

MojoSort™ Human CD3 T Cell Isolation Kit

Catalog# / Size	480021 / 10 tests 480022 / 100 tests 480131 / 200 tests
Regulatory Status	RUO
Description	Non CD3 ⁺ T cells are depleted by incubating your 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 CD3 ⁺ T cells are collected by decanting the liquid in a clean tube. These are your 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# 480021: <ul style="list-style-type: none">• 100 µl Biotin-Antibody Cocktail• 100 µl Streptavidin Nanobeads For Cat# 480022 <ul style="list-style-type: none">• 1 ml Biotin-Antibody Cocktail• 1 ml Streptavidin Nanobeads
---------------------	--

Product Details

Verified Reactivity	Human
Formulation	Cocktail: Phosphate-buffered 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 1x10 ⁷ cells in 100 µl of buffer. 10 µl Streptavidin Nanobeads for 1x10 ⁷ cells in 100 µl of buffer. For this kit, incubation with antibody cocktail and Streptavidin Nanoparticles can be done for 5 to 15 minutes, without loss in purity and yield.
Application Notes	This kit is designed for the isolation of untouched CD3 ⁺ T 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: 2.5X Nanobead dilution to use in columns: 4X
Additional Product Notes	View more applications data for this product in our Scientific Poster Library . Watch a Scientific Poster video of this product.

Product Citations

1. Rennie K, *et al.* 2020. Clin Cancer Res. 26:5019. [PubMed](#)

Antigen Details

Biology Area	Immunology
Molecular Family	CD Molecules, TCRs
Gene ID	NA

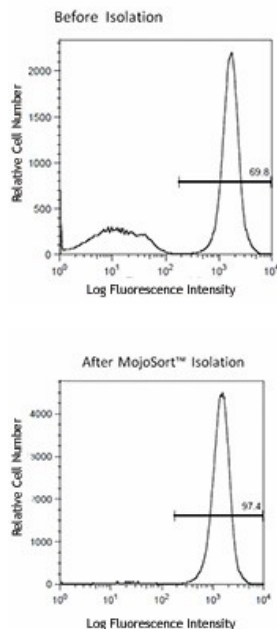
Related Protocols

[MojoSort™ Isolation Kits Protocol - 1](#)

[MojoSort™ Isolation Kits Column Protocol - 1](#)

[MojoSort™ General Protocol - Video](#)

Product Data



PBMCs were prepared to isolate CD3⁺ T cells using the MojoSort™ Human CD3 T Cell Isolation Kit. Cells were stained with CD3ε (UCHT1) APC.

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.

BioLegend Inc., 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