

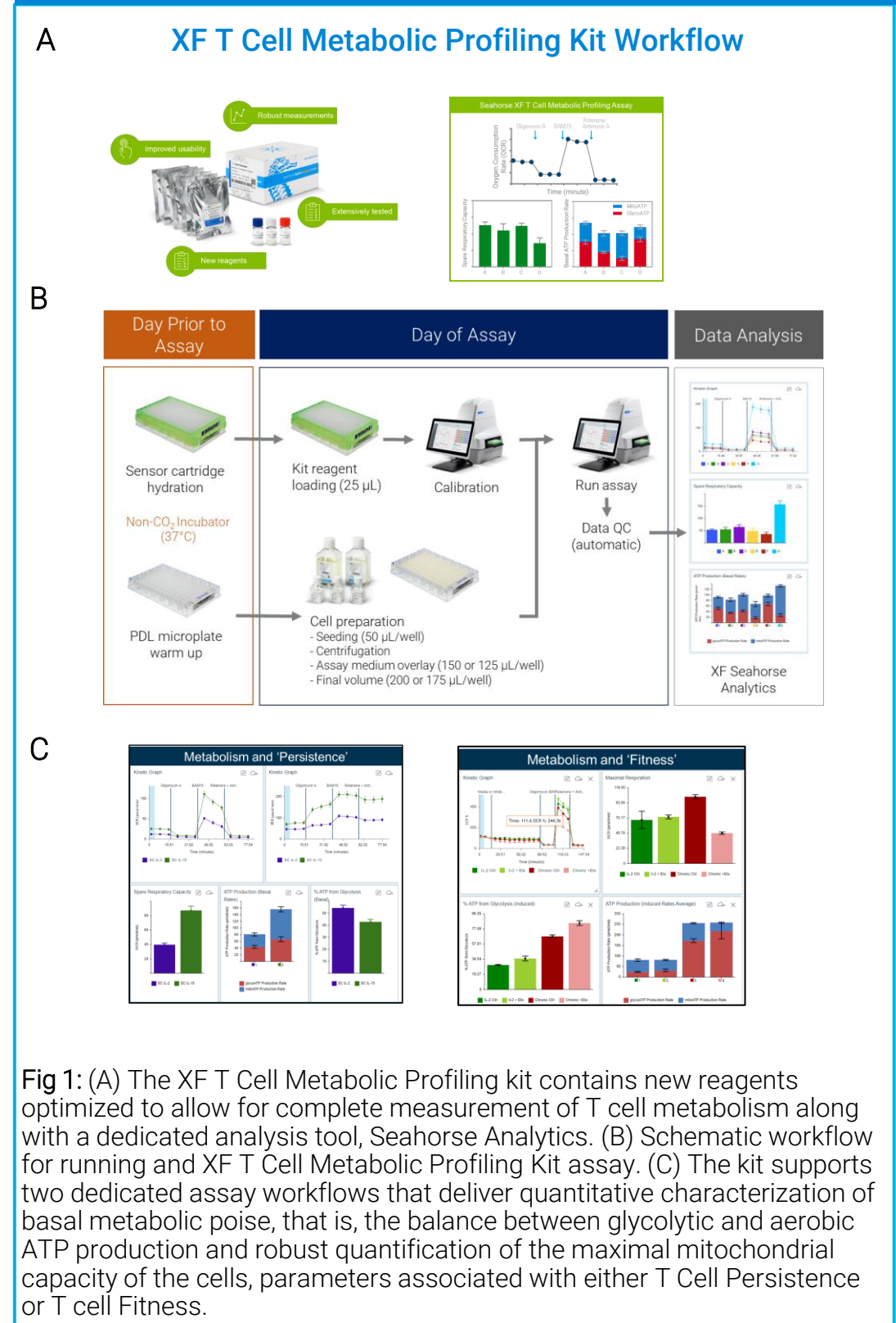
Introduction

Cellular metabolism is a known driver of T cell fate and function. For this reason, it represents a key attribute to be assessed during T cell therapy optimization and production.

Agilent Seahorse XF T Cell Metabolic Profiling kit is a robust new solution that allows for the simultaneous measurement of basal metabolic requirements, metabolic poise, and mitochondrial bioenergetic capacity in T cells. These parameters have previously been shown to correlate with increased T cell persistence and improved metabolic fitness.

In this study, we used the assay for comprehensive assessment of the bioenergetic profile at different time points of human T cells when expanded in different cell culture media and cytokine formulations that were reported to increase T cell persistence and improved metabolic fitness.

Experimental



Results and Discussion

BAM15 is an Improved Uncoupler for T cells

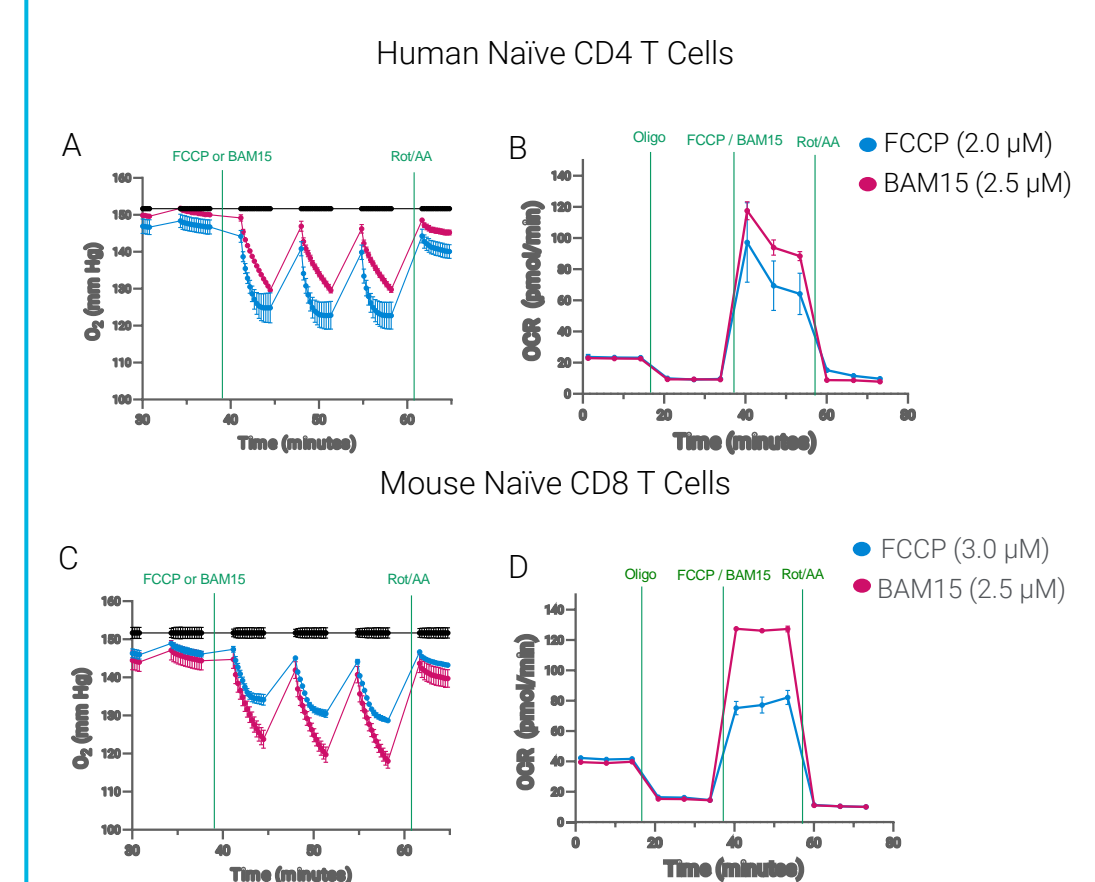


Fig 2: BAM15 addition results in a more robust uncoupling response in T cells compared with FCCP. BAM15 or FCCP were titrated in either human naïve CD4 T cells or mouse naïve CD8 T cells. (A), (C): Changes in extracellular oxygen levels after uncoupler addition, highlighting the more consistent rate during the 3 minutes of instrument measurement obtained with FCCP. (B), (D): OCR kinetic profile in naïve human CD4 T cells, and mouse naïve CD8 T cells, respectively illustrating underestimation of Max respiration obtained when FCCP is used as uncoupler in the assay.

BAM15 Results in a Broader Effective Range and Simplified Concentration Optimization

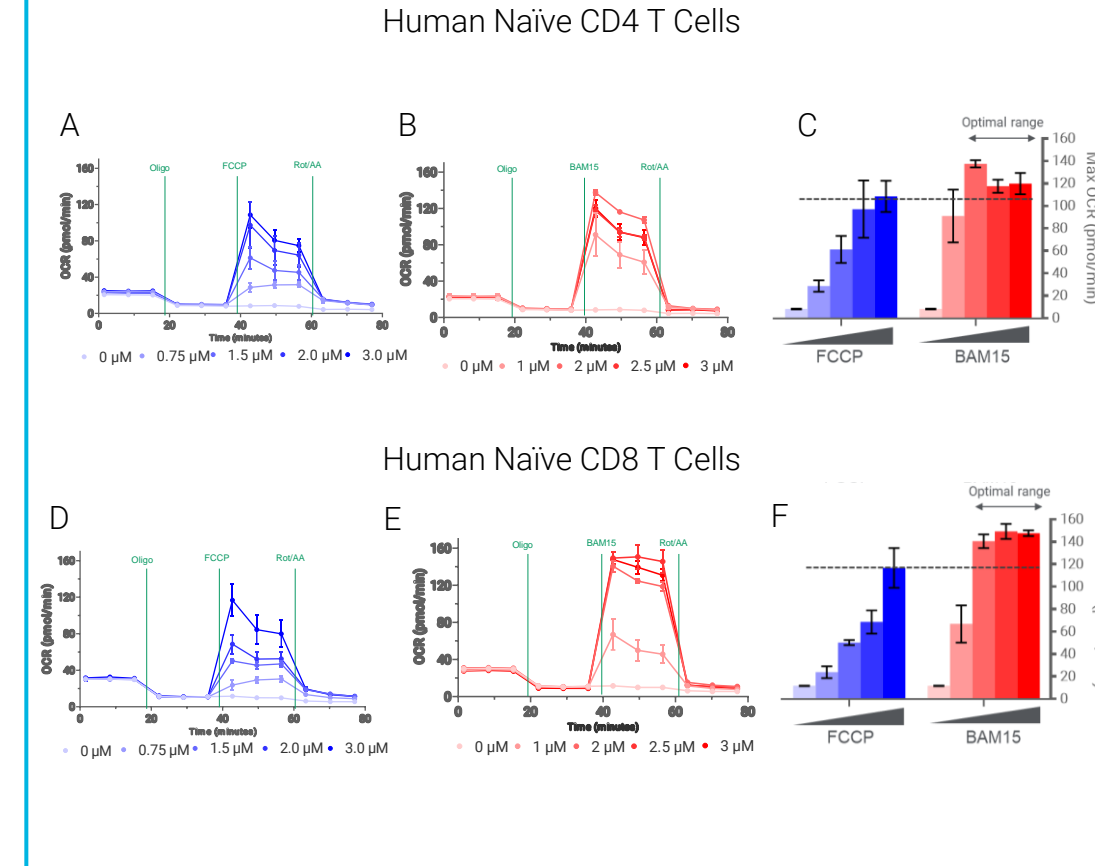
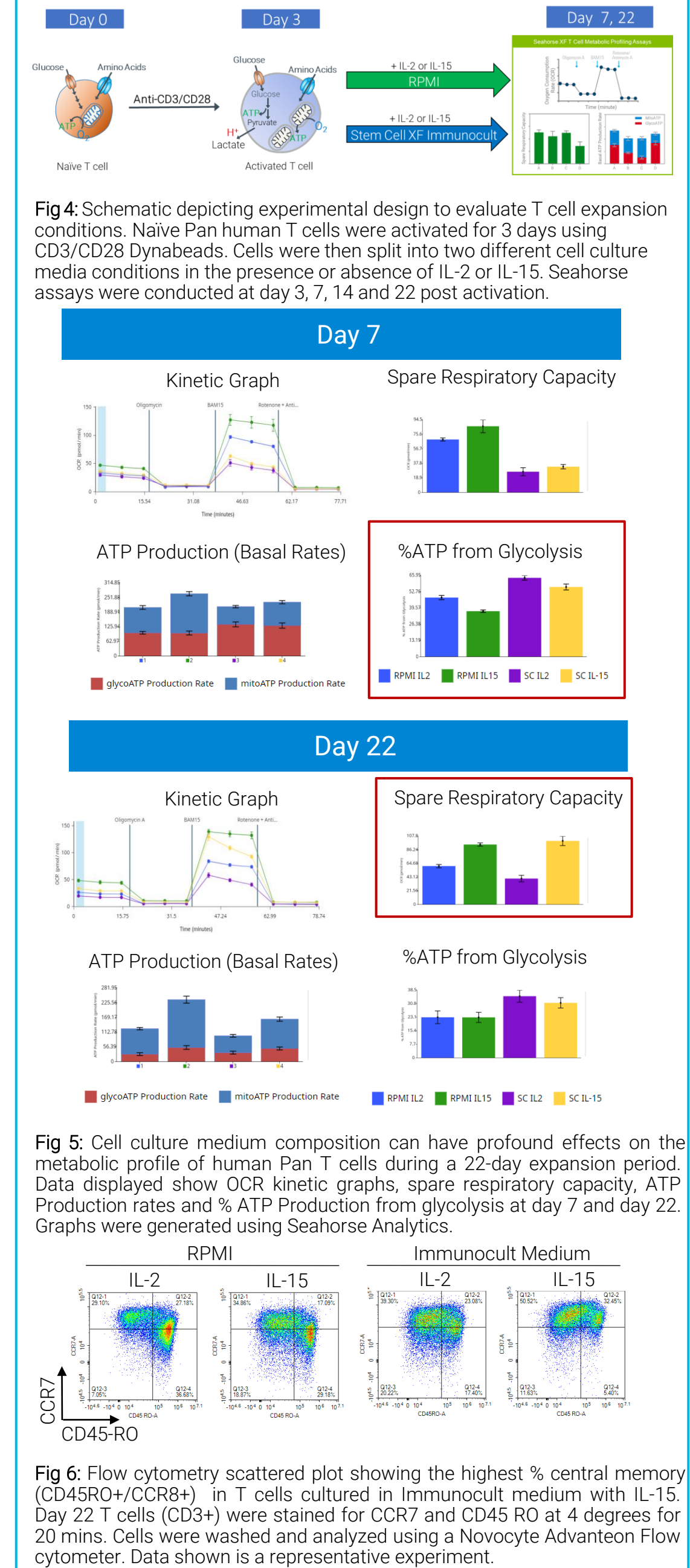


Fig 3: BAM15 at 2.5 µM shows the optimal response in naïve T cells, minimizing requirement of concentration optimization for each cell sample: (A) Kinetic OCR trace of FCCP titration (0 – 3 µM) in naïve CD4 T cells. (B) Kinetic OCR trace of BAM15 titration (0 – 3 µM) in naïve CD4 T cells. (C) Bar graph comparing maximum OCR obtained with FCCP or BAM15 titration. (D) Kinetic OCR trace of FCCP titration (0 – 3 µM) (E) Kinetic OCR trace of BAM15 titration (0 – 3 µM) in naïve CD8 T cells. (F) Bar graph showing maximum OCR obtained with FCCP or BAM15 titration.

Metabolic evaluation of T cell expansion conditions



Conclusions

- BAM15 demonstrates significant benefit over the traditional uncoupler, FCCP, when assessing T cell oxidative metabolism. It supports more robust oxygen consumption rates over measurement time and allows for accurate measurement of ATP production rates.
- Higher spare respiratory capacity has often been associated with better T cell persistence *in vivo*. At day 7 post-activation, IL-15 cultured cells have a lower % of ATP production from glycolysis, indicating a more aerobic poise. By day 22, this results in IL-15 cultured cells having significantly higher spare respiratory capacity than IL-2 cultured cells. This parameter may offer unique early insight into T cell metabolism during expansion.
- The Agilent XF T Cell Metabolic Profiling kit allows for simultaneous evaluation of glycolytic and mitochondrial activity and capacity, providing a robust and simple assay for comprehensive profiling of T cell populations.
- The data presented show the impact of cell culture conditions during expansion on the resulting T cell profile and fate, highlighting the importance of including metabolic profiling to monitor and tune culture conditions during T cell therapy development.

Methods

Cell Culture
Human T cells (CD4 and Pan T cells) were obtained from Stem Cell Technologies. For naïve T cell experiments, T cells were thawed and rested overnight in Stem Cell XF Immunocult Medium. For T cell expansion experiments, Pan T cells were rested overnight in Stem Cell Technologies Immunocult XF medium. On day 1 cells were activated using CD3/CD28 dynabeads (Thermo Fisher). After 3 days, dynabeads were magnetically removed and cells were put into either RPMI + 10% FBS or Immunocult XF culture medium with either IL-2 or IL-15 and maintained over 22 days. Mouse CD8 T cells were magnetically purified from splenocytes obtained at Hooke Laboratories and XF assays were run the same day.

Seahorse XF Assays
Naïve human T cells, expanded T cells and mouse CD8 T cells were resuspended in Seahorse XF RPMI, pH 7.4 medium and seeded at 200,000 cells, 100,000 cells and 150,000 cells per well respectively in PDL-coated microplates. Assays were carried out using a Seahorse XF Pro. For detailed information see XF T Cell Metabolic Profiling Kit user guide at Agilent.com

References

Select publications relating to the utility of T cell metabolic measurements for T cell therapy

Alizadeh et al. (2019). IL15 Enhances CAR-T Cell Antitumor Activity by Reducing mTORC1 Activity and Preserving Their Stem Cell Memory Phenotype. *Cancer Immunol Res* 7(5): 759-766. DOI: 10.1158/2326-6066.CCR-18-0466

Buck et al. (2017). Metabolic Instruction of Immunity. *Cell* 169(4): 570-86. doi: 10.1016/j.cell.2017.04.004

Rostamian et al. (2021). A metabolic switch to memory CAR T cells; Implications for cancer treatment *Cancer Letters* 500, 107–118. <https://doi.org/10.1016/j.canlet.2020.12.004>

Schärping et al. (2021). Mitochondrial stress induced by continuous stimulation under hypoxia rapidly drives T cell exhaustion. *Nature Immunology* 22(2): 205-215. doi: 10.1038/s41590-020-00834-9

Song et al. (2018). IRE1α-XBP1 controls T cell function in ovarian cancer by regulating mitochondrial activity *Nature* 562(7727): 423–428. <https://doi.org/10.1038/s41586-018-0597-x>

J. Walls, N. Romero (2022). Assessing T Cell Bioenergetic Poise and Spare Respiratory Capacity Using Extracellular Flux Analysis. Application Note, Agilent.com. 5994-4494EN

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