

## Objective

Accumulation of reactive oxygen species (ROS) in dermal fibroblasts can cause oxidative stress and premature skin aging. To identify compounds which can prevent free radical buildup and/or treat oxidative stress inducers, Mattek's Normal Human Dermal Fibroblasts (NHDFs) were used to screen potential antioxidants.

## Methods

Mattek's NHDFs (NHDF-CRY-NEO) were cultured in DMEM-10 according to manufacturer's protocol (Figure 1). Cells were pre-treated with increasing concentrations of potential antioxidant for 24 hours. Cells were incubated with 100  $\mu$ M of 2'-7'-dichlorodihydrofluorescein diacetate (DCFH-DA, 10 $\mu$ M) for 1 hour. Intracellular ROS was induced by a 1 hour incubation with 10mM H<sub>2</sub>O<sub>2</sub>. Levels of 2'-7'-dichlorodihydrofluorescein (DCF), generated by ROS, were measured and changes in 485/528 nm ROS signal were compared to control (Figure 2).



Figure 1. Human small intestine model cultured on PermaCell Insert, CCI24-PET-0.4, showing a highly differentiated structure with villi and brush border formation: A) H&E cross section, B) TEM micrograph.

## Results

NHDFs treated with Antioxidant showed dose-dependent decreases in intracellular reactive oxygen species following induction by H<sub>2</sub>O<sub>2</sub> (Figure 2).

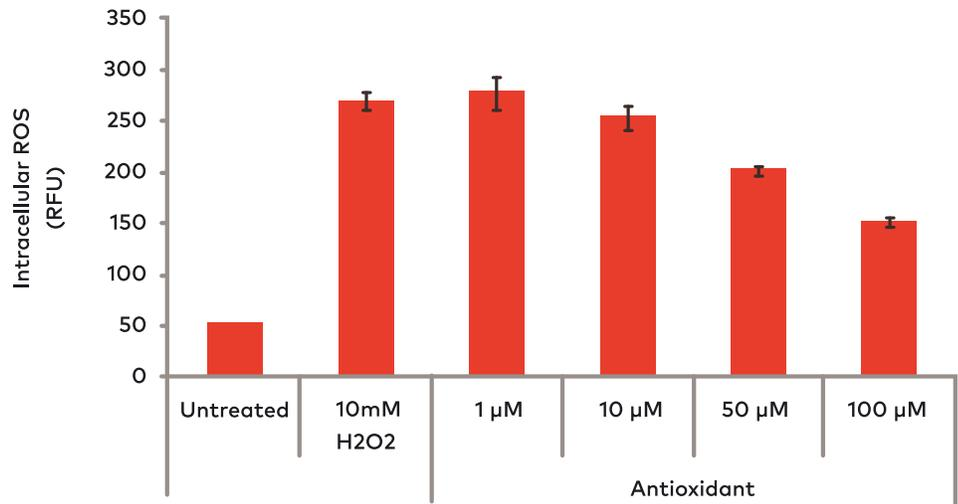


Figure 2. Mattek's Normal Human Dermal Fibroblasts (NHDF-CRY-NEO) were treated with increasing concentrations of antioxidant test compound for 24 hours and then exposed to 10mM H<sub>2</sub>O<sub>2</sub>. Significant decreases in intracellular ROS levels were observed with increased concentrations of Antioxidant.

## Conclusion

Mattek's Normal Human Dermal Fibroblasts can be used to identify antioxidant compounds for personal care and cosmetic product development.