

Background

Protein biomarker assays are well-established tools for interrogating biological responses to treatment. Modulation of context-relevant biomarkers allows researchers to gain insights on treatment-relevant dynamics. Herein, we discuss the development of four multiplex biomarker assays designed to monitor critical attributes associated with chimeric antigen receptor T cells (CAR-T). These U-PLEX® Combos include markers useful in research regarding efficacy, safety, persistence and exhaustion in cultured cell models and biofluids.

Engineered cellular therapies such as CAR-T have revolutionized the management of hematological malignancies and solid tumors. Multiple CAR-T therapies (including Kymriah, Yescarta, Breyanzi, Carteyva, and Tecartus) were approved by the FDA between 2017 and 2023. Patients treated with CAR-T often experience remarkable, relapse-free survival. These immune cell therapies demonstrate prolonged CAR-T cell persistence, enhanced CAR-T effector function, limited T cell exhaustion, and good safety profiles in many responding patients. The most common adverse reactions in patients receiving CAR-T therapy are cytokine release syndrome (CRS) and immune effector cell-associated neurotoxicity syndrome (ICANS).¹⁻⁵ These adverse events may lead to patients being treated with highly immunosuppressive corticosteroids to mitigate effects, hospitalization, treatment discontinuation, and in rare instances, death.

Table 1. Definitions of CAR-T attributes.

Term	Definition
Efficacy	CAR-T cell engagement, activation, and subsequent lysis of tumor cells
Safety	Avoidance of immune-related adverse events such as cytokine release syndrome and immune effector cell-associated neurotoxicity syndrome
Persistence	Proliferation, expansion, or presence of CAR-T cells
Exhaustion	Chronic antigen stimulation of T cells resulting in upregulation of co-inhibitory molecules

CAR-T therapies must be optimized for persistence and efficacy while minimizing T cell exhaustion and adverse events (AE) that can result from excessive immune system stimulation (Figure 1). Recent studies have identified biomarkers associated with all of these key attributes. For example, the persistence of CAR-T cells is required to promote effective anti-tumor responses and lasting remission, and can be monitored by measuring cytokines such as IL-7, IL-15, IL-18, and IFN- γ . These markers correlate well with the number of CAR-T cells present in circulation.^{4,6-8} CAR-T persistence is associated with a durable increase in T cell activity and secretion of proteins including IFN- γ , Perforin, and Granzyme B.⁶⁻⁸

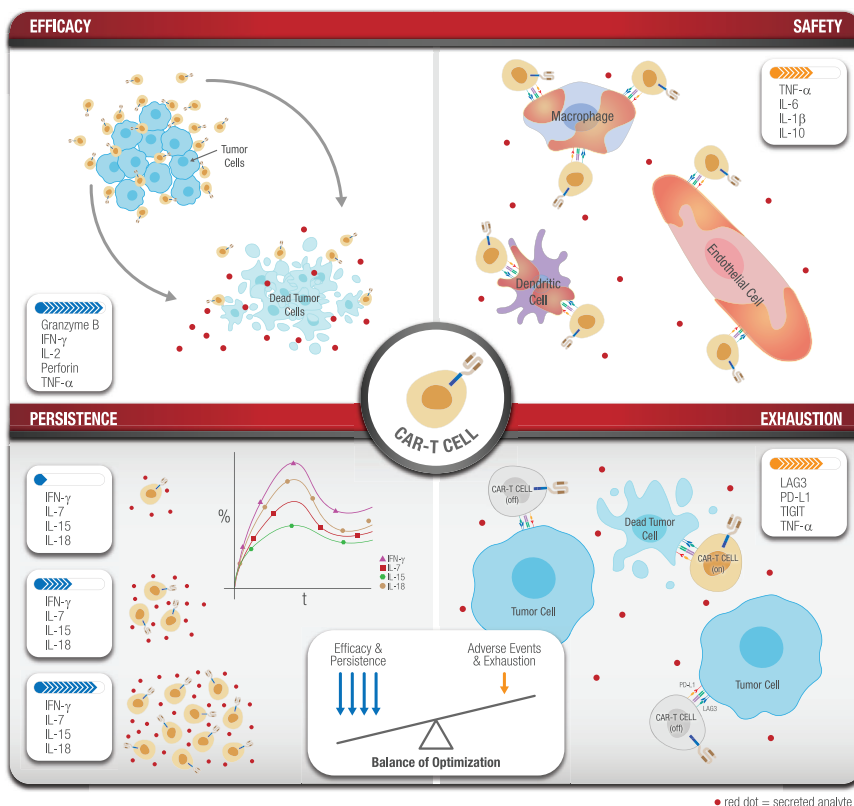


Figure 1. Efficacy biomarkers are surrogate indicators of tumor lysis and T cell function. Safety biomarkers are surrogates of AEs of CAR-T, such as CRS, ICANS, or on-target/off-tumor effects. Persistence and exhaustion biomarkers provide surrogate measures of T cell activity.

In parallel with efficacy, safety must also be tuned. Several markers associated with pro-inflammatory responses, including IL-6, IL-1 β , and MCP-1 are commonly secreted following CAR-T treatment in subjects with Grade 2 or Grade 3 AEs or greater tumor burden.⁸⁻¹⁰ A balance between the safety and efficacy of CAR-T therapy is desirable and specific for each patient. In a retrospective phase 2 study by Neelapu and colleagues⁹, a biomarker signature associated with CRS and ICANS in patients dosed with axi-cel (Yescarta) was found. Specifically, peak levels of cytokines such as IL-6, GM-CSF, IFN- γ , and others were detected in the serum of patients within 8 days of dosing. The presence and magnitude were at least two times greater in patients who experienced AEs \geq Grade 3 than those who did not. Interestingly, this study demonstrated an overall complete response rate of 78%, indicating that CAR-T therapy removed the tumor burden despite most patients experiencing AEs such as CRS. These findings indicate that efficacy and safety may be counterparts in determining response rates in current CAR-T clinical trials.

Lastly, T cell exhaustion limits the expansion, persistence, and efficacy of CAR-T products. Surface and shed biomarkers such as CTLA-4, PD-L1, and TIGIT are common inhibitory signals linked to the exhaustion of T cells following chronic antigen stimulation.¹¹ Patients in clinical trials with upregulation of these biomarkers typically relapse.¹² Deletion of the extracellular domain of surface markers such as PD-L1 or CTLA-4 on CAR-T cells has limited the exhaustion profile in preclinical models with increased persistence and expansion of CAR-T and improved efficacy,¹³ indicating that monitoring and appropriately modifying the markers associated with exhaustion may improve clinical outcomes.

Product Design and Characterization

Four U-PLEX Combos (multiplex assays) were designed to support research regarding the functions and mechanisms associated with CAR-T cell therapies. The sensitivity and specificity of individual U-PLEX assays are characterized using assay-specific reagents as well as serum, plasma, and cell culture supernatant samples. The four Combos noted in Table 2 combine multiple assays for translationally relevant biomarkers. Like all Combos, these assays were characterized for non-specific interactions between each assay at the component level. Non-specific binding (NSB) for each assay was calculated as a percentage of non-specific signal to specific signal. The observed NSB was less than 0.2% for all assays across all Combos, with the exception of CD28 and LAG3 in the CAR-T Exhaustion Combo. That pair-wise %NSB was 2.27%.

For further characterization, peripheral blood mononuclear cells (PBMC) from a healthy donor were cultured and stimulated using various conditions for 24-72 hours. Cell culture supernatants were collected and the Efficacy (Fig. 2A), Persistence (Fig. 2B), and Exhaustion (Fig. 2C) Combos were evaluated for their functional readouts. Stimulation with various Toll-like receptor (TLR) and T cell agonists resulted in the upregulation of interferon and checkpoint molecule secretion from immune cells over multiple logs of concentration range. These data emphasize the sensitivity and dynamic range integral to MSD's U-PLEX platform.

Potential research applications for these Combos include:

- Examination of CAR-T cell target-specific engagement and tumor cell killing in relevant human cell culture models.
- Elucidation of persistence and exhaustion biomarker profiles of CAR-T cell products in development.
- Linking exploratory biomarker profiles to adverse event findings.
- Rapid measurement of biomarkers commonly associated with efficacy in CAR-T therapy in relevant humanized animal models and translational studies.
- Justification of lymphodepletion regimens in pre-clinical applications.

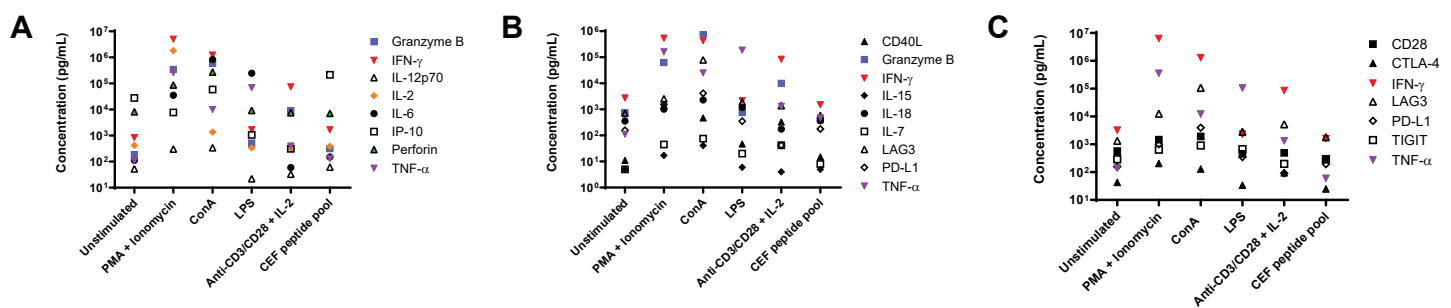


Figure 2. The concentration of analytes in the Efficacy (A), Persistence (B), and Exhaustion (C) U-PLEX Combos were tested from cell culture supernatants that were diluted 40-fold. PBMCs from an individual donor were cultured and stimulated under various conditions for 24-72 hours. Concentrations of each analyte are reported for each stimulation. (PMA = phorbol-12-myristate-13-acetate; LPS = lipopolysaccharide; IL = interleukin; ConA = Concanavalin A; CEF = cytomegalovirus (C), Epstein-Barr virus (E), and influenza virus (F)).

Conclusion

These assays are designed to simultaneously measure multiple biomarkers relevant to CAR-T cell research and development. They perform robustly with minimal non-specific binding with broad dynamic range and sensitivity.

Table 2. List of biomarkers in U-PLEX CAR-T cell Combos and their relevance to CAR-T cell research.

Panel	Assay	Description
Persistence	CD40L	Co-stimulatory molecule to prime and expand CAR-T cells
	Granzyme B	Effector molecules secreted by T cells upon target engagement, induces apoptosis in target cells
	IFN- γ	Secreted by proliferating CAR-T cells
	IL-7	Promotes CAR-T cell survival through up-regulation of anti-apoptotic genes
	IL-15	Regulates CAR-T cell effector function
	IL-18	Enhances the proliferative capacity of CAR-T cells
	LAG3	Inhibitory ligand resulting in reduced CAR-T cell activation and regulation of immune cell homeostasis
	PD-L1	Promotes T cell exhaustion and self-tolerance
	TNF- α	Pro-inflammatory cytokine secreted by numerous cells upon inflammatory stimuli, promoting activation and proliferation of naïve and effector CAR-T cells
Efficacy	Granzyme B	Effector molecule secreted by T cells upon target engagement, induces apoptosis in target cells
	IFN- γ	Secreted by proliferating CAR-T cells
	IL-2	Promotes CAR-T cell proliferation and effector function
	IL-6	Pro-inflammatory cytokine secreted by macrophages in response to microbial molecules following engagement with T cells
	IL-12p70	Stimulates CAR-T cells to improve cytotoxic capacity
	IP-10	Stimulates chemotaxis of CAR-T cells to tumors
	Perforin	Pore-forming cytolytic protein expressed in granules of cytotoxic CAR-T cells
	TNF- α	Pro-inflammatory cytokine secreted by numerous cells upon inflammatory stimuli promoting activation and proliferation of naïve and effector CAR-T cells
Exhaustion	CD28	Activates CAR-T cells by increasing their sensitivity to the antigen of interest
	CTLA-4	Inhibits CAR-T activation and upregulates co-inhibitory molecules
	IFN- γ	Secreted by proliferating CAR-T cells
	LAG3	Inhibitory ligand resulting in reduced CAR-T cell activation and regulation of immune cell homeostasis
	PD-L1	Promotes CAR-T cell exhaustion and self-tolerance
	TIGIT	Co-inhibitory ligand for T cells impacting CAR-T function
	TNF- α	Pro-inflammatory cytokine secreted by numerous cells upon inflammatory stimuli promoting activation and proliferation of naïve and effector CAR-T cells
Safety	GM-CSF	Affects myeloid cell proliferation and activation, activates CD4+ T cells
	IFN- γ	Secreted by proliferating CAR-T cells
	IL-1 β	Pro-inflammatory cytokine that is secreted following CAR-T engagement with endothelial cells
	IL-6	Pro-inflammatory cytokine secreted by macrophages in response to stimulus following engagement with CAR-T cells
	IL-10	Inhibits tumor-specific function while upregulating pro-inflammatory effects following on-target, off-tumor CAR-T cell engagement
	MCP-1	Enhances macrophage responses to pro-inflammatory stimuli following CAR-T cell engagement
	TNF- α	Pro-inflammatory cytokine secreted by numerous cells upon inflammatory stimuli promoting activation and proliferation of naïve and effector CAR-T cells

References

1. Siegler EL, Kenderian SS. Neurotoxicity and cytokine release syndrome after chimeric antigen receptor T cell therapy: Insights into mechanisms and novel therapies. *Front Immunol.* 2020 Aug 28;11:1973.
2. Sievers S, Watson G, Johny S, Adkins S. Recognizing and grading CAR T-cell toxicities: An advanced practitioner perspective. *Front Oncol.* 2020; 10:885.
3. Sheth S, Gauthier J. Taming the beast: CRS and ICANS after T-cell therapy for ALL. *Bone Marrow Transplantation.* 2021; 56:552-566.
4. Pietrobon V, Todd LA, Goswami A, Stefanson O, Yang Z, Marincola F. Improving CAR T-Cell persistence. *Int J Mol Sci.* 2021; 22(19):10828.
5. Gumber D, Wang LD. Improving CAR-T immunotherapy: Overcoming the challenges of T cell exhaustion. *eBioMedicine.* 2022; 77:103941.
6. Guedan S, Posey Jr AD, Shaw C, Wing A, Da T, Patel PR, McGettigan SE, Casado-Medrano V, Kawalekar OU, Uribe-Herranz M, Song D, Melenhorst JJ, Lacey SF, Scholler J, Keith B, Young RM, June CH. Enhancing CAR T cell persistence through ICOS and 4-1BB costimulation. *JCI Insight.* 2018; 3(1): e96976.
7. Locke FL, Neelapu SS, Bartlett NL, Siddiqi T, Chavez JC, Hosing CM, Ghonadi A, Budde LE, Bot A, Rossi JM, Jiang Y, Xue AX, Elias M, Aycock J, Wiezorek J, Go WY. Phase 1 Results of Zuma-1: A multicenter study of KTE-C19 Anti-CD19 CAR T cell therapy in refractory aggressive lymphoma. *Mol Ther.* 2017;25(1):285-295.
8. Kockenderfer JN, Somerville RP, Lu T, Yang JC, Sherry RM, Feldman SAA, McIntyre L, Bot A, Rossi J, Lam N, Rosenberg SA. Long-duration complete remissions of diffuse large B cell lymphoma after Anti-CD19 chimeric antigen receptor T cell therapy. *Mol Ther.* 2017; 25(10):2245-2253.
9. Neelapu SS, Dickinson M, Munoz J, Ulrichson ML, Thieblemont C, Oluwole OO, Herrera AF, Ujjani CS, Lin Y, Riedell PA, Kekre N, de Vos S, Lui C, Milletti F, Dong J, Xu H, Chavez JC. Axicabtagene ciloleucel as first-line therapy in high-risk large B-cell lymphoma: the phase 2 ZUMA-12 trial. *Nat Med.* 2022; 28:735-742.
10. Locke FL, Rosi JM, Neelapu SS, Jacobson CA, Miklos DB, Ghobadi A, Oluwole OO, Reagan PM, Lekakis LJ, Lin Y, Sherman M, Better M, Go WY, Wiezorek JS, Wue A, Bot A. Tumor burden, inflammation, and product attributes determine outcomes of axicabtagene ciloleucel in large B-cell lymphoma. *Blood Adv.* 2020; 4(19):4898-4911.
11. Wherry EJ. T Cell Exhaustion. *Nat Rev Immunol.* 2011;12(6): 492-499.
12. Gargett T, Yu W, Dotti G, Yvon ES, Christo SN, Hayball JD, Lewis ID, Brenner MK, Brown MP. GD2-specific CAR T cells undergo potent activation and deletion following antigen encounter but can be protected from activation-induced cell death by PD-1 blockade. *Mol Ther.* 2016;24(6):1135-1149.
13. Agarwal S, Aznar MA, Rech AJ, Good CR, Kuramitsu S, Da T, Gohil M, Chen L, Hong S-J A, Ravikumar P, Rennels AK, Salas-Mckee J, Kong W, Ruella M, Davis MM, Piesa G, Fraietta JA, Porter DL, Young RM, June CH. Deletion of the inhibitory co-receptor CTLA-4 enhances and invigorates chimeric antigen receptor T cells. *Immunity.* 2023;56(10):2388-2407.

MESO SCALE DISCOVERY®

1601 Research Boulevard

Rockville, Maryland 20850-3173

Phone: 1-240-314-2600 Fax: 1-240-632-2219

For Research Use Only. Not for use in diagnostic procedures.

MESO SCALE DISCOVERY, Meso Scale Diagnostics, MSD, mesoscale.com, www.mesoscale.com, methodicalmind.com, www.methodicalmind.com, DISCOVERY WORKBENCH, InstrumentLink, MESO, MesoSphere, Methodical Mind, MSD GOLD, MULTI-ARRAY, MULTI-SPOT, QuickPlex, ProductLink, SECTOR, SECTOR HTS, SECTOR PR, SULFO-TAG, TeamLink, TrueSensitivity, TURBO-BOOST, TURBO-TAG, N-PLEX, R-PLEX, S-PLEX, T-PLEX, U-PLEX, V-PLEX, MSD (design), MSD (luminous design), Methodical Mind (head logo), 96 WELL SMALL-SPOT (design), 96 WELL 1-, 4-, 7-, 9-, & 10-SPOT (designs), 384 WELL 1- & 4-SPOT (designs), N-PLEX (design), R-PLEX (design), S-PLEX (design), T-PLEX (design), U-PLEX (design), V-PLEX (design), It's All About U, Spot the Difference, The Biomarker Company, and The Methodical Mind Experience are trademarks and/or service marks owned by or licensed to Meso Scale Diagnostics, LLC. All other trademarks and service marks are the property of their respective owners. ©2024 Meso Scale Diagnostics, LLC. All rights reserved.

