

## Abstract

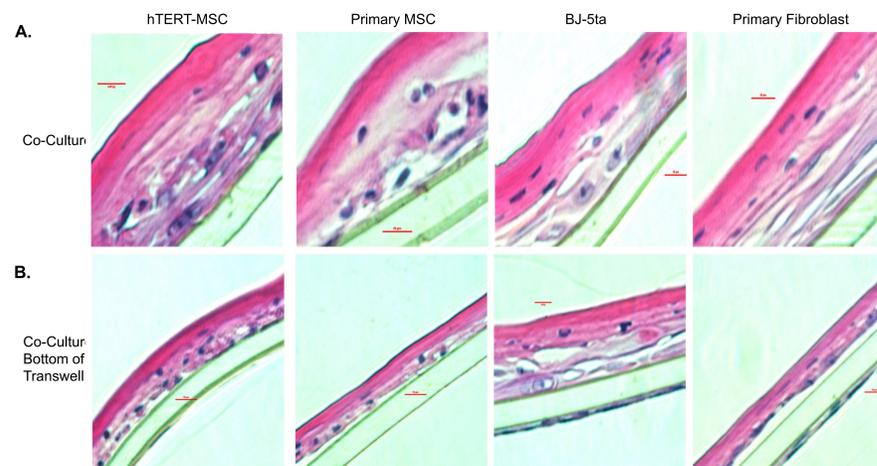
In this study, we compared primary keratinocytes (ATCC® No. PCS-200-010) to hTERT immortalized keratinocytes (Ker-CT; ATCC® No. CRL-4048™), co-cultured with either primary fibroblasts (ATCC® No. PCS-201-010), primary adipose-derived mesenchymal stem cells (MSCs; ATCC® No. PCS-500-011), hTERT-immortalized fibroblasts (BJ-5ta; ATCC® No. CRL-4001™), or hTERT-immortalized MSCs (hTERT-MSCs; ATCC® No. SCRC-4000™). We confirmed that both primary keratinocytes and Ker-CT are able to fully differentiate into skin equivalents in a 3D culture model when co-cultured with primary fibroblasts, primary MSCs, BJ-5ta, or hTERT-MSCs. To confirm the functionality of the co-culture models, both the primary keratinocytes and the Ker-CT air-liquid interface (ALI) co-cultures were subjected to a scratch assay. Re-epithelialization occurred in both cell lines, and interleukin 8 (IL-8) showed an increase in expression from day 0 to day 1 and 3, corresponding to migration of cells into the wound. The continuous nature of the Ker-CT cell line makes it an invaluable model for the research of keratinocyte biology, as it eliminates the issue of short life span and donor variation seen with primary cells.

## Introduction

- Keratinocytes undergo a program of terminal differentiation that results in a potent barrier against microbial infection, water loss, and chemical attack.
- Primary keratinocyte cultures are important cell models for the study of normal and pathological biology of the cutaneous epithelia.
- Keratinocyte differentiation is promoted by factors secreted by cells in the dermis.
- Primary keratinocytes can form skin equivalents that mimic the architectural features and behavior of normal skin in a 3D organotypic culture model in an ALI.
- Primary keratinocytes have finite lifespan in culture, which greatly restricts their use as an *in-vitro* cellular model.
- The constitutive overexpression of the catalytic subunit of telomerase, hTERT, allows for extended passaging of cells with the benefit of little loss of normal cellular physiology.
- hTERT-immortalized keratinocytes cultured with primary fibroblasts, primary MSCs, BJ-5ta, or hTERT-immortalized MSCs comprise a 3D model of the epidermis with high value for investigating drug delivery, toxicity of cosmetics, and the pathology of psoriasis.

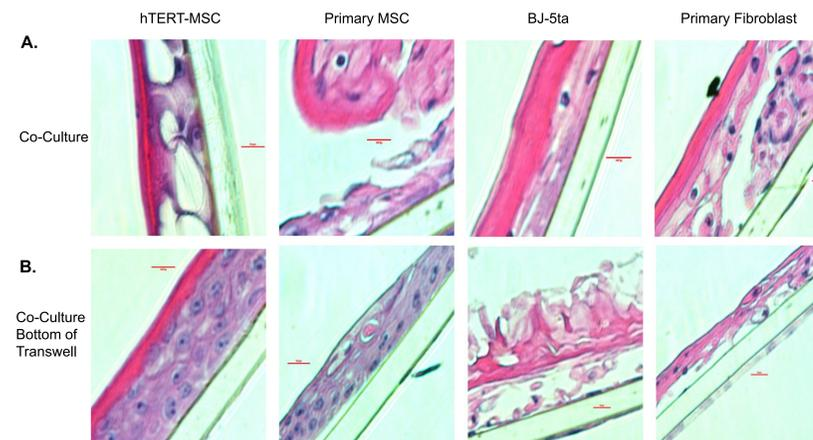
## Results

### hTERT-immortalized and primary stroma-secreting cells directly and indirectly promote primary keratinocytes to form a stratified, 3D, differentiated epidermis

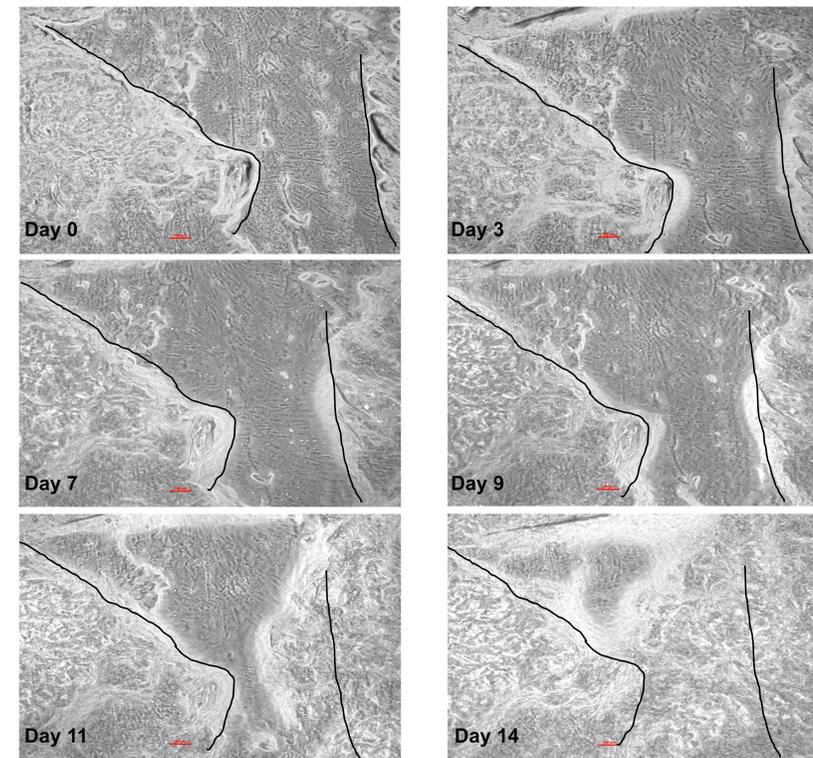


**Figure 1. Hematoxylin and Eosin (HE)-stained cross sections of primary keratinocytes co-cultured with hTERT-immortalized and primary MSCs or fibroblasts differentiated for 21 days.** Primary fibroblasts, MSCs, or their hTERT-immortalized counterparts were seeded onto A) the Transwell™ (Corning) apical chamber or on B) the underside of the Transwell. The fibroblasts and MSCs were allowed to adhere for 24 hours in Mesenchymal Stem Cell Basal Medium (ATCC® PCS-500-030) supplemented with Mesenchymal Stem Cell Growth Kit (ATCC® PCS-500-040). BJ-5ta and primary dermal fibroblasts were grown in Eagle's Minimal Essential Medium (ATCC® 30-2003) supplemented with 10% Fetal Bovine Serum (ATCC® 30-2020). The keratinocytes were then seeded into the apical chamber and allowed to adhere for 4 hours. Cultures were then maintained according to Kalabis *et al.*: Media was changed to epidermalization medium 1 (EPM1). After 3-5 days media was removed and EPM2 was added to the basal chamber only, with media changes every 2-3 days<sup>1</sup>. The cells remained in culture for 21 days, were cross-sectioned, and then subjected to HE staining. All images were taken at 20 x magnification.

### hTERT-immortalized and primary stroma-secreting cells directly and indirectly promote Ker-CT to form a stratified, 3D, differentiated epidermis

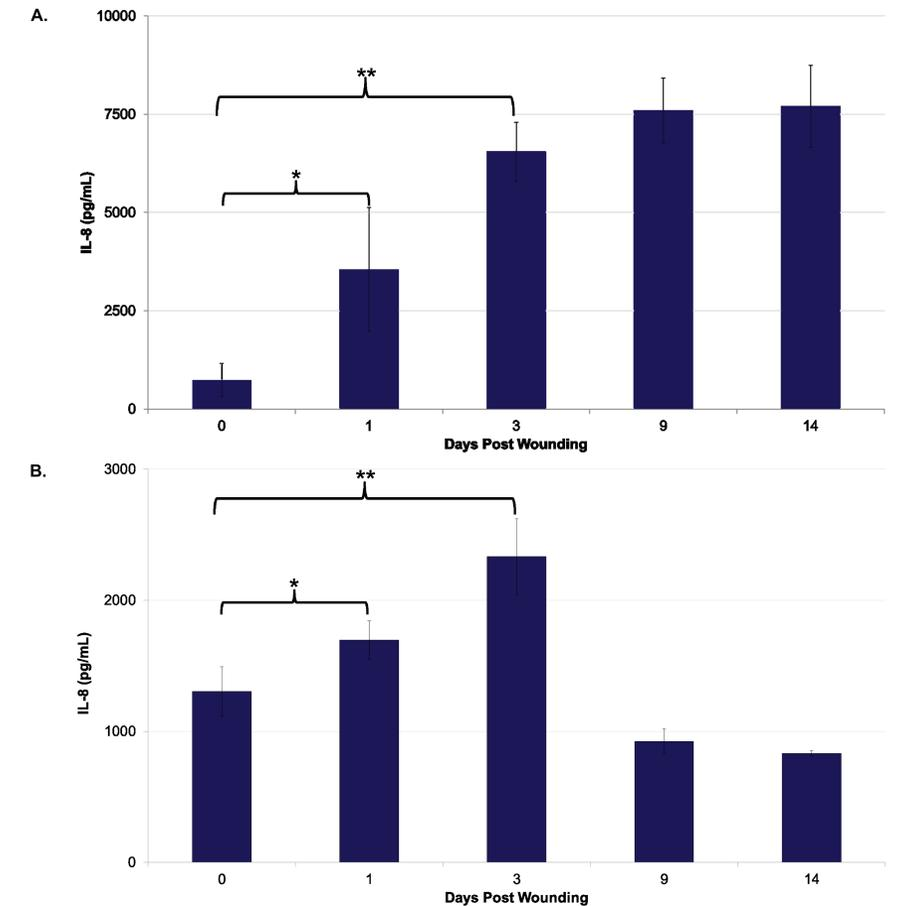


**Figure 2. HE-stained cross sections of Ker-CT co-cultured with hTERT and primary MSCs or fibroblasts, then differentiated for 21 days.** Primary fibroblasts, MSCs, or their h-TERT-immortalized counterparts were seeded onto A) the Transwell apical chamber or B) the underside of the Transwell, then co-cultured with Ker-CT and processed as in Figure 1.



**Figure 3. Ker-CT co-cultured with hTERT-MSC healed after scratch test.** Ker-CT cells were seeded onto Transwells containing hTERT-MSCs in the basal chamber as described in figure 1. After 21 days of differentiation, the apical chamber of the Transwells were scratched using a P1000 tip. Media (EPM2) was changed every 2-3 days for a total of 14 days. Phase images at a 10 x magnification were taken at the indicated days to observe the wound healing. Black lines indicate the size of the wound at day 0.

### Primary keratinocytes and Ker-CT provide an *in vitro* model of epidermal inflammation



**Figure 4. Concentration of IL-8 secreted into the basal media of the primary and immortalized keratinocytes during wound healing.** Basal medium of organotypic A) primary keratinocytes and B) Ker-CT at day 0 (before wounding) and after wounding were collected. Concentrations were then analyzed using the Luminex™100 (Luminex) platform (n=2, performed in triplicate, \*represents  $p < 0.005$  vs day 0, \*\* represents  $p < 0.00001$  vs day 0).

## Summary

- *In vitro* epidermal assay models were established using an ALI Transwell system, primary epidermal keratinocyte cells or Ker-CT, and primary fibroblasts, MSCs, BJ-5ta, or hTERT-MSCs.
- Primary keratinocytes and Ker-CT displayed stratification and terminal differentiation.
- Ker-CT filled in wounds after being subjected to a scratch assay.
- Ker-CT displayed a similar inflammatory response as compared to the primary keratinocytes after wounding.
- Immortalized keratinocyte co-culture model provides a consistent and robust *in vitro* system which increases consistency between assays by decreasing the lot to lot variability of primary cells.

## References

1. Kalabis J, *et al.* Isolation and characterization of mouse and human esophageal epithelial cells in 3D organotypic culture. *Nature Protocols* 7:235-247, 2012.
2. Safferling K, *et al.* Wound healing revisited: A novel reepithelialization mechanism revealed by *in vitro* and *in silico* models. *J Cell Biol* 203(4):691-709, 2013.
3. Sasaki M, *et al.* Mesenchymal stem cells are recruited into wounded skin and contribute to wound repair by transdifferentiation into multiple skin cell type. *J Immunol* 180; 2581-2587, 2008.
4. Lai Y, *et al.* Interleukin-8 Induces the endothelial cell migration through the activation of phosphoinositide 3-kinase-rac1/rhoA pathway. *Int J Biol Sci* 7(6):782-791, 2011.

© 2014 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. Transwell and Corning are registered trademarks of Corning Incorporated. Luminex is a trademark of Luminex Corporation.