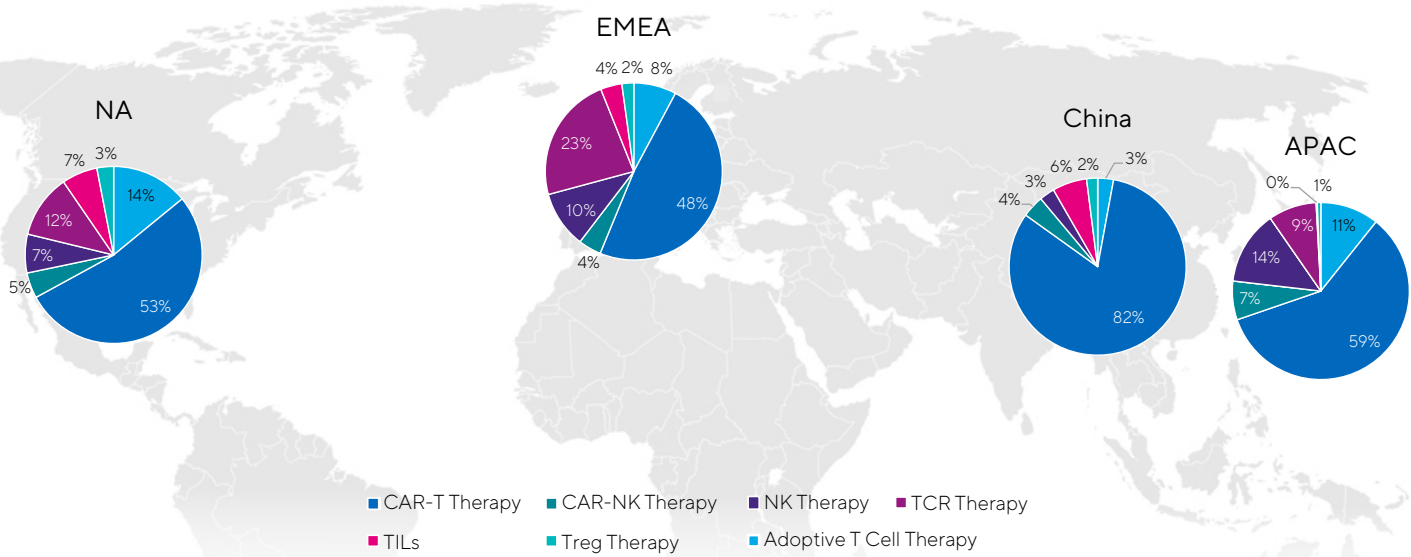


Enhancing Cell Growth With Cytokines

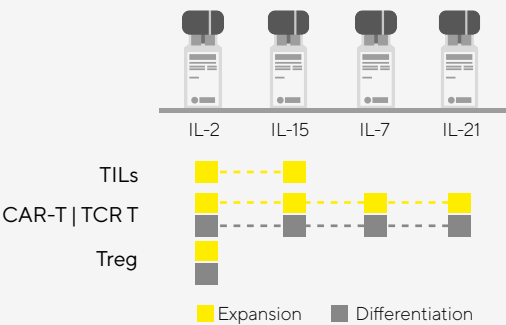
With nearly 4,800 immune cell therapy products in development today, scientists around the world depend on cytokines to support the culture of a variety of immune cell types.



A breakdown of projects per immune cell type globally. Source: Global Trends, February 2022

Cytokine Combinations to Tailor Your Immune Cell Culture

Tune cytokines to help promote successful expansion and differentiation of a variety of immune cells.



Combining CellGenix® cytokines makes T cell production more efficient and robust while reducing handling steps.

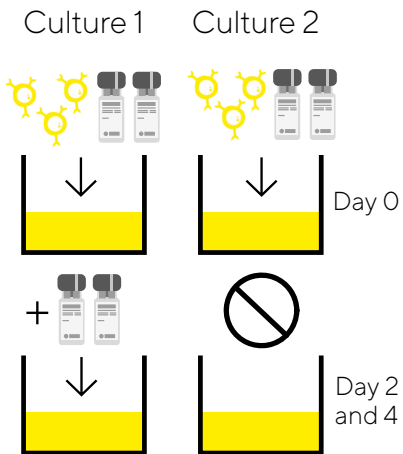


Data Points:

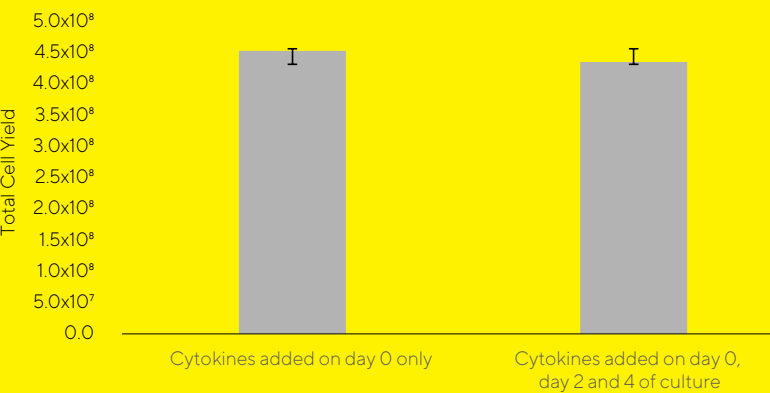
1. CellGenix® IL-7, CellGenix® IL-15, along with anti-CD3/anti-CD28-coupled magnetics beads were added to pre-warmed media containing 5×10^6 purified CD3+ T cells/well.
2. The media cultured over an 11-day expansion, with one culture receiving CellGenix® IL-7 and CellGenix® IL-15 on days 2 and 4.

Results:

At day 11 of culture, no significant differences in expansion and viability were observed between the two conditions. This demonstrates that **CellGenix® IL-7 and IL-15 support T cell culture over 11 days** — meaning you could **reduce open handling steps and save up to 15% of aliquot volume**.



Day 11 Cell Yield



Day 11 Cell Viability

