



Characterization of hTERT-immortalized Prostate-derived Stromal and Epithelial Cells: an Authentic *in vitro* Model for Tumor Microenvironment Studies

Luis G. Rodriguez¹, Russell E. McDaniel¹, Xiangshan Zhao¹, Elizabeth Turner², Christopher Annesi¹, and Chaozhong Zou¹

1. ATCC Cell Systems, 217 Perry Parkway, Gaithersburg, MD 20877, USA
2. ATCC, 217 Perry Parkway, Gaithersburg, MD 20877, USA

Abstract

Tumor development begins with mutational changes to the genetic makeup of a cell; tumor progression is not solely determined by the mutated cell, but also by the tumor's microenvironment. Prostate cancer, a leading cancer diagnosed in men, has been determined to be highly influenced by its surrounding stroma, particularly fibroblasts. It has been demonstrated that cancer-associated prostate fibroblasts (CAFs) differ from normal-associated prostate fibroblasts (NAFs). However, human prostate cancer model systems have focused largely on prostate cancer epithelial cells. Currently, a need exists for a more physiologically relevant human cell model system to study prostate cancer progression within the context of its tumor microenvironment. In this study, we characterized three prostate-derived cells: CAFs, NAFs, and prostate epithelial cells (PrEs); all three lines were immortalized by human telomerase reverse transcriptase (hTERT) alone, and have been continuously passaged for more than 40 PDL in our hands. Our data shows that the hTERT-immortalized CAFs proliferate faster than the NAFs; in addition, both CAFs and NAFs express fibroblast markers such as TE7 and alpha smooth muscle actin (α -SMA), while neither cell line expresses epithelial markers such as CK14. Both CAFs and NAFs also express elevated levels of α -SMA upon TGF- β stimulation. All three prostate-derived cells weakly express the prostate specific marker AR, and show similar markers staining after long time passaging. Importantly, conditioned media collected from CAFs promotes tumor cell growth better than NAF conditioned media. In conclusion, CAFs, NAFs, and immortalized PrEs may provide a very valuable model system for the study of prostate cancer cell progression and tumor microenvironment studies.

Introduction

Prostate cancer remains a common cancer with an estimated 180,000 new cases each year, representing about 11% of all new cancer cases in the US in 2016. Over two million men are estimated to be living with prostate cancer and men have a 13 percent chance of being diagnosed with prostate cancer at some point in their lifetime¹. Mortality rates for prostate cancer have remained consistent over the past decade². Further, treatment options have not changed, although prostate cancer animal models are available for use to aid in development of new treatments. In fact, a comprehensive analysis of genetically engineered mouse models identified that invasive prostate cancers are associated with the presence of fibroblast or myofibroblast cells. Additionally, these cells were later observed within a subset of human cancers³. It is known that prostate cancer progression results not only from genetic changes within a cell, but also by the tumor microenvironment⁴, which includes the underlying connective tissue and fibroblasts, myofibroblasts, endothelial cells, and immune cells. However, the mechanisms behind prostate cancer progression remain poorly understood, due to a lack of human cellular models. Developing human cell models is difficult, particularly human prostate cells that are difficult to grow continuously in culture. Thus, an *in vitro* human cell model that incorporates normal prostate cells and their surrounding stromal cells is vital to understanding prostate cancer progression mechanisms.

Results

1. The hTERT Immortalized Prostate Derived Cells Grow Continuously

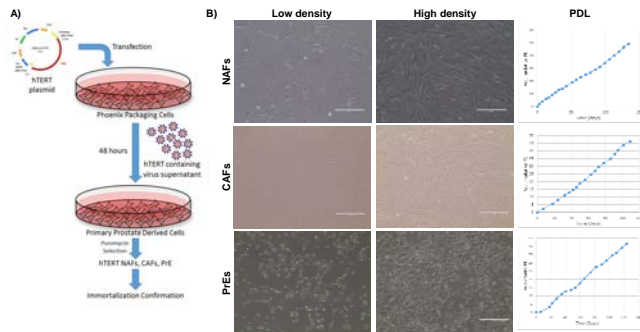


Figure 1. Prostate-derived fibroblast and epithelial cells grow continuously in culture. (A) NAFs, CAFs, and PrEs were immortalized by introducing hTERT alone⁵, selected with puromycin. Successful introduction of hTERT was confirmed by TeloTAGGG[™] telomere length assay and telomeric repeat amplification protocol (TRAP) assay. (B) The selected clones grew continuously in culture for more than forty population doublings (PD). As depicted in the graphs, CAFs grow faster than NAFs, reaching forty population doublings in under 100 days. (n = 1). Scale bar, 400 μ m.

2. Prostate-associated Fibroblasts Express Fibroblast-specific Markers

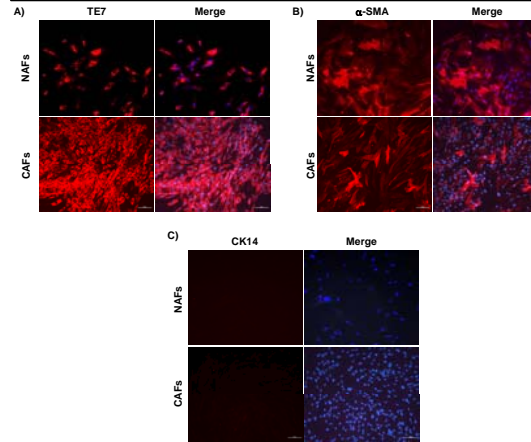


Figure 2. NAFs and CAFs express fibroblast specific markers. NAFs and CAFs cells were fixed using 4% paraformaldehyde, then immunostained with A) an anti-human fibroblast monoclonal antibody (TE7), B) alpha-smooth muscle actin antibody (α -SMA), and C) a cytokeratin 14 antibody (CK14). The nucleus was stained with DAPI, then imaged with a Nikon fluorescent microscope. Both cell lines positively stained for TE7 and α -SMA, but not for CK14, suggesting CAFs and NAFs are mesodermal derived human connective tissue. (n=5). Scale bar, 100 μ m.

3. Normal Prostate Epithelial Cells Express Epithelial-specific Markers

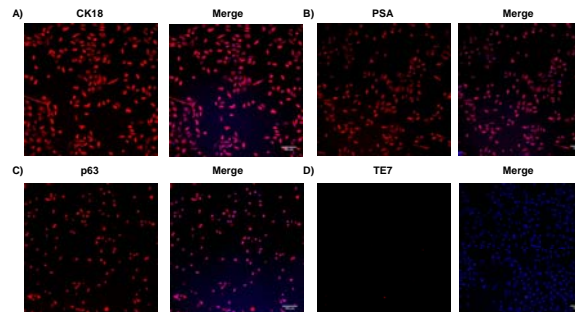


Figure 3. PrEs express epithelial markers, but do not express a fibroblast marker. PrEs were fixed using 4% paraformaldehyde, then immunostained with A) Cytokeratin 18 antibody (CK18), B) Prostate Specific Antigen (PSA) antibody, C) p63 antibody and D) the anti-human fibroblast monoclonal antibody (TE7). Nuclei were stained with DAPI and imaged at 10X using a CellSight CX7 HCS instrument. PrE cells show expression of CK18, PSA and p63, but do not express TE7, suggesting PrE cells are of prostate origin. (n=2). Scale bar, 100 μ m.

References

- National Cancer Institute. Cancer Stat Facts: Prostate Cancer, 2016; Available from: <https://seer.cancer.gov/statfacts/html/prost1.html>.
- American Cancer Society. Survival Rates for prostate cancer, 2017; available from <https://www.cancer.org/cancer/prostate-cancer/detection-diagnosis-staging/survival-rates.html>
- Ittmann M, et al. Animal Models of Human Prostate Cancer: The Consensus Report of the New York Meeting of the Mouse Models of Human Cancers Consortium Prostate Pathology Committee. *Cancer Res* 73(9): 2718-36, 2013.
- Krušlin B, Ullamec M, Tomas D. Prostate cancer stroma: an important factor in cancer growth and progression. *Bosn J Basic Med Sci* 15(2):1-8, 2015.
- Kogan I, et al. hTERT-immortalized prostate epithelial and stromal-derived cells: an authentic *in vitro* model for differentiation and carcinogenesis. *Cancer Res* 66(7):3531-40, 2006.

©2017 American Type Culture Collection. The ATCC trademark and trade name, and any other trademarks listed in this publication are trademarks owned by the American Type Culture Collection unless indicated otherwise.

4. α -Smooth Muscle Actin is Upregulated in Prostate-derived Fibroblasts

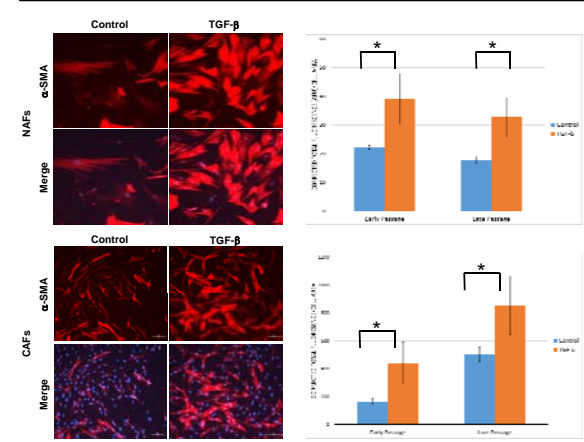


Figure 4. α -SMA upregulation in NAFs and CAFs cells when stimulated with TGF- β . NAFs and CAFs were untreated (control) or treated with 1ng/ml TGF- β . After 24 hours, cells were fixed with 4% paraformaldehyde and immunostained with an α -SMA antibody, counterstained with DAPI, then imaged with a Nikon fluorescence microscope. Multiple fields of view were quantified for total cell fluorescence and normalized by the total area of the cells using ImageJ software. α -SMA levels increased in both early passage and late passage NAFs and CAFs cells after TGF- β stimulation. (n=3, Students t-test, * P<0.05). Scale bar, 100 μ m.

5. CAF-Conditioned Media Promotes Tumor Cell Growth.

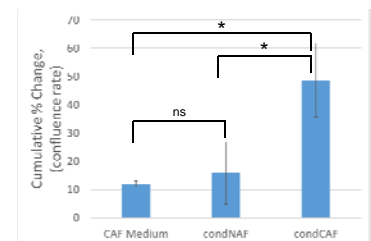


Figure 5. Human prostate carcinoma cells grow faster when cultured in conditioned CAF medium than in conditioned NAF medium. LnCap clone GFC cells (ATCC[®] CRL-1740[™]) were plated onto a 12-well plate at subconfluent levels, allowed to adhere, and then incubated in an Incucyte FLR[®] imaging system in the presence of four distinct growth mediums, including: standard growth medium, conditioned normal-associated fibroblast (condNAF) medium, unconditioned cancer-associated fibroblast (CAF) medium, and conditioned cancer-associated fibroblast medium (condCAF). Growth was then monitored by confluence. The confluence rate was determined by the slope of the line during the exponential phase of growth. Percent change was calculated by normalization with growth in the standard growth medium. The data demonstrated that LnCap clone GFC cells have a faster growth rate in conditioned condCAF, suggesting that cancer-associated fibroblasts, from which the medium was collected, may promote cancer cell growth. (n=3, Students t-test, ns = not significant, * P<0.05).

Summary

- Prostate hTERT-immortalized fibroblast and epithelial cells have been grown continuously for over 100 days without a change in growth rate or senescence.
- Normal-associated and cancer-associated fibroblasts express fibroblast specific markers, but do not express an epithelial specific marker.
- Prostate-derived fibroblasts respond to stimulation with TGF- β , by increasing expression of α -SMA.
- Normal prostate epithelial cells express epithelial specific markers, but do not express a fibroblast specific marker.
- Medium collected from cancer-associated fibroblast cells made cancer cells grow faster than medium collected from normal-associated fibroblast cells.
- These three characterized cell lines provide a human cell model for the study of prostate cancer progression.