

CTS™ Dynabeads™ CD3/CD28 enables simultaneous one-step naïve and early memory T cell isolation and activation

Tor Espen Stav-Noraas¹, Reidun Hartberg¹, Anette Kullmann¹, Katrine Alfnes¹, Øystein Åmellem¹, Lauren Miller², Tim Looney², Geoffrey M. Lowman² and Hilde Almåsbak¹

¹Thermo Fisher Scientific, Ullernchaussen 52, Oslo, Norway ²5781 Van Allen Way, Carlsbad, CA 92008 USA

ABSTRACT

CTS™ Dynabeads™ CD3/CD28 has been used successfully in advancing T cell therapies through >88 clinical trials and lately also in commercial T cell drug manufacturing.

CTS™ Dynabeads™ CD3/CD28 requires no upstream T cell selection as this technology simultaneously isolates and activates naïve and early memory T cells based on CD3 and CD28 co-expression. T cells are isolated with high recovery (>90%) and purity (>95%) and uniformly stimulated (>95% CD25⁺ day 3 post-activation). Activated and expanded T cells preserve a young phenotype (CD28⁺CD62L⁺) with low PD-1 expression day 7-10, as well as demonstrate increased clonal diversity compared to starting T cells.

Early removal of Dynabeads™ (days 2, 3, and 5 post-activation) is feasible but comes with some cell loss days 2 and 3 (T cell recovery 50-80%). Nearly all T cells are recovered following bead removal day 5. Transduction efficiency was higher when beads were removed and cells transduced day 2 compared to day 3 post-activation.

MATERIALS AND METHODS

- T cells from healthy donor peripheral blood mononuclear cells (PBMCs) were isolated and activated using CTS™ Dynabeads™ CD3/CD28 (bead: cell ratio 3:1) (Thermo Fisher Scientific) in Permalife cell culture bags (Origen) with CTS™ OptiMizer™ T Cell Expansion SFM supplemented with CTS™ Immune Cell Serum Replacement, L-Glutamine and Gentamicin (incomplete media) (all Thermo Fisher Scientific).
- Post bead removal T cells were expanded in G-rex plates (Wilson Wolf) or Permalife cell culture bags (Origen)
- Isolated and activated T cells were expanded in complete media (incomplete media + 100 U/mL IL-2) for 7-10 days. Dynabeads were magnetically removed at day 3 post stimulation or day 2 and 5 as described (Figure 6).
- T cells were transduced with γ-retrovirus encoding EGFP (BionTech AGH) in retronectin (TaKaRa) coated corning microplates (Sigma-Aldrich) post bead removal.
- Flow cytometry analysis of T cells were performed on a BD Fortessa (BD bioscience) instrument using FlowJo
- RNA from T cells was isolated using Dynabeads™ mRNA DIRECT™ Purification Kit (Thermo Fisher Scientific) and analyzed with AmpliSeq™ TCRβ sequencing (Thermo Fisher Scientific)

TYPICAL T CELL MANUFACTURING WORKFLOW

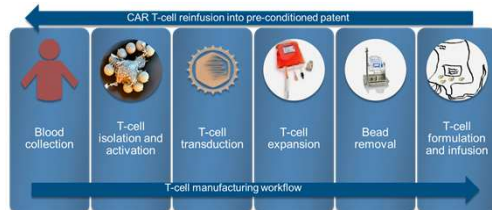


Figure 1: Example of manufacturing and delivery pipeline of CAR T cell therapies (modified from ref 1). PBMCs are harvested from the patient (or a T cell donor) and transferred to a GMP facility where the T cells are isolated and activated in the presence of CTS™ Dynabeads™ CD3/CD28 and genetically engineered by viral transduction to express the CAR. The activated T cells are expanded in CTS™ OptiMizer™ T Cell Expansion SFM supplemented with CTS™ Immune Cell Serum Replacement for a period, typically 7-10 days, to reach a therapeutic relevant number before magnetic bead removal and drug formulation, either for freezing or for adoptive transfer.

RESULTS

Nearly 100% recovery of naïve and early memory T cells with one-step simultaneous T cell isolation and activation

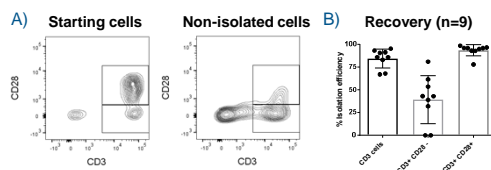


Figure 2. CTS™ Dynabeads™ CD3/CD28 preferentially isolates CD3⁺ CD28⁺ positive cells. PBMCs were incubated for 30 min with CTS™ Dynabeads™ CD3/CD28 at a ratio of 3:1 beads:cells and the negative (non isolated) fraction was analyzed and used to calculate isolation efficiency. Representative cytogram demonstrating efficiency of CD3⁺CD28⁺ T cells isolation (A). Recovery (n=9) experiments (B)

One-step isolation and activation yields pure and uniformly activated T cells

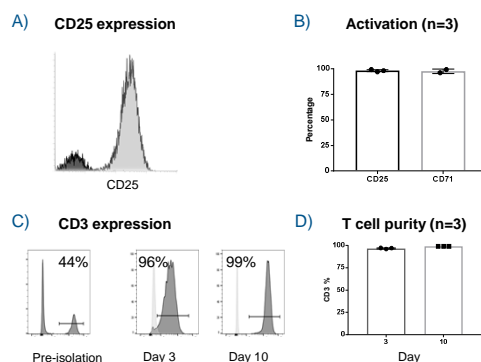


Figure 3. CTS™ Dynabeads™ CD3/CD28 isolated T cells are pure and uniformly stimulated. Isolated T cells were expanded for 3 days before analyzing activation markers (A,B). Overlay histogram of activated T cells and a non-activated control (A). T cell purity was analyzed at days 0 (starting material), 3 and 10 post-activation (C, D).

CTS™ Dynabeads™ CD3/CD28 expanded CD4⁺ and CD8⁺ T cells maintains an early memory phenotype (day 10)

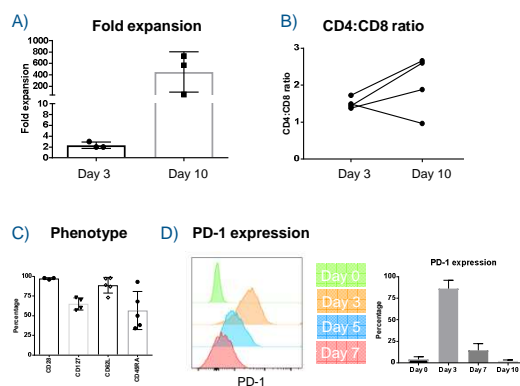


Figure 4. Activated T cells typically expands >100 fold in 10 days. Isolated and activated T cells were expanded for up to 10 days (A) and both CD4 and CD8 T cells proliferate (B). Day 10 post-activation, cells were mainly CD45RA⁺CD127⁺CD28⁺CD62L⁺ (C) with low PD-1 expression (D).

Dynabeads expanded T cells have increased clonal diversity

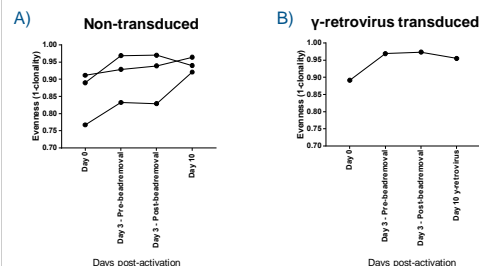


Figure 5. Increased clonal diversity after T cell expansion. T cells were isolated and stimulated with CTS™ Dynabeads™ CD3/CD28 (3 donors). At days 0, 3 and 10, samples were harvested for TCR-sequencing (A). In one experiment, T cells were also transduced with γ-retrovirus day 3 post stimulation (B).

Early bead removal is feasible and day 2 transduction improves γ-retroviral transduction

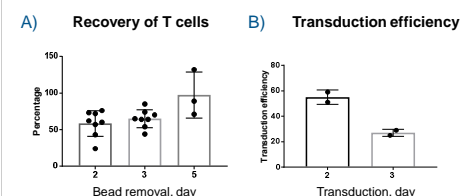


Figure 6. Shorter stimulation of T cells improves γ-retrovirus transduction efficiency. T cells were isolated with CTS™ Dynabeads™ CD3/CD28 in Origen cell culturing bags and stimulated for 2 or 3 days. The beads were magnetically removed day 2 or 3 before transducing the cells with γ-retrovirus encoding an EGFP reporter in retronectin coated wells (2.5 MO). The recovery of T cells were lower after early bead removal (A) compared to 5 days stimulation. Transduction efficiency was assessed three days post-transduction (B)

CONCLUSIONS

- CTS Dynabeads CD3/CD28 isolates CD3⁺CD28⁺ T cells with high efficiency and purity
- Isolated & activated CD4 and CD8 T cells expand 100-800 fold in 10 days while retaining a young phenotype (CD45RA⁺CD62L⁺CD127⁺PD-1⁻)
- Expanded T cells have increased clonal diversity (TCR sequencing)
- Bead removal day 2 improves transduction efficiency of γ-retroviral vector

REFERENCE

(1) Modified from Journal of Immunology Research (2016), CAR T Cell Therapy: A Game Changer in Cancer Treatment. Almåsbak, Aarvak, and Vemuri

TRADEMARKS/LICENSING

© 2018 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

For Research Use or Manufacturing of Cell, Gene, or Tissue-Based Products. CAUTION: Not intended for direct administration into humans or animals.

ThermoFisher
SCIENTIFIC