

## Objective

To evaluate the anti-aging efficacy of topically applied cosmetic ingredients and formulations by measuring the expression of ECM components in the EpiDermFT™ *in vitro* human skin model.

## Methods

EpiDermFT tissues were treated with 25 µL of each formulation topically and then cultured at 5% CO<sub>2</sub>, 37 degrees, 95% humidity for 24hrs. After treatment, EpiDermFT tissues were processed for total RNA isolation. Total RNA was utilized for gene expression analysis by quantitative PCR.

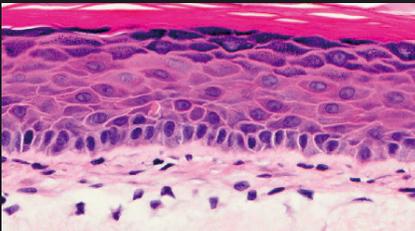


Figure 1. Histology of EpiDermFT™ H&E stained cross-section showing that the tissue morphology of EpiDermFT™ closely parallels that of normal human skin. The epidermis contains basal, spinous, granular and stratum corneum layers, and the dermis contains viable fibroblasts (400X).

## Results

EpiDermFT™ tissues treated with Formulation A showed significant increases in Collagen 1A1, Collagen 3A1 and Elastin gene expression. Tissues treated with Formulation B showed significant increases in COL3A 1 expression (Figure 2).

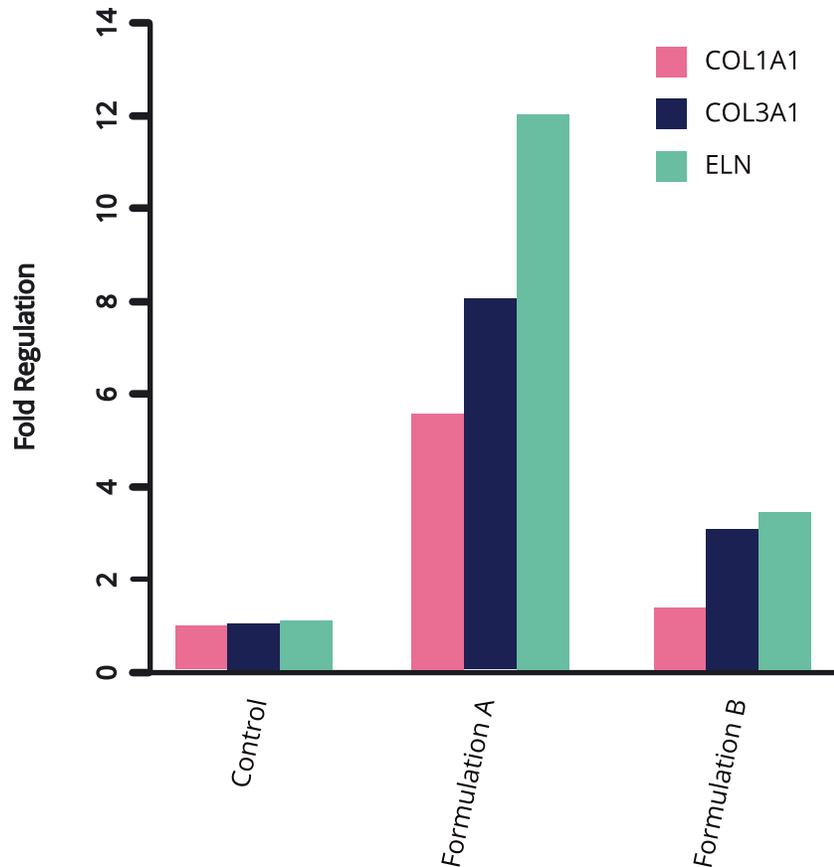


Figure 2. Gene Expression of EpiDermFT™. Genes of interest are compared to untreated controls. Data are presented as the average fold regulation of experimental replicates.

## Conclusion

Evaluation of ECM components by quantitative PCR in the EpiDermFT™ *in vitro* human skin model can be used in efficacy and claims substantiation studies.