

A Sendai viral reprogramming kit for completely xeno-free generation of iPSC

Chad C. MacArthur and Uma Lakshmiopathy

Thermo Fisher Scientific, Cell Biology, 5781 Van Allen Way, Carlsbad, CA, USA 92008

ABSTRACT

The induced pluripotent stem cell (iPSC) field has grown immensely in the past ten years, starting with basic research, and now moving more toward clinical applications. As iPSC move towards the clinic, it is vital that high quality cells are used. These cells should be free of reprogramming transgenes, free of adventitious agents, genetically stable, and functionally pluripotent. Typical reprogramming workflows often include animal origin components, such as fetal bovine serum (FBS) or bovine serum albumin (BSA), which can be potential sources of adventitious agents. iPSC that are generated in a xeno-free, clinically relevant workflow will be free from any such potential contaminants, and much more readily applicable to clinical uses. Here, we show the use of a Sendai reprogramming kit manufactured under a certified ISO 9001 Quality Management System, which is also free of animal origin components, to generate high quality iPSC. The kit was used to generate iPSC derived from both blood-derived cells, and skin fibroblasts. The reprogramming workflows used were xeno-free, starting from initial cell isolation, all the way through to iPSC expansion and banking. The iPSC generated with this method were shown to be of high quality; they had a normal karyotype, were free of Sendai viral vectors, expressed the standard self-renewal markers, and demonstrated functional pluripotency for all three germ layers. The ability to create high quality iPSC in clinically relevant, xeno-free workflows, will allow researchers to more smoothly transition cells to clinical applications.

INTRODUCTION

For iPSC to be used as starting material for clinical applications, it is important that the iPSC are created using clinically relevant workflows. This includes removing animal origin components, and using reagents which are manufactured reliably. In order to achieve this, the CTS CytoTune-iPS 2.1 Reprogramming Kit was created by applying three major changes to the existing CytoTune-iPS 2.0 Sendai Reprogramming Kit:

- 1.) Removal of animal-origin components (e.g. BSA) from the viral dilution buffer.
 - 2.) Replacing the C-Myc transgene with L-Myc, which is shown to be a less oncogenic member of the same gene family. (1)
 - 3.) Production of the viral vectors under a certified ISO 9001 Quality Management System.
- Given these changes, the new reprogramming kit was used to reprogram both fibroblasts and blood-derived cells in xeno-free workflows, and both reprogramming efficiency and quality of the resulting iPSC clones were examined.

MATERIALS AND METHODS

Fibroblast Reprogramming

Fibroblasts were cultured in either standard serum-based fibroblast medium (DMEM, 10% FBS, and 1% Non-essential amino acids), or xeno-free fibroblast medium, and transduced with the CTS™ CytoTune™-iPS 2.1 Sendai Reprogramming Kit. Virus was removed after 24 hours, and cells were harvested and plated onto rh-Laminin521 (LN521) at 7 days after transduction. At 8 days after transduction, medium was changed to Essential 8™ Medium, and cells were fed daily. At 21 days after transduction, cells were either isolated for clonal expansion, or reprogramming efficiency was determined by using the Vector® Red Alkaline Phosphatase stain.

PBMC Reprogramming

Peripheral Blood Mononuclear Cells (PBMCs) were cultured in StemPro™34 SFM, plus SCF (100ng/ml), FLT3 (100ng/ml), IL-3 (20 ng/ml), IL-6 (20ng/ml) for 4 days after thaw. Cells were then transduced with CTS™ CytoTune™-iPS 2.1 Sendai Reprogramming Kit. Virus was removed after 24 hours, and cells were plated onto Vitronectin (VTN) or rh-Laminin521 (LN521) at 3 days after transduction. At 7 days after transduction, medium was changed to Essential 8™ Medium, and cells were fed daily. At 18-21 days after transduction, cells were either isolated for clonal expansion, or reprogramming efficiency was determined by using the Vector® Red Alkaline Phosphatase stain.

T-Cell Reprogramming

CD3+ cells were purified from whole blood using the Dynabeads™ FlowComp™ Human CD3 Kit, according to manufacturer instructions. Purified CD3+ cells were cultured with Dynabeads™ Human T-Expander CD3/CD28 and 100U/mL recombinant human IL-2 CTS™ OptMizer™ T Cell Expansion SFM for seven days. Cells were then reprogrammed with a protocol modified from Seki et al. (2). Cells were plated onto CD3 antibody coated plates, then transduced with CytoTune. After 48 hours, cells were plated onto either VTN, or rhLaminin-521 in Essential 8™ Medium. At 18 days after transduction, cells were analyzed for AP activity.

RESULTS

Figure 1. CTS CytoTune 2.1 successfully reprograms human dermal fibroblasts

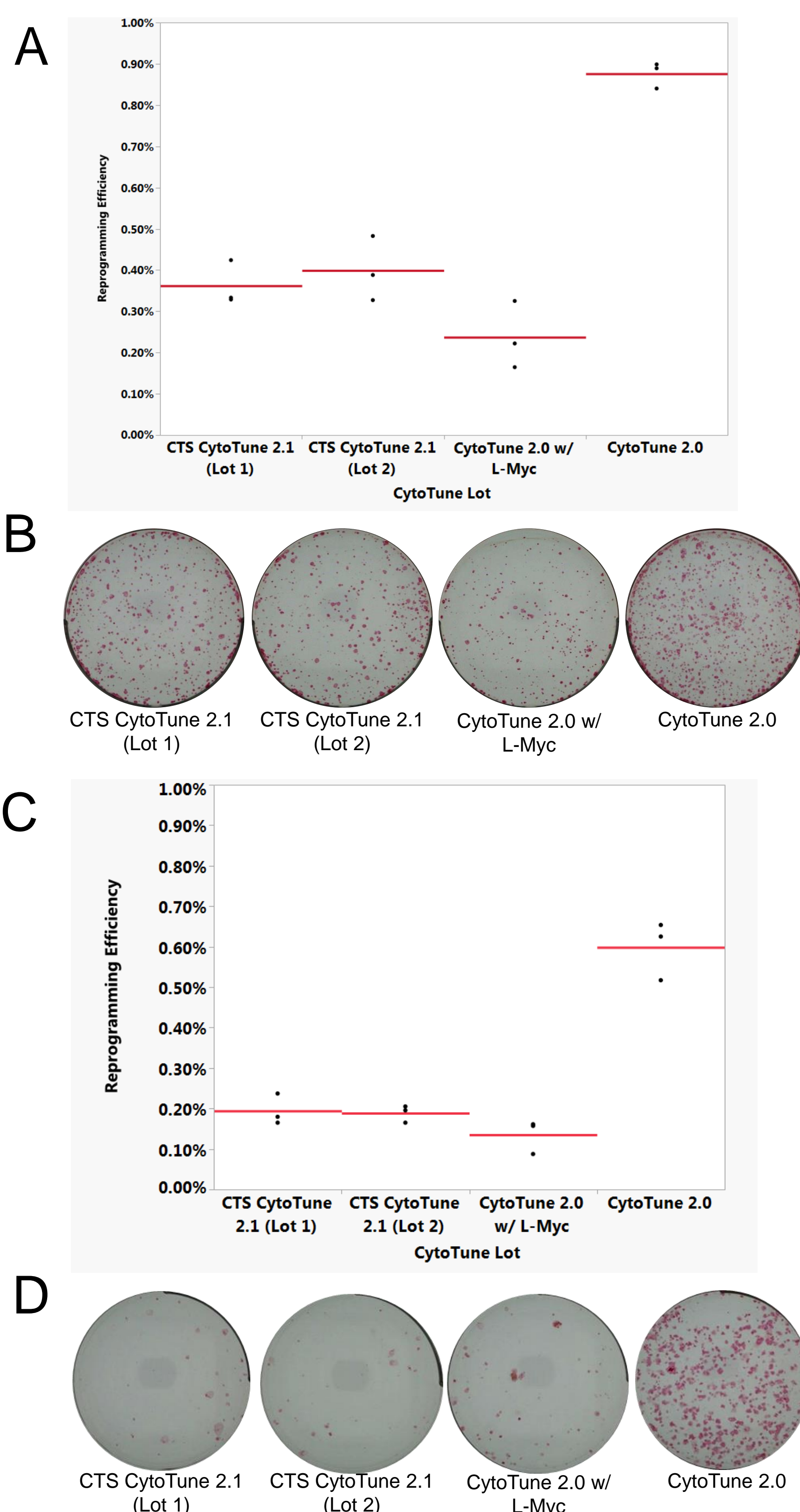
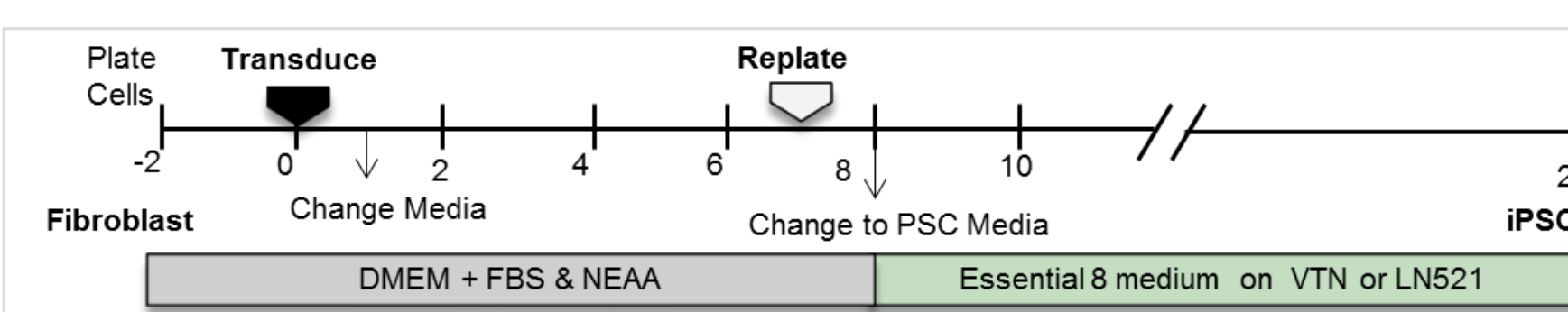


Figure 2. CTS CytoTune 2.1 vectors clears from iPSC clones by Passage 8

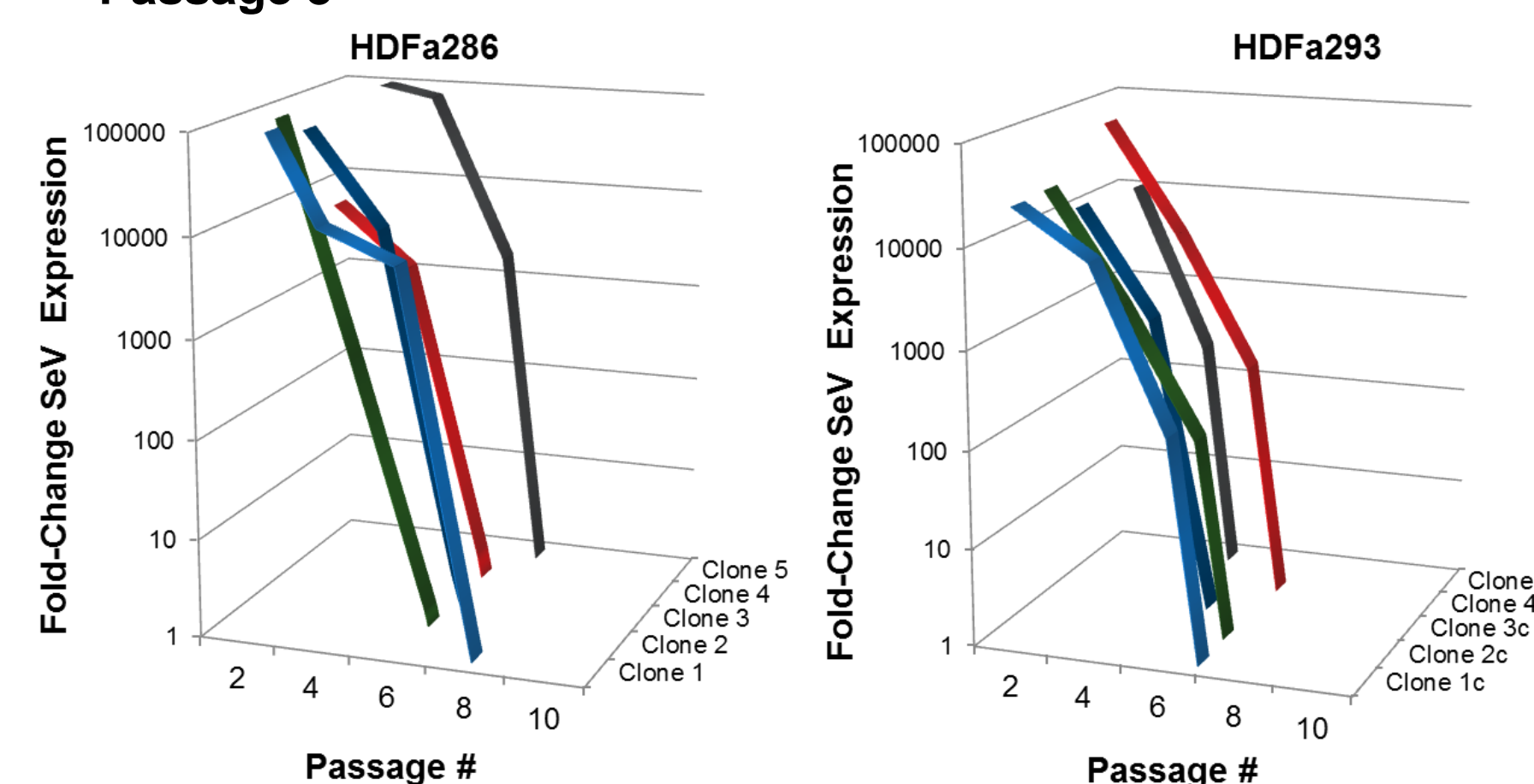


Figure 3. Human Dermal Fibroblasts isolated and expanded in xeno-free medium have typical morphology and a normal karyotype

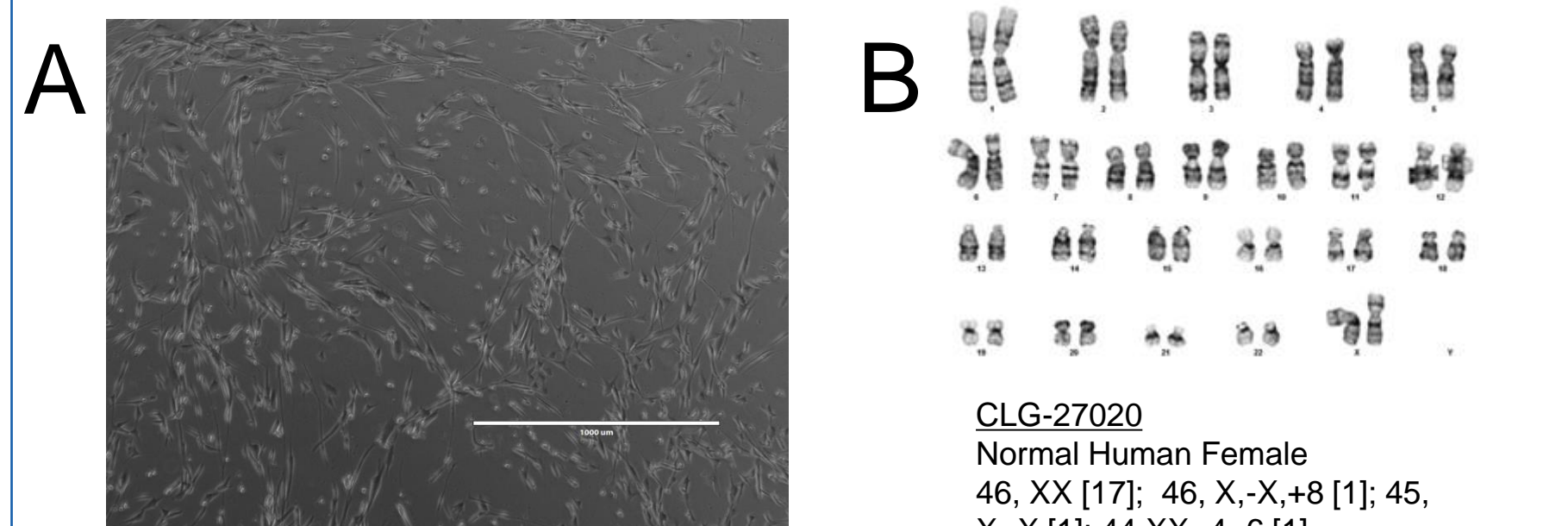
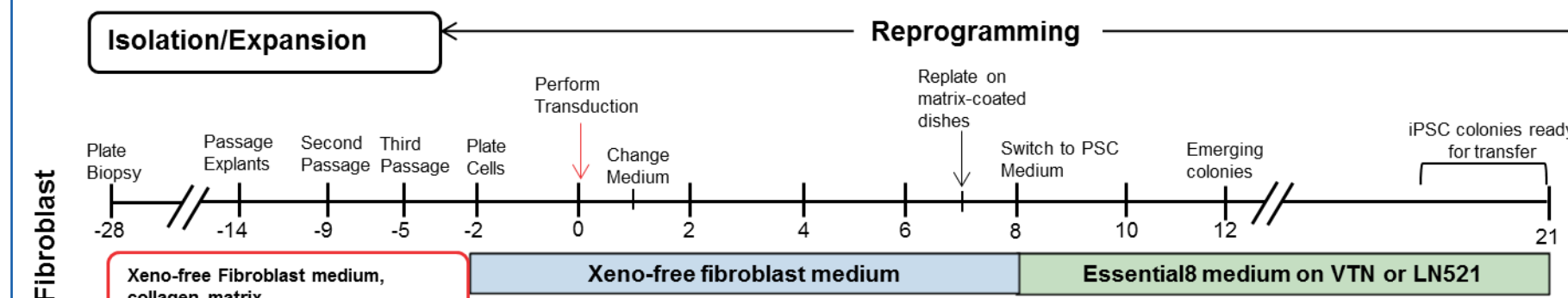


Figure 4. iPSCs derived in a completely xeno-free workflow have a normal karyotype, expression of self-renewal markers, and display functional pluripotency

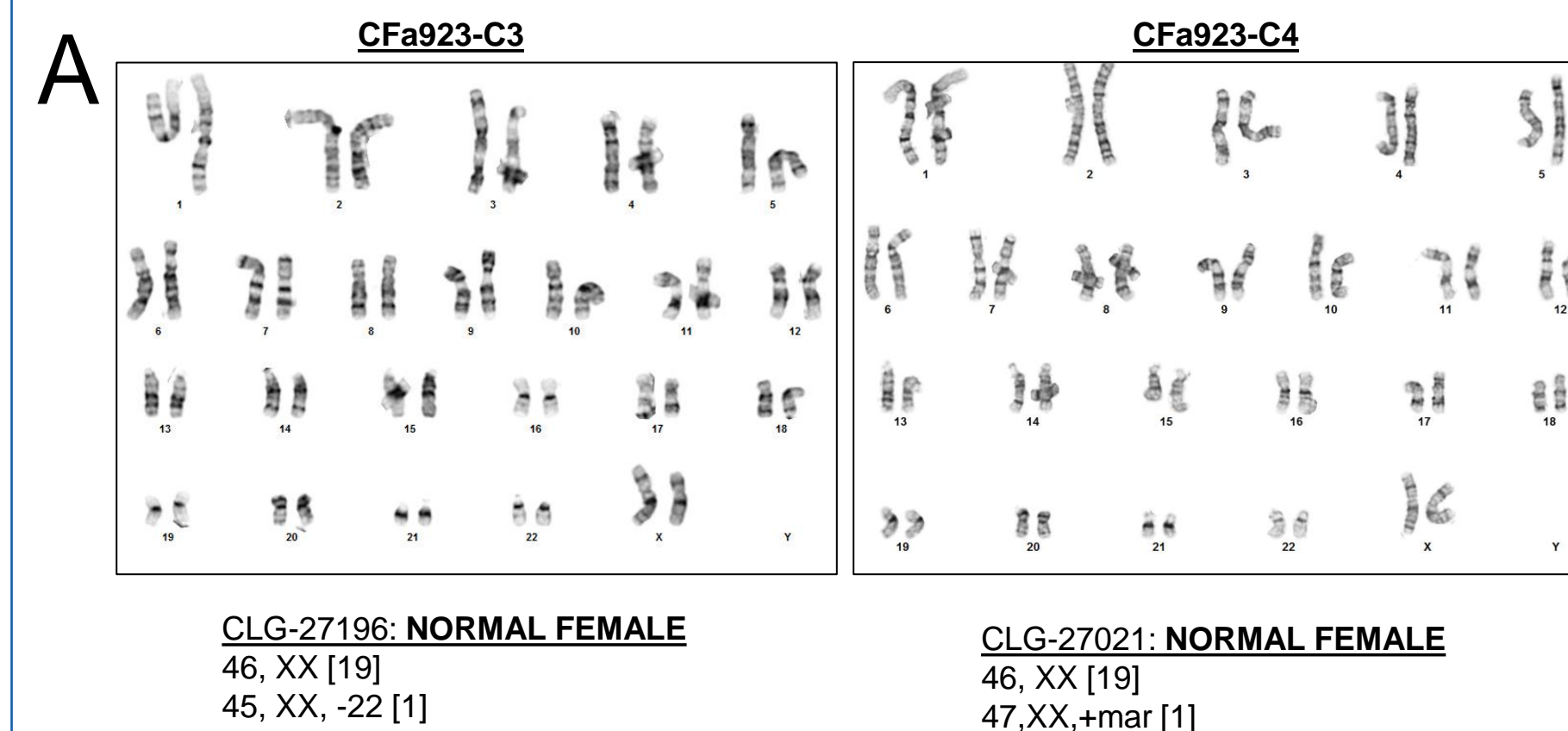
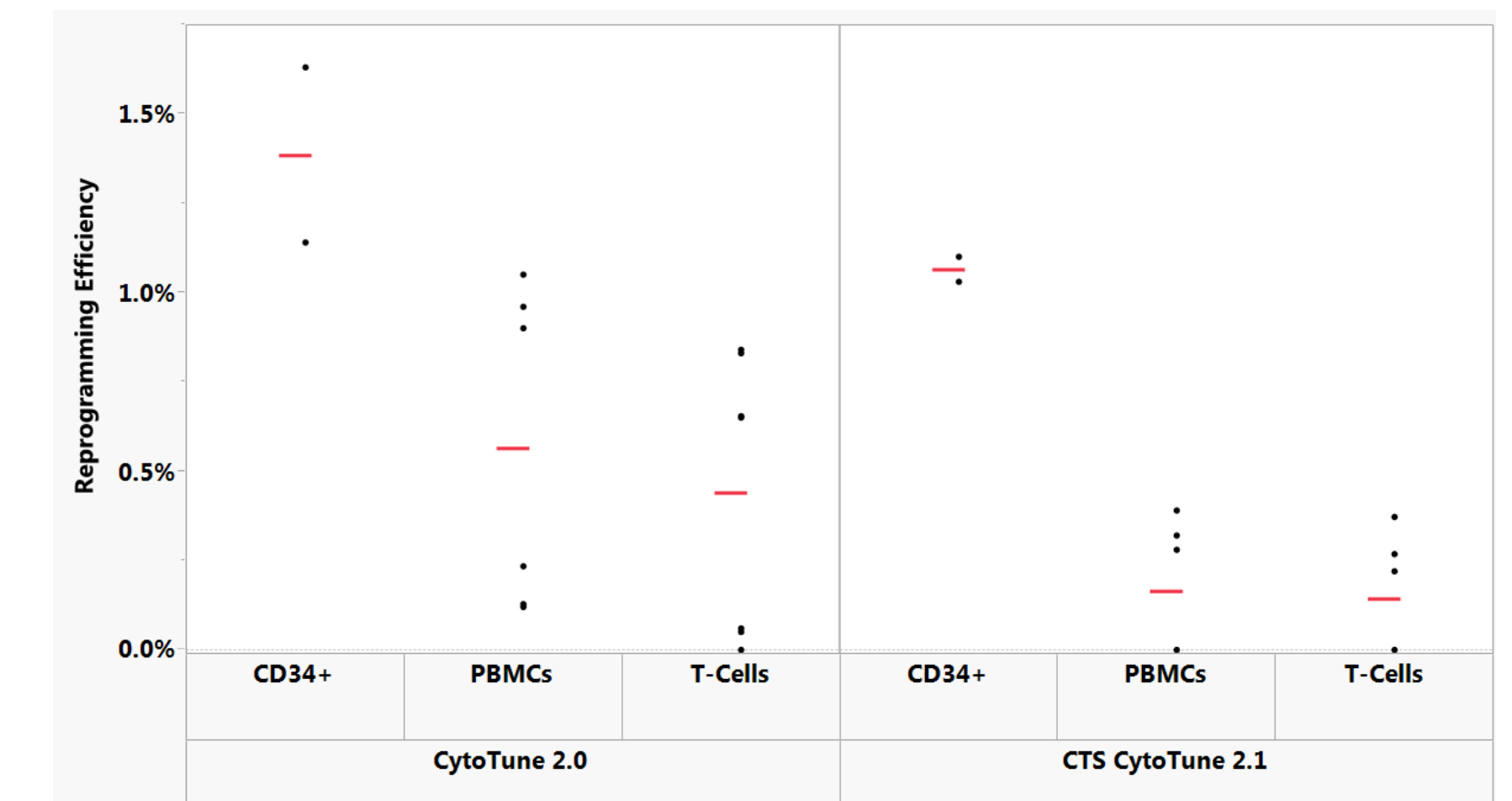
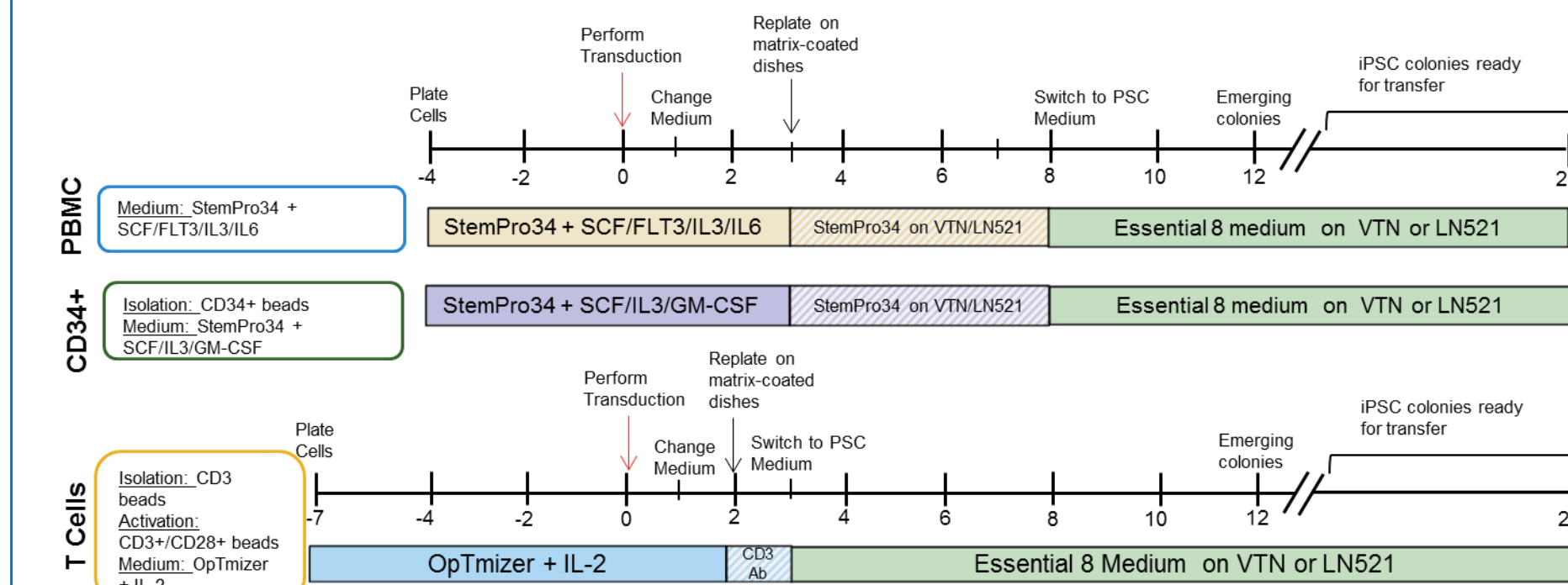


Figure 5. CTS CytoTune 2.1 successfully reprograms blood-derived cells



Three different types of blood-derived cells (CD34+ Cells, PBMCs, and T-Cells) were reprogrammed with CTS CytoTune 2.1, according to the schematic above. At 21 days after transduction, cells were stained for AP, and reprogramming efficiency was calculated as the number of AP positive colonies, relative to the number of cells replated.

CONCLUSIONS

CTS CytoTune 2.1 can successfully reprogram HDF and blood derived cells (CD34+, PBMCs, T-Cells); though efficiency is about 2 fold lower than reprogramming with CytoTune 2.0.

The reduction in reprogramming efficiency is likely due to replacement of C-Myc with L-Myc, and not due to the xeno-free formulation, or GMP manufacturing.

Ten CTS CytoTune 2.1-derived iPSC clones from two different donors showed no detectable virus by Passage 8.

Human dermal fibroblasts derived and expanded in completely xeno-free conditions showed a normal karyotype.

iPSC generated with CTS CytoTune 2.1 in completely xeno-free conditions - from fibroblast derivation through iPSC expansion - displayed all of the standard characteristics of iPSC (i.e. normal karyotype, self-renewal marker expression, functional pluripotency).

REFERENCES

1. Nakagawa M, Takizawa N, Narita M, Ichisaka T, Yamanaka S (2010). Promotion of direct reprogramming by transformation-deficient Myc. Proc Natl Acad Sci U S A. 2010 Aug 10;107(32):14152-7.
2. Seki T, Yuasa S, Fukuda K (2012) Generation of induced pluripotent stem cells from a small amount of human peripheral blood using a combination of activated T cells and Sendai virus. Nat Protoc;7(4):718-28.

TRADEMARKS/LICENSING

© 2016 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. CytoTune is a trademark of ID Pharma Co., Ltd. Essential 8 is a trademark of Cellular Dynamics International, Inc. TaqMan is a registered trademark of Roche Molecular Systems, Inc., used under permission and license. Vector Red® is a registered trademark of Vector Laboratories.

ThermoFisher
SCIENTIFIC