Usage	10 µl of antibody cocktail for 1 x 10 ⁷ cells in 100 µl of buffer. 10 µl of Streptavidin Nanobeads for 1 x 10 ⁷ cells in 100 µl of buffer.
Application Notes	This kit is designed for the isolation of untouched Natural Killer (NK) Cells from peripheral blood mononuclear cells (PBMCs). Each lot has been individually optimized. Do not mix and match components from different lots or different kits. Antibody or cocktail dilution to use in column: 4X Nanobead dilution to use in columns: 4X
Product Citations	1. Harmon C, <i>et al.</i> 2019. Front Immunol. 10:1502. PubMed

Antigen Details

Biology Area	Immunology, Innate Immunity
---------------------	-----------------------------

Molecular Family CD Molecules
Gene ID NA

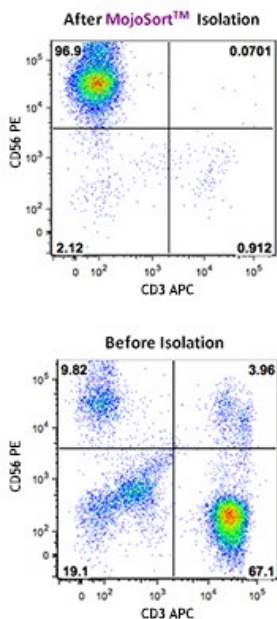
Related Protocols

[MojoSort™ Human NK Cell Isolation Kit Protocol](#)

[MojoSort™ Human NK Cell Isolation Kit Column Protocol](#)

[MojoSort™ General Protocol - Video](#)

Product Data



PBMCs were prepared to isolate CD56⁺ CD3⁻ cells using the MojoSort™ Human NK Cell Isolation Kit. Cells were stained with CD56 (clone hCD56) PE, CD3ε (UCHT1) APC. Dead cells were excluded by 7-AAD.

For research use only. Not for diagnostic use. Not for resale. BioLegend will not be held responsible for patent infringement or other violations that may occur with the use of our products.

*These products may be covered by one or more Limited Use Label Licenses (see the BioLegend Catalog or our website, www.biolegend.com/ordering#license). BioLegend products may not be transferred to third parties, resold, modified for resale, or used to manufacture commercial products, reverse engineer functionally similar materials, or to provide a service to third parties without written approval of BioLegend. By use of these products you accept the terms and conditions of all applicable Limited Use Label Licenses. Unless otherwise indicated, these products are for research use only and are not intended for human or animal diagnostic, therapeutic or commercial use.

8999 BioLegend Way, San Diego, CA 92121 www.biolegend.com
Toll-Free Phone: 1-877-Bio-Legend (246-5343) Phone: (858) 768-5800 Fax: (877) 455-9587