



# Advanced Products for Cell Therapy Overview

Application Note

# CONTENTS

Cell Therapy Overview	03
Workflow Example and Applicable Solutions	04
Cryopreservation and Thawing	05
Benz-Neburase™ Nuclease	
Cell Isolation	05
CytoSinct™ System	
T Cell Isolation	
NK Isolation	
Stem Cell Isolation	
TCR Removal	
Automated Cell Isolation with Data Integrity	
T Cell Activation	08
Optimized T Cell Activation with Enceed™	
Transduction	08
Accurate & Fast Lentiviral Titration with p24 ELISA Kit Pro	
Expansion	09
Catalog Cytokines and Growth Factors	
Analytical/Quality Control in Cell Therapy	09
Flow Cytometry Analysis of CAR+ Cells with MonoRab™ anti-scFv or anti-VHH Antibodies	
Endotoxin Detection with ToxinSensor™ Chromogenic LAL Endotoxin Assay Kit	
BSA Detection ELISA Kit, 2G	
Regulatory Support for cGMP Products	11
References	11

# OVERVIEW

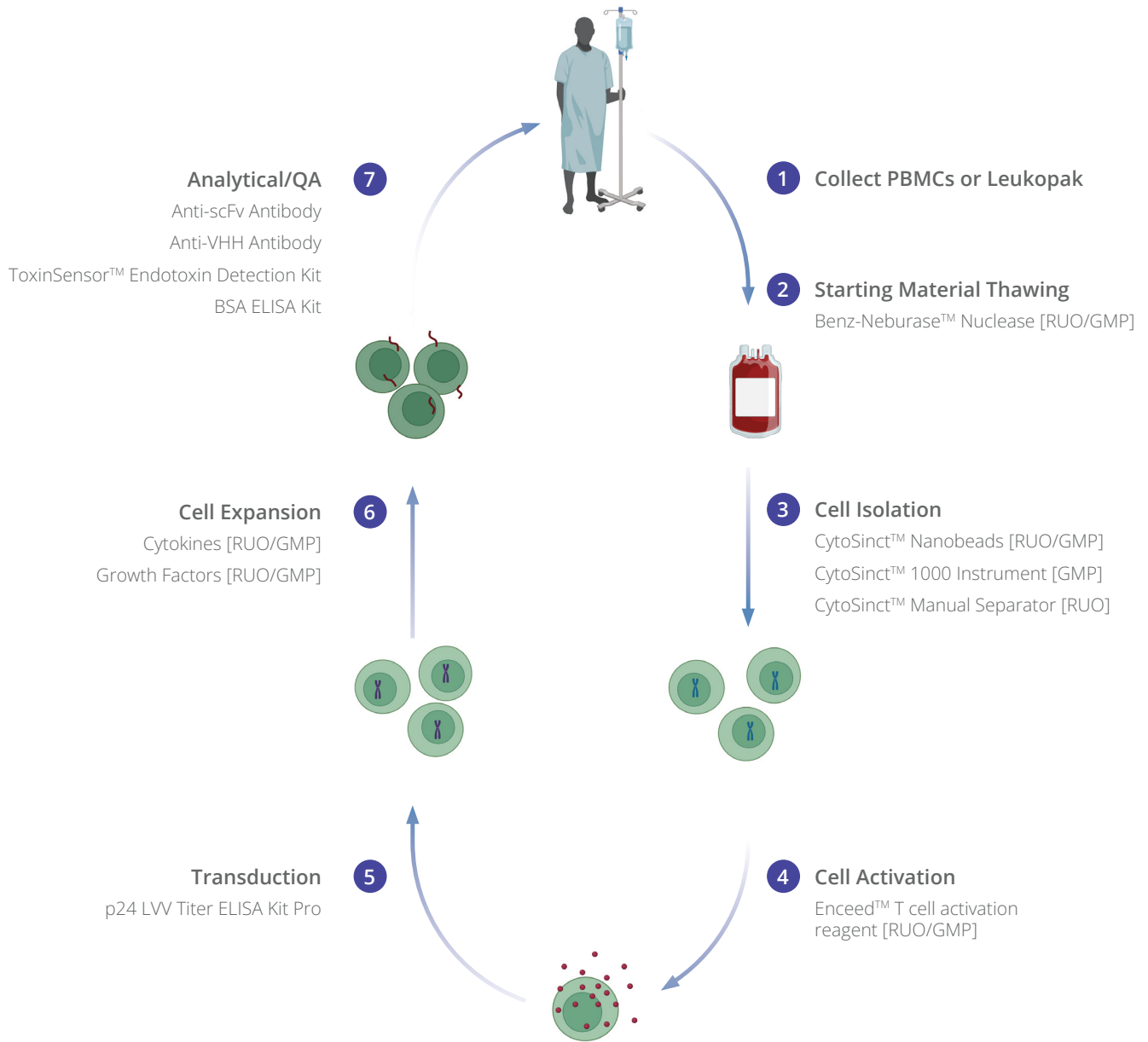
Cell based therapies have emerged as one of the most promising approaches in modern medicine, offering novel solutions for a wide range of diseases, from cancer to autoimmune disorders and regenerative medicine. These therapies harness the power of the body's immune and regenerative systems, with various types of cells, including T cells, NK cells, macrophages, and stem cells, playing pivotal roles in targeted treatments. Advances in genetic engineering and in vitro expansion techniques have enhanced the efficacy of these therapies, with some already receiving regulatory approval and others showing promise in clinical trials. Below is an overview of key cell based therapies currently shaping the field, their mechanisms, and recent developments.

- CAR T cell therapy has shown great promise in recent years. As of 2024, six CAR T cell therapies targeting hematological malignancies have been approved by the FDA since the first approval in 2017. Some of these therapies are now being used in earlier lines of treatment, expanding their target populations. Recently, CAR T cell therapy has also demonstrated encouraging clinical outcomes in autoimmune diseases<sup>[1]</sup>.
- TCR T cell therapy represents another promising T cell based approach, particularly in the treatment of solid tumors. This is due to certain intrinsic advantages of TCRs, such as a target repertoire not limited to membrane proteins and their high antigen sensitivity<sup>[2]</sup>.
- $\gamma\delta$  T cells, while accounting for less than 5% of the total T cell population, have emerged as a highly promising effector cell type for cancer immunotherapy. Their ability to recognize and kill transformed cells independently of HLA antigen presentation is key to their therapeutic potential. Current strategies for  $\gamma\delta$  T cell therapies involve developing both autologous and allogeneic approaches, typically including in vitro expansion using cytokines and CD3 monoclonal antibodies, with potential incorporation of CAR engineering or gene modification to enhance chemotherapy resistance<sup>[3]</sup>.
- Natural killer (NK) cells, a vital component of the innate immune system, play a key role in identifying and eliminating distressed cells, including tumor and virus infected cells. NK cells exert direct effector functions against these targets and support the broader immune response. Advances in receptor engineering have led to significant progress in NK cell based cancer immunotherapies<sup>[4],[5]</sup>.
- Monocyte derived, ex vivo generated macrophages have also gained attention in recent years, showing superior performance over stem cells in certain clinical trials. The development of chimeric antigen receptor (CAR) macrophages has further highlighted the potential of genetically engineered macrophages in advancing cell therapy<sup>[6]</sup>.
- Tumor infiltrating lymphocyte (TIL) therapy is an adoptive cell transfer treatment, uses a patient's own immune cells to fight cancer. This involves harvesting TILs from the patient's tumor, expanding those with high specificity for tumor antigens in vitro, and reinfusing them to enhance antitumor responses. In 2024, TIL therapy received FDA approval for the treatment of relapsed or refractory melanoma and is currently being tested in clinical trials for other cancers<sup>[7]</sup>.
- Stem cell therapy is a offers regenerative treatment options for various diseases and conditions. Hematopoietic stem cell transplantation (HSCT), commonly known as a bone marrow transplant, involves isolating stem cells from a healthy donor and infusing them into a patient with dysfunctional or depleted bone marrow to treat hematologic disorders or malignancies. Mesenchymal stem cell therapy has also gained worldwide attention due to the cells' multipotency in tissue regeneration and their strong immunosuppressive properties<sup>[8]</sup>. This approach typically involves isolating and expanding cells in vitro before infusion back into the patient.

Despite the promising advancements in cell based therapies, several challenges remain. Manufacturing scalability and high production costs are significant hurdles, particularly in personalized therapies like CAR T and TIL, which require complex in vitro expansion processes. Ensuring long term persistence of therapeutic cells and managing potential off target effects or immune related toxicities are also critical concerns.

GenScript is addressing these challenges by offering comprehensive solutions that streamline the manufacturing process, improve scalability, and reduce costs. Through innovations in automated cell processing systems, gene editing technologies, and robust quality control protocols, GenScript is helping to optimize cell therapy workflows and improve the accessibility and efficiency of these life life-saving treatments.

# Cell Therapy Workflow Example and Applicable Solutions



## Cryopreservation and Thawing

Cell clumping is a major challenge in cell therapy manufacturing, impacting cryopreservation, thawing, and expansion by reducing viability and therapeutic efficacy. It complicates downstream processing, affecting cell isolation, quality, and scalability, while increasing contamination risks.

### Benz-Neburase™ Nuclease For Safely Reducing Cell Clumping

Benz-Neburase™ *S. marcescens* Nuclease is a genetically engineered endonuclease derived from *Serratia marcescens*, designed for efficient digestion of all forms of DNA and RNA into 3-5 base nucleotides. It is highly effective in reducing nucleic acids without impacting cell viability, even during extended treatments. Available in both RUO and GMP grades, the tag-free Benz-Neburase™ offers a safe and efficient solution for reducing cell clumping, particularly when thawing frozen cell samples.

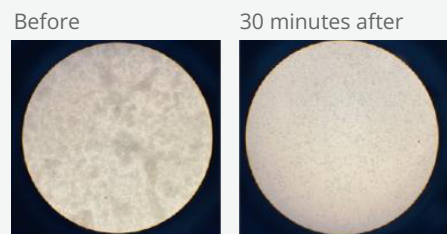


Figure 1: Benz-Neburase efficiently reduced cell clumping from SUP-T1 cell culture 30 minutes after treatment at 50 units/mL, 37°C.

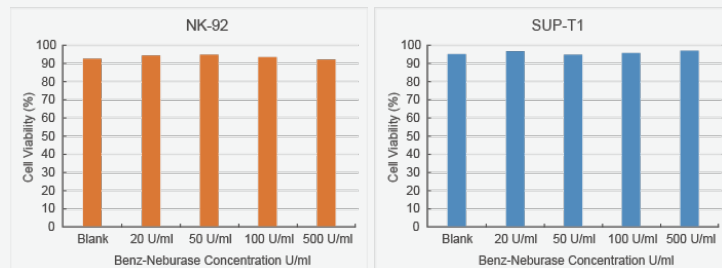


Figure 2: Benz-Neburase has minimal to no impact on the cell viability after incubation with each time of cell line over night at 20-500 units/mL, 37°C.

## Cell Isolation

Cell isolation is a critical step in the development and manufacturing of cell-based therapies. Precise isolation of target cells based on surface markers ensures the purity and potency of the therapeutic product while minimizing contamination risks. Traditional methods, such as density gradient centrifugation and FACS, are labor-intensive, time-consuming, and may not scale well or achieve high specificity. Advanced immuno-magnetic technologies, using magnetic beads coated with antibodies, offer more efficient, scalable, and targeted cell isolation.

### Cell Isolation with CytoSinct™ System

Leveraging GenScript's expertise in antibody development and magnetic bead production, CytoSinct™ Nanobeads are nano-sized, biodegradable, column-compatible, and do not require de-beading, ensuring efficient isolation of both common and rare cell populations. Available in RUO and GMP grades, CytoSinct™ Nanobeads, along with manual and automated instruments, support seamless scaling from research to manufacturing.

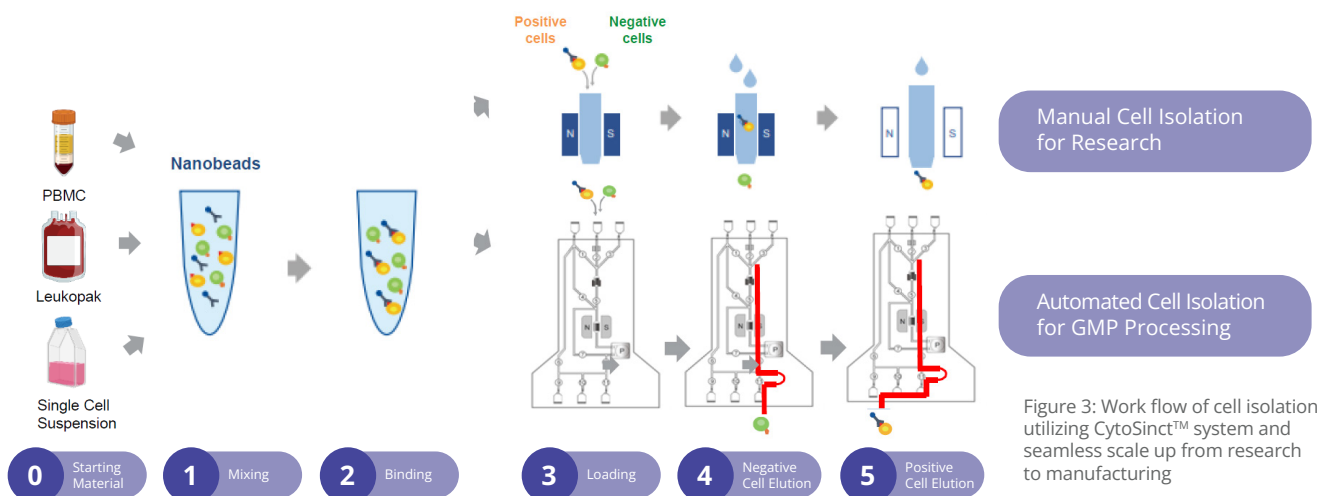


Figure 3: Work flow of cell isolation utilizing CytoSinct™ system and seamless scale up from research to manufacturing

## CD4+/CD8+ T Cell Isolation

Purified T cell subsets can enhance the biological activity of the final CAR T cell therapy product, in line with FDA recommendations<sup>[9]</sup>. CD4+ and CD8+ combined isolation is one of the most commonly used strategies for T cell enrichment due to its ability to avoid T cell activation. The co-incubation and single-step elution of both CD4+ and CD8+ cells using CytoSinct™ Nanobeads significantly reduces processing time compared to sequential isolation (CD4+ followed by CD8+), while maintaining the same high level of purity. In addition, CD4+ cell isolation with CytoSinct™ Nanobeads has shown a trend of higher recovery rates compared to the current standard method.

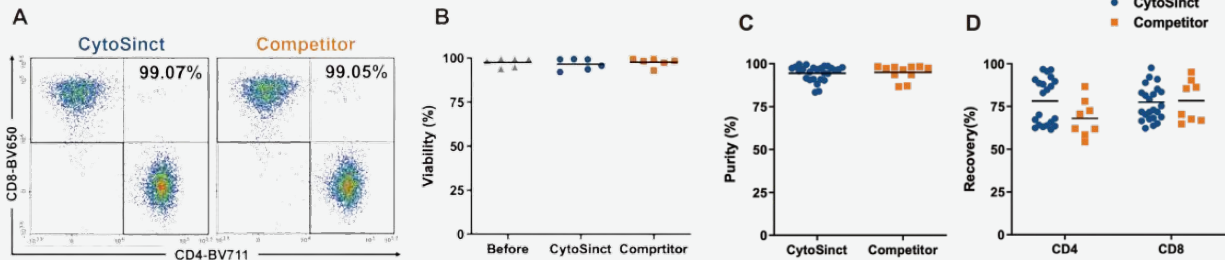


Figure 4. T cells isolation from various healthy donor leukopheresis samples using CytoSinct™ CD4 and CD8 Nanobeads and gM columns or the current standard method. (A) Representative flow analysis of CD4 and CD8 population and combined percentage after isolation. (B) Viability before and after isolation. (C) Purity of total T cells (CD4+ and CD8+ combined) after isolation (D) Recovery after isolation.

## NK Cell Isolation

Efficient enrichment of NK cells is critical for producing NK cell based therapies, such as CAR NK. However, isolating NK cells is more challenging than isolating T cells due to their relatively low abundance, typically around 10% in starting materials like leukopak or PBMCs. This process can be optimized by first depleting CD3+ cells, followed by enrichment of CD56+ cells, ensuring high purity without inducing NK cell activation.

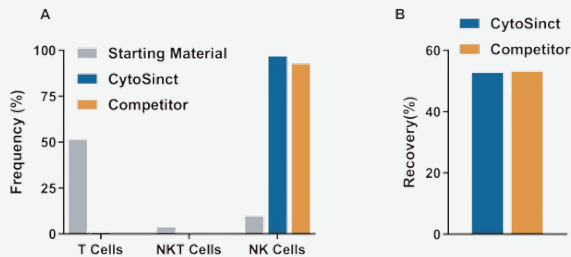


Figure 5. Representative flow analysis of NK (CD3-CD56+), NKT (CD3+ CD56+) and T cells (CD3+) following CD3+ cell removal and CD56+ cell enrichment from a leukapheresis sample using CytoSinct™ CD3 and CD56 Nanobeads and gM columns. (A) High purity of NK cells and sufficient removal of other cell types after isolation. (B) Recovery of NK cells (CD56+ CD3 -) after isolation.

## CD34+ Stem Cell Isolation

Isolating CD34+ stem cells presents a significant challenge due to their extremely low percentage, typically around 1%, in starting materials such as cord blood and mobilized peripheral blood. These sources also contain high levels of red blood cells and platelets, which can interfere with the isolation process. To achieve optimal results, PBMC enrichment is often required before magnetic isolation. Following this step, high purity stem cells can be efficiently isolated using CytoSinct™ CD34 Nanobeads.

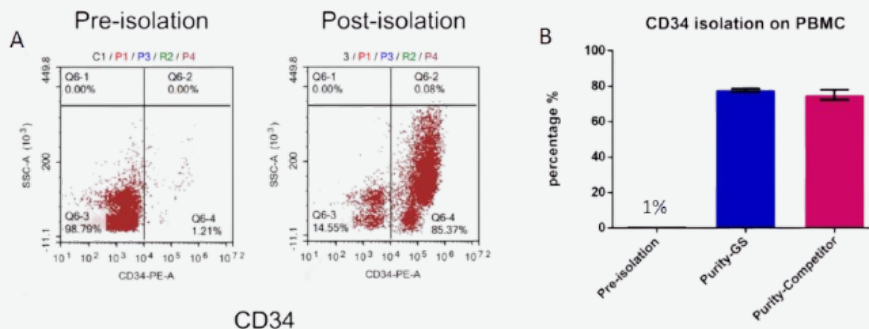


Figure 6. CD34+ cell line enrichment from a mix with PBCMs using CytoSinct™ CD34 Nanobeads and gL columns (A) Representative flow cytometry analysis before and after isolation (B) purity before and after isolation using CytoSinct™ CD34 Nanobeads or the current standard method.

## One-Step TCR αβ Depletion For Allogeneic Cell Therapy

In allogeneic cell therapy, a key challenge is the risk of graft versus host disease (GvHD), where the transplanted cells attack the host, leading to adverse reactions. αβ T cell is the primary immune cell type that cause acute GvHD, and its removal at the end of the production process is an effective measure for allogeneic cell therapy to alleviate GvHD. Therefore, αβ T cell removal is a regulatory mandate by FDA for allogeneic CAR T cell therapy production. Additionally, αβ T cell depletion is being increasingly studied in allogeneic hematopoietic stem cell transplantation (allo HCT) to overcome the human leukocyte antigen (HLA) barrier in haploidentical (haplo) HCT, showing promising results<sup>[10]</sup>.

CytoSinct™ TCR αβ Nanobeads provide an efficient solution for removing αβ T cells from various allogeneic cell products, including but not limited to allogeneic CAR T, γδ T cells, stem cells and CAR NK samples, while preserving the remaining cell components. Unlike standard methods, CytoSinct™ TCR αβ Nanobeads come pre conjugated and ready to use, eliminating the need for additional steps such as labeling, washing, and resuspension. This simplified process saves 30 minutes to 2 hours of processing time, potentially reduces the use of one tubing set, and enhances cell recovery by 10-20% (depending on whether using RUO or GMP process and the choice of cell washing instrument).

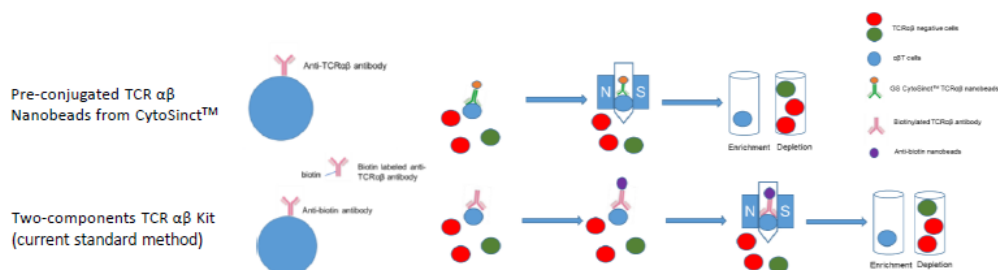


Figure 7. Component(s) and workflow of CytoSinct™ TCR αβ beads or the current standard method

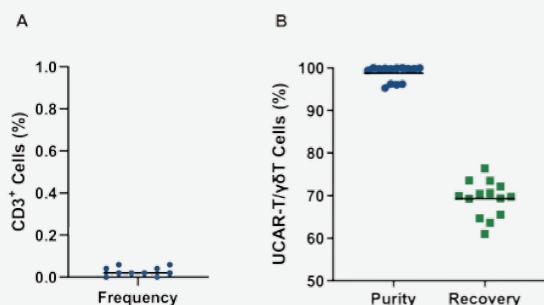


Figure 8. Flow analysis of various donor derived allogeneic CAR T cells or γδ T cells after TCR αβ depletion using CytoSinct™ TCR αβ Nanobeads and gL columns. (A) Percentage of CD3+ cells below 0.1% after TCR αβ depletion. (B) Purity and recovery of CD3 cells after TCR αβ depletion.

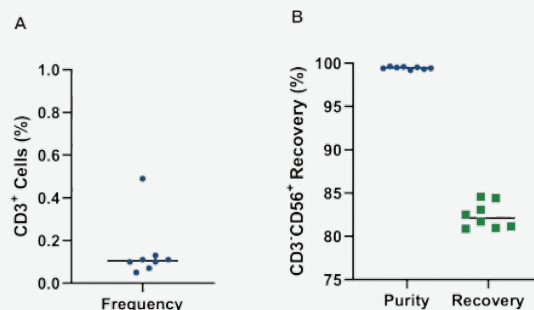
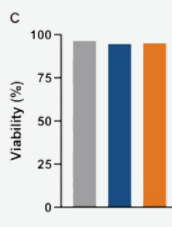
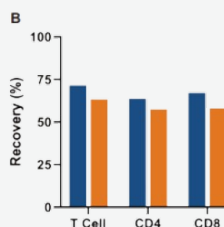
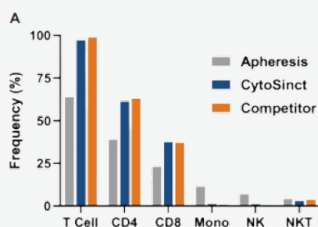


Figure 9. Flow analysis of various donor derived allogeneic CAR NK cells after TCR α/β depletion using CytoSinct™ TCR αβ Nanobeads and gL columns. (A) Percentage of CD3+ cells mostly below 0.1% after TCR αβ depletion. (B) Purity and recovery of CD3 cells after TCR αβ depletion.

## Closed, Automated Cell Isolation for Cell Therapy Manufacturing with Data Integrity

Automated, large scale cell isolation in a closed system is essential for successful GMP manufacturing and clinical application of cell therapy products. The CytoSinct™ 1000 is designed with 21 CFR Part 11 compliance, adhering to FDA regulations for electronic records and e-signatures, ensuring data integrity throughout the cell isolation process. Its customizable programming and alert functions streamline the operation, providing an efficient and reliable solution for automated cell isolation with peace of mind.



Test Groups	CytoSinct	Competitor
Instrument	CytoSinct™ 1000	Current Standard Method
Tubing Set	CytoSinct™ 1000 Tubing Set	Current Standard Method
Reagent	CytoSinct™ CD4 and CD8 Nanobeads	CytoSinct™ CD4 and CD8 Nanobeads

Figure 10. T cells isolation from various donor leukapheresis samples using CytoSinct™ 1000 instrument and tubing set and similar instrument and tubing set from another provider, both with CytoSinct™ CD4 and CD8 Nanobeads. (A) Subpopulation analysis. (B) Recovery after isolation. (C) Viability before and after isolation.

## T Cell Activation

Effective and gentle activation of T cells is critical for achieving high cell count in final product while also keeping young phenotype for persistence and clinical efficacy after infusion. Selected stem like or “young” T cells can significantly influence the biological activity of the final therapy, according to FDA guidelines<sup>[9]</sup>.

### Optimized T Cell Activation with Enceed™ for Enhanced Memorial Phenotype & Consistency

Pre-conjugated with both anti-CD3 and anti-CD28 antibodies, Enceed™ T cell activation reagent is designed to activate T cells by engaging the T cell receptor (TCR) and the CD28 co-stimulatory module. Thanks to its nanometer size, unbound Enceed™ can be easily removed through cell washing or a medium change without the need for a magnet. T cells activated by Enceed™ present enhanced memory phenotype and greater expansion compared to those activated using standard methods. The consistent performance of Enceed™ across both RUO and GMP grades enables seamless scale up from research to manufacturing, ensuring its utility for large-scale T cell based therapies.

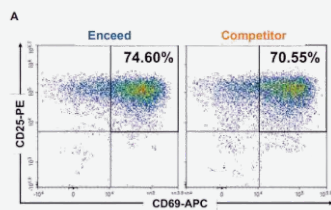


Figure 11. Flow analysis of cell activation markers CD25 and CD69 expression on T cells enriched from one (A) or multiple (B) healthy donor PBMC samples 2 days after treated with Enceed™ T cell activation reagent or using the current standard method.

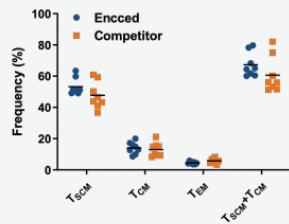


Figure 13. Enhanced memorial phenotype T cells from healthy donor PBMC samples 10 days after treated with Enceed™ and cultured with IL-2.

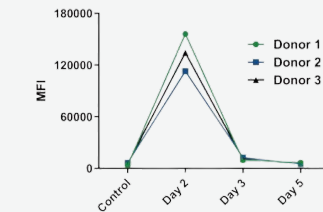
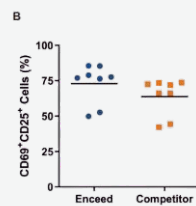


Figure 14. Flow cytometry analysis showing removal of cell surface Enceed™ by centrifugation 2 days after treatment.

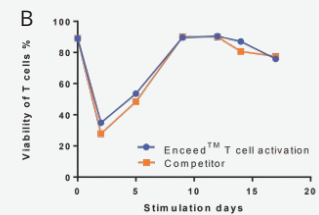
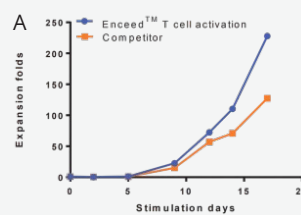


Figure 12. (A) Fold expansion and (B) viability of T cells enriched from a healthy donor PBMC after activated by Enceed™ or the current stand method and cultured with IL-2 for 17 days.

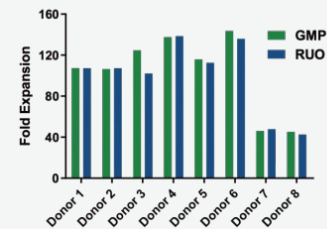


Figure 15. Consistent T cell expansion between GMP & RUO grade Enceed™ from various health donor derived PBMCs and after 10 days of culture with IL-2.

## Transduction

Lentiviral transduction is a crucial technique in cell therapy that involves using lentivirus vectors (LVV) to introduce genetic material into target cells. This method allows for stable and long term expression of therapeutic genes, making it a powerful tool for developing treatments for various genetic disorders and cancers.

### Accurate and Fast Lentiviral Titration with p24 ELISA Pro Kit

ELISA is a widely used method for the physical titration of lentiviral particles by quantifying p24 protein, offering a quicker and simpler alternative to measuring lentiviral RNA via qPCR. However, traditional ELISA may lack accuracy as it measures all p24 in the sample, including free p24 not incorporated into viral particles. The p24 ELISA Pro kit overcomes this limitation by using proprietary magnetic bead technology to remove free p24, enhancing specificity and accuracy. It supports a broad linear range of 30 to 2,000 pg/mL, reducing dilution requirements, and can be completed in just 2.5 hours, allowing for multiple runs per day as needed.

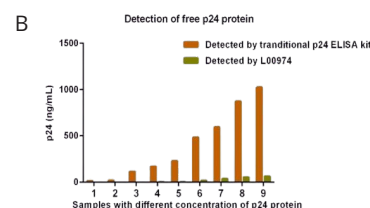
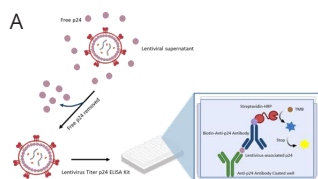


Figure 16. Highly accurate lentiviral titration by p24 ELISA assay through removal of free p24 protein from sample using magnetic beads. (A) schematic diagram of p24 Pro ELISA kit workflow. (B) Free p24 added to dilution buffer was detected by traditional P24 ELISA kit but not p24 Pro kit.



## Cell Expansion

In cell therapy manufacturing, the in vitro expansion of therapeutic cells is a critical step that relies heavily on the use of cytokines and growth factors. These essential ancillary materials drive the growth and proliferation of various cell types. Cytokines like IL-2, IL-15, and IL-7 are indispensable for lymphocyte based therapies, such as CAR T, TCR T, NK cells, and tumor infiltrating lymphocytes (TILs)<sup>[7]</sup>. Additionally, growth factors such as basic fibroblast growth factor (bFGF) have been shown to significantly enhance mesenchymal stem cell (MSC) expansion in vitro<sup>[8]</sup>.

### Optimized Cell Expansion with GenScript Catalog Cytokines and Growth Factors

GenScript offers a comprehensive catalog of recombinant cytokines and growth factors at both RUO and GMP grades, designed to support efficient cell expansion for both research and large scale therapeutic manufacturing. These products are characterized by superior activity, excellent lot to lot consistency, and remarkably low endotoxin levels, ensuring reliable and effective cell growth in any application.

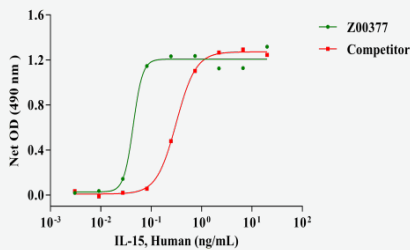


Figure 17. GenScript IL-15 (Cat.No. Z00377) ED50 < 0.5 ng/ml as determined by the dose-dependent stimulation of the proliferation of CTL2-2 cells, corresponding to a specific activity of > 2.0 × 10<sup>6</sup> IU/mg.

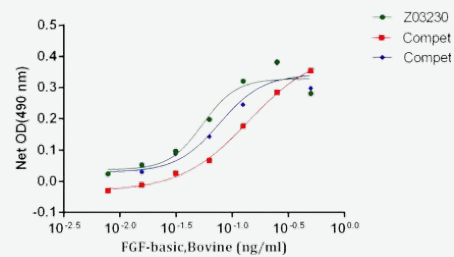


Figure 18. GenScript FGF-basic (Cat.No. Z03230) ED50 < 1.0 ng/ml as determined by the dose-dependent proliferation assay using 3T3 cells, corresponding to a specific activity of > 1.0 × 10<sup>6</sup> units/mg.

## Analytical/Quality Control in Cell Therapy

### Flow Cytometry Analysis of CAR+ cells with MonoRab™ Anti-scFv and Anti-VHH Antibodies

Characterization of CAR expression, CAR+ cell count, and percentage is critical for CAR engineered cell therapy research, GMP manufacturing in process analysis, and lot release. The FDA recommends directly detecting CAR to accurately determine the percentage of CAR positive cells. Additionally, if surrogate proteins (co expressed genes) or broad specificity reagents (e.g., protein L) are used, their correlation with CAR expression should be evaluated<sup>[9]</sup>. Single chain variable fragments (scFv) and variable heavy chain domains (VHH) are commonly used CAR constructs in both research and clinical settings. MonoRab™ Anti-scFv and Anti-VHH antibodies are specifically developed for the detection of scFv and VHH CAR constructs, offering high specificity, affinity, and sensitivity across a variety of constructs. These antibodies provide non-blocking detection, ensuring that they do not interfere with the CAR's ability to bind its target cells, making them ideal for downstream application like functional assays.

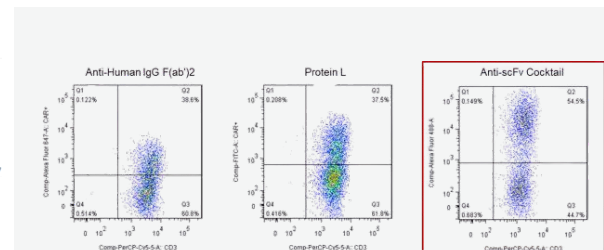
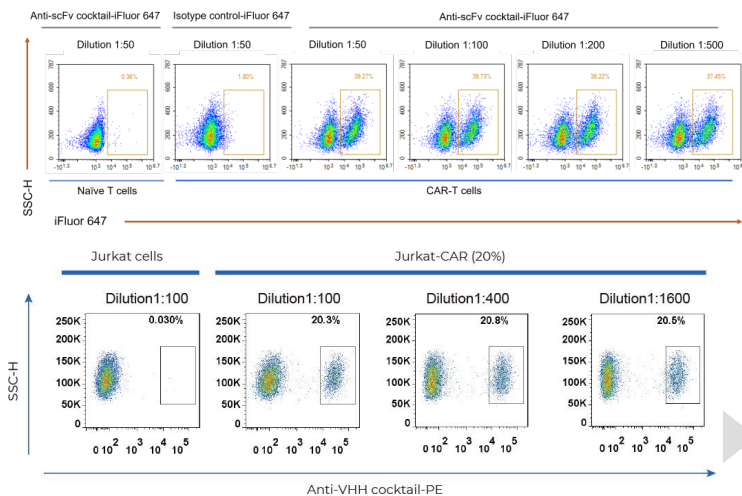


Figure 19. Flow analysis in comparison of scFv CAR T cell detection using anti-IgG, protein-L, and MonoRab™ anti-scFv antibody cocktail.

Figure 20. Flow analysis shows scFv CAR+ T cells and VHH CAR+ Jurkat cells detected by MonoRab™ Rabbit Anti-scFv Cocktail [iFluor 647] (Cat. No. A02288) and Anti-Camelid VHH Cocktail [PE] (Cat. No. A02018) with high accuracy at various dilution ratios.

## Highly Sensitive Endotoxin Detection with ToxinSensor™ Chromogenic LAL Endotoxin Assay Kit

Endotoxins, which are lipopolysaccharides from gram negative bacteria, can cause severe adverse reactions in patients. Therefore, it is crucial that cell therapy products and viral vectors used for gene editing are tested for endotoxin levels according to current good manufacturing practices (cGMP) and FDA regulations<sup>[9]</sup>. This can be done by the limulus amoebocyte lysate (LAL) test, an FDA approved method, with highly sensitive detection of bacterial endotoxins.

The ToxinSensor™ Chromogenic LAL Endotoxin Assay Kit (Cat. No. L00350) allows for quantitative detection of endotoxins over a wide range, with high sensitivity (0.01 1 EU/ml) and strong reproducibility. It also includes ready-to-use, endotoxin-free reagents and materials, ensuring reliable and efficient testing for endotoxins.

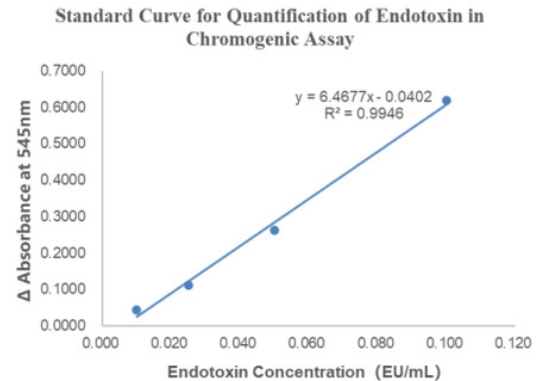


Figure 21. Representative standard curve showing wide linear range. Note the OD values of standards may vary depending on the assay.

## Highly Sensitive BSA Detection with GenScript BSA ELISA Kit 2G

Highly Sensitive BSA Detection with GenScript BSA ELISA Kit 2G Fetal bovine serum (FBS) is widely used in large-scale mammalian cell culture and serves as a key raw material in the production of biological products, including cell and gene therapies. Many commercial serum-free media formulations contain biological macromolecules, such as bovine serum albumin (BSA), transferrin, and insulin, to replace the functions of serum. However, as a heterologous protein, BSA can trigger adverse immune reactions in humans if present above a certain threshold, impacting the safety of biological products. Therefore, detecting BSA residues is a critical quality control measure in the production of biologics.

The GenScript BSA ELISA Kit 2G (Cat. No. L00976) is a highly sensitive sandwich ELISA kit designed for the quantitative measurement of BSA in test samples. With a limit of blank (LoB) of 0.14 ng/mL and a limit of quantification (LoQ) of 0.5 ng/mL, this kit offers exceptional sensitivity, accuracy, and stability. Its traceable standard and calibration make it an ideal solution for routine BSA impurity detection, quality control, lot release, and the optimization of purification processes.

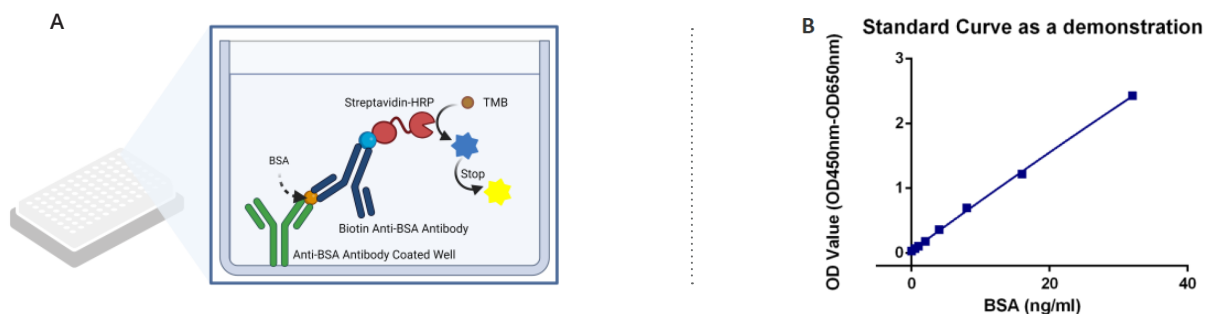


Figure 22.(A) Schematic diagram of BSA ELISA Kit 2G working mechanism. (B) Representative standard curve showing wide linear range of 0.5 35 ng/mL.

## Regulatory Support

---



### GenScript Biologics Manufacturing Center for GMP Grade Products

The GenScript Manufacturing Center in Zhenjiang China, spans over 100,000 square meters. The center is designed and constructed to fully comply with the good manufacturing practice (GMP) standards set by the FDA, EMA, PMDA, and NMPA. This state of the art facility is equipped to meet the demands of Phase I to III clinical sample production, as well as commercial scale manufacturing.

GenScript's quality management system is certified to ISO 13485 standard, ensuring rigorous processes for the design, development, manufacture, and distribution of GMP products. Drug Master Files (DMF) for GenScript's GMP grade products have been submitted to the FDA, supporting Investigational New Drug (IND) applications in the U.S.

## References

---

- [1] : The changing landscape of cancer cell therapies: clinical trials and real world data. *Nature Reviews Drug Discovery* , 2024 May 31.
- [2] : TCR engineered T cell therapy in solid tumors: State of the art and perspectives. *Science Advances* , 2023 Feb 15, Vol 9, Issue 7.
- [3] : Gamma Delta T Cell Based Cancer Immunotherapy: Past Present Future. *Frontiers in Immunology* , 2022 Jun 16:13:915837.
- [4] : Natural killer cell therapies. *Nature*, volume 626, pages727 736 (2024).
- [5] : Development of NK cell based cancer immunotherapies through receptor engineering. *Cellular & Molecular Immunology*, volume 21, pages315 331 (2024).
- [6] : A new era of macrophage based cell therapy. *Experimental & Molecular Medicine*, volume 55, pages 1945 1954
- [7] : Expansion of tumor infiltrating lymphocytes (TIL) from human pancreatic tumors. *J Immunother Cancer* , 2016 Oct 18:4:61.
- [8] : Large Scale Expansion of Human Mesenchymal Stem Cells. *Stem Cells Int*, 2020: 9529465.
- [9] : Considerations for the Development of Chimeric Antigen Receptor (CAR) T Cell Products Guidance for Industry. FDA, 2024 Jan.
- [10] : Role of  $\alpha\beta$  T Cell Depletion in Prevention of Graft versus Host Disease. *Biomedicines*. 2017 Sep; 5(3): 35.

## CONTACT US

### USA

Tel: +1 732 885 9188 Toll Free: +1 877 436 7274

Fax: +1 732 210 0262

Email: [product.techsupport@genscript.com](mailto:product.techsupport@genscript.com)

Address: 860 Centennial Ave. Piscataway NJ 08854

### China

Tel: +86-025-58897288 ext 5810

Fax: +86-025-58897288-5815

Email: [product@genscript.com.cn](mailto:product@genscript.com.cn)

Address: 28 YongXi Road, JiangNing District, Nanjing, China

### Asia/Pacific

Tel: +65 3159 1898

Email: [sales.AP@genscript.com](mailto:sales.AP@genscript.com)

Fax: +65 6491 5073

Address: 164 Kallang Way, #06-12, Singapore 349248

### Europe

Tel: +31 (0) 71 569 0120

Email: [support.EU@genscript.com](mailto:support.EU@genscript.com)

Address: Treubstraat 1, 2288Eg Rijswijk Zh Netherlands

