

# Comparison of EMT biomarker expression in 2D monolayer and 3D spheroid cultures in a prostate cancer cell model

#5793

Jen Carlstrom, Jeanine Hinterneder, Lindsay Nelson, and Stephen Hurt

## 1 Introduction

Epithelial-mesenchymal transition (EMT) is characterized by rearrangement of the extracellular matrix (ECM) and differential regulation of ECM proteins. We induced EMT in DU 145 cells using TGF- $\beta$  and phorbol-12-myristate-13-acetate (PMA) and compared expression levels of specific biomarkers, including E-cadherin, fibronectin, and IL-6, using AlphaLISA® and LANCE® (TR-FRET) assay technologies. We confirmed that treatment with TGF- $\beta$  is sufficient for inducing changes in both EMT biomarker expression and in promoting development of characteristic mesenchymal stromal cell morphology in monolayer cultures. However, in 3D spheroid cultures, we only observed a partial EMT response to the same TGF $\beta$  treatment as demonstrated by changes in the expected biomarker expression pattern. Cellular proliferation, growth and vitality were assessed using ATPlite luminescence assays and confocal microscopy of live-stained cells with a high content imaging system. Though we observe increased proliferation in monolayer cultures compared to 3D spheroids, the changes observed in protein expression patterns cannot be sufficiently explained by differences in cell number or viability. These data illustrate the differences in protein expression levels and in cellular tolerance for compound treatment between a human prostate cancer cell line grown in monolayers and those same cells grown in 3D spheroids.

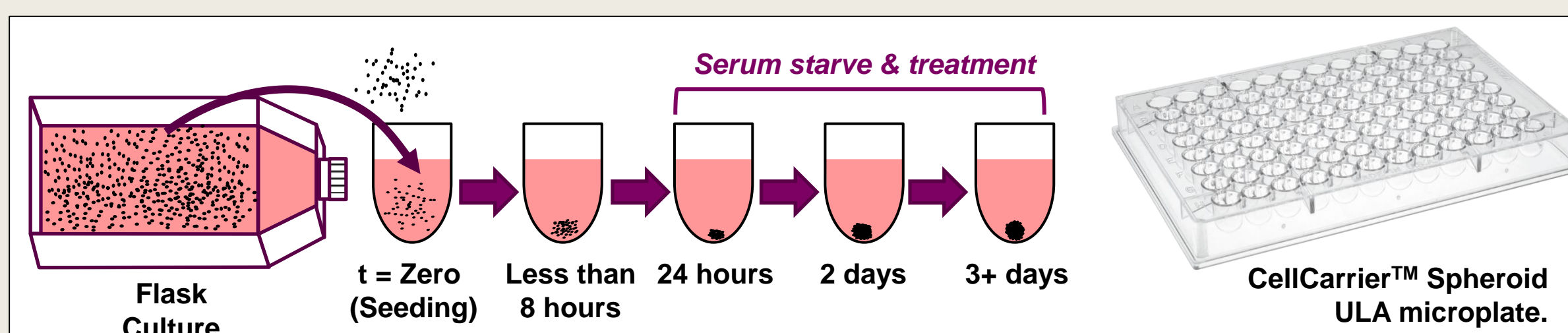
## 2 Materials and Methods

**Cell culture and treatment:** DU 145 cells (ATCC® HTB-81™) were seeded (100  $\mu$ L/well) into PerkinElmer 96-well CellCarrier™ (6005550) or CellCarrier Spheroid ULA 96-well microplate (3D) (6055330) and grown for at least 18 hours (see spheroid growth illustration below). Cells were serum starved for 24 hours prior to treatment with recombinant human TGF- $\beta$ 1 (BioLegend, 580702) or PMA (Sigma cat# P1585) to induce EMT. In some experiments, cells were treated with 2.5  $\mu$ M of TGF- $\beta$  inhibitor SD 208 (Sigma, S7071) for two hours followed by 48 hours with TGF- $\beta$  (5 ng/mL) in a total volume of 100  $\mu$ L.

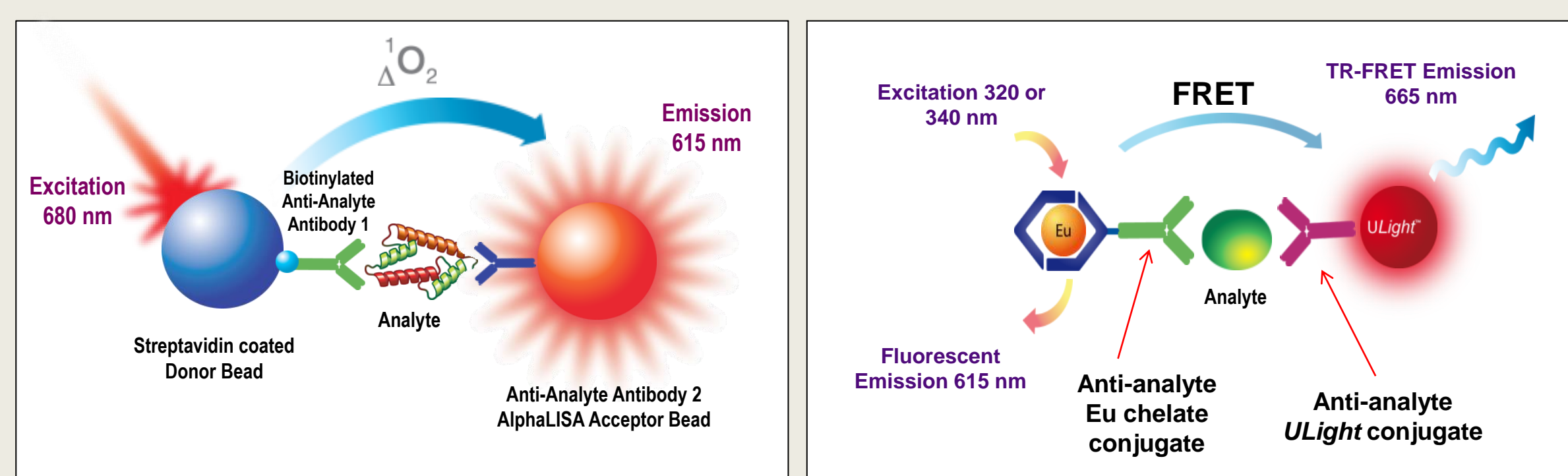
**Biomarker quantification assays:** For biomarker detection assays, 50  $\mu$ L of media was removed (for testing supernatants). Cells were then lysed with 50  $\mu$ L of 2X Alpha SureFire® Ultra lysis buffer (ALSU-LB-10mL) for 10 minutes. Lysates and supernatants were frozen at -80°C and later thawed for testing in AlphaLISA and LANCE assays (see models for each assay type below). For **AlphaLISA assays** detecting E-cadherin (AL370C), fibronectin (AL351C), and human IL-6 (AL223C), 5  $\mu$ L of each lysate or supernatant sample was added to a 384-well white OptiPlate (6007290) and assays were performed according to the manual. For Fibronectin **LANCE Ultra TR-FRET assay** (TRF1351C), 15  $\mu$ L of lysate was added to a 384-well white OptiPlate and assays were performed according to the manual. AlphaLISA and LANCE assays were measured on a standard EnVision® multilabel plate reader using standard Alpha Settings and reading TR-FRET with the Laser excitation.

**ATPlite™ 1step and ATPlite™ 1step 3D assays:** Cellular proliferation, growth & vitality were measured by assessing the concentration of ATP using ATPlite 1step and ATPlite 1step 3D luminescence-based assays following kit protocols. Luminescence was measured using the EnSight™ multimode plate reader.

**Cellular imaging:** For cellular imaging, cultures were first labeled with Hoechst 33342 (Life Technologies, #H3570), Tetramethylrhodamine (TMRM; Life Technologies, #T-668) and CellTox Green (Promega, #G8742). Monolayer cultures were imaged using the cellular imaging module of the EnSight plate reader using Brightfield and UV fluorescence filters and cell numbers were quantified using the Count Nuclei function. 3D Spheroid cultures were imaged with the 10X long WD objectives on the Operetta® and Opera Phenix™ High Content Imaging systems using Brightfield and appropriate fluorescence optics. Cross-sectional spheroid area was measured with Harmony software using an intensity cutoff in the UV channel (Hoechst).



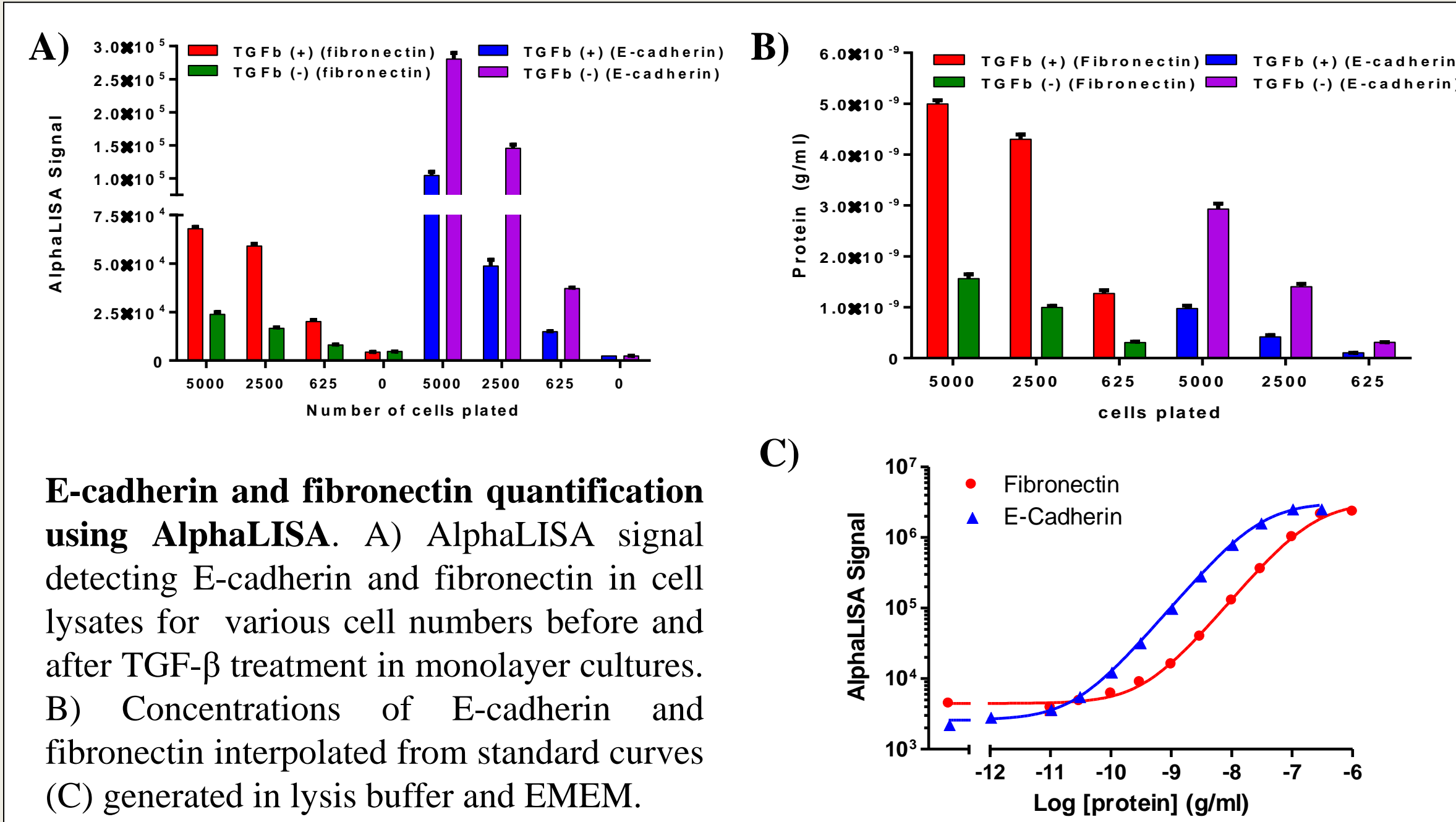
Model of Spheroid culture growth.



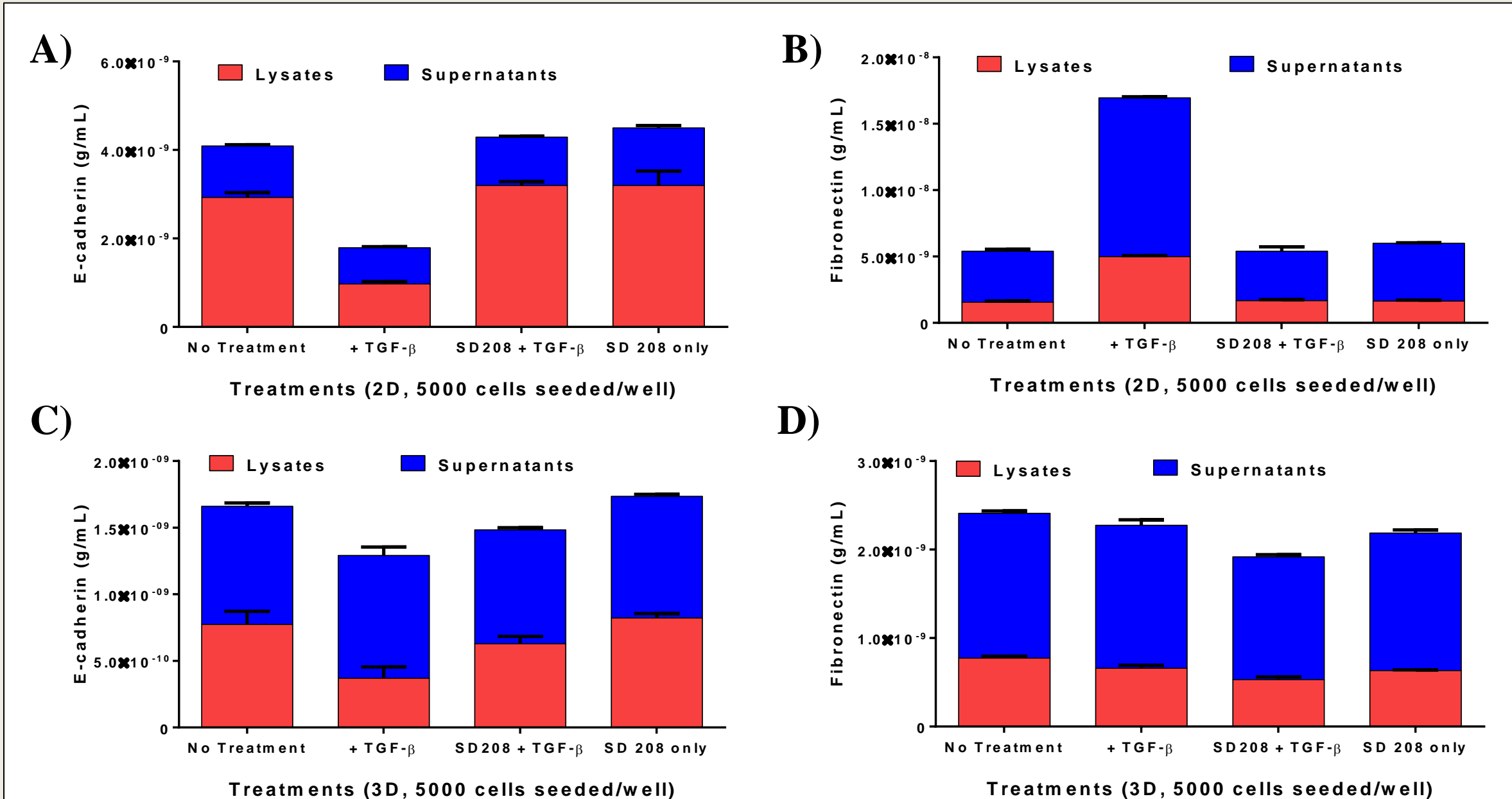
AlphaLISA assay

LANCE TR-FRET assay

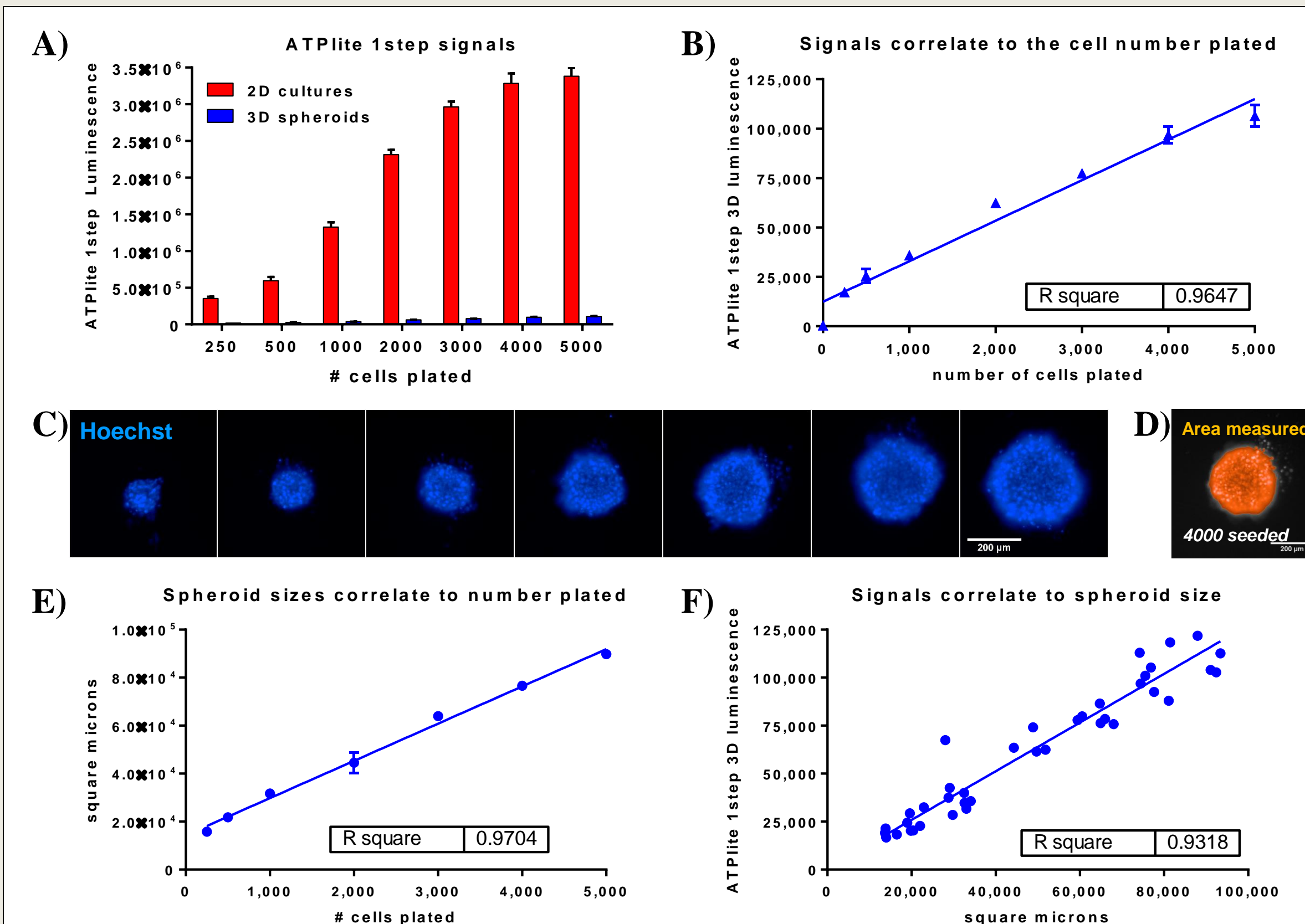
## 3 E-cadherin and Fibronectin modulated by TGF- $\beta$ measured with AlphaLISA in monolayer cultures



## 4 Modulation of E-cadherin & Fibronectin Levels in 2D and 3D cultures

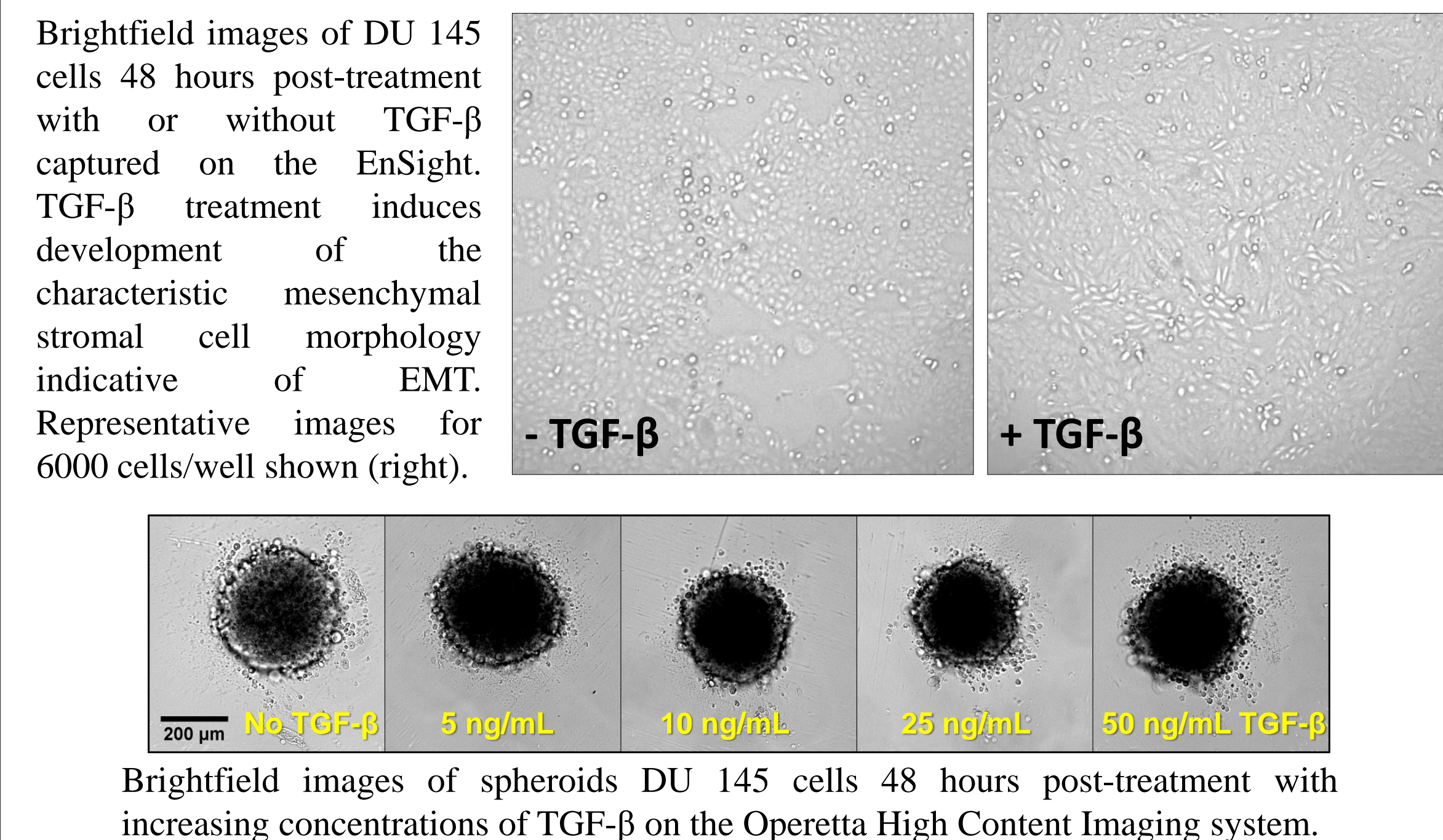


## 5 Assessing cell proliferation in 2D & 3D

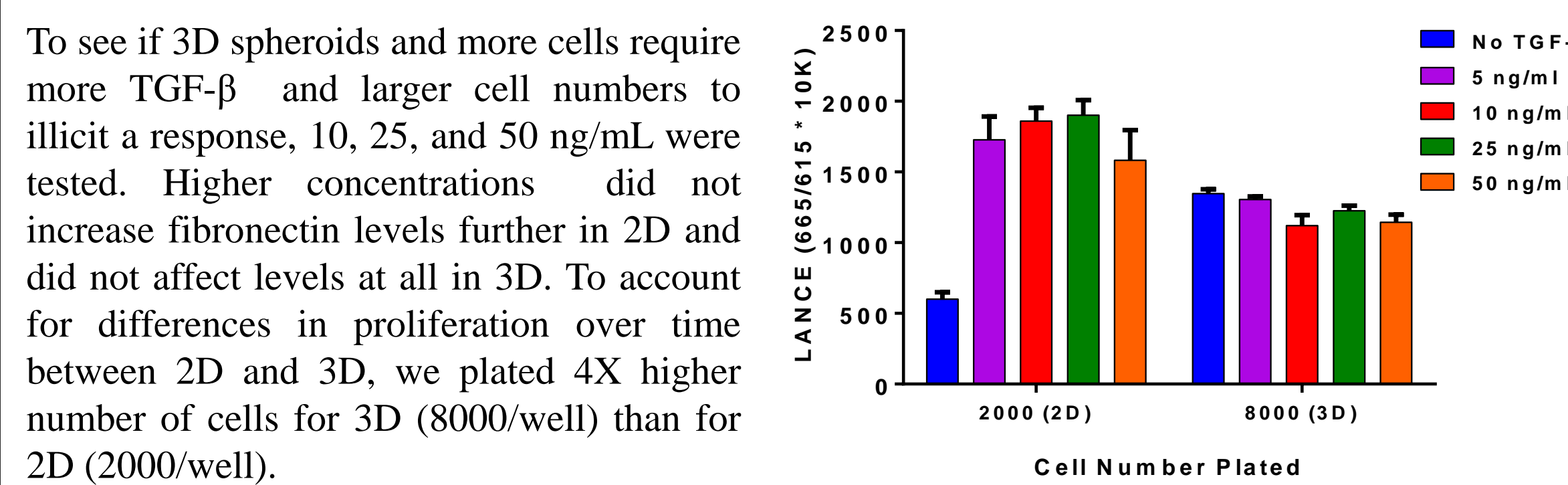


Monolayer (2D) cultures proliferate much more than spheroid (3D) cultures. 2D and 3D cultures were plated at 7 concentrations, grown for 4 days and assessed with ATPlite 1step and ATPlite 1step 3D assays (with the final step being to transfer 50  $\mu$ L of the reaction to a gray OptiPlate for reading luminescence on the EnSight (n=3). 3D cultures were labeled with Hoechst, spheres were imaged on the Operetta (C), and cross-sectional spheroid area measured using an intensity cutoff (D). Spheroid area correlates to number plated (E) and luminescence signals (F).

## 6 TGF- $\beta$ effects on morphology

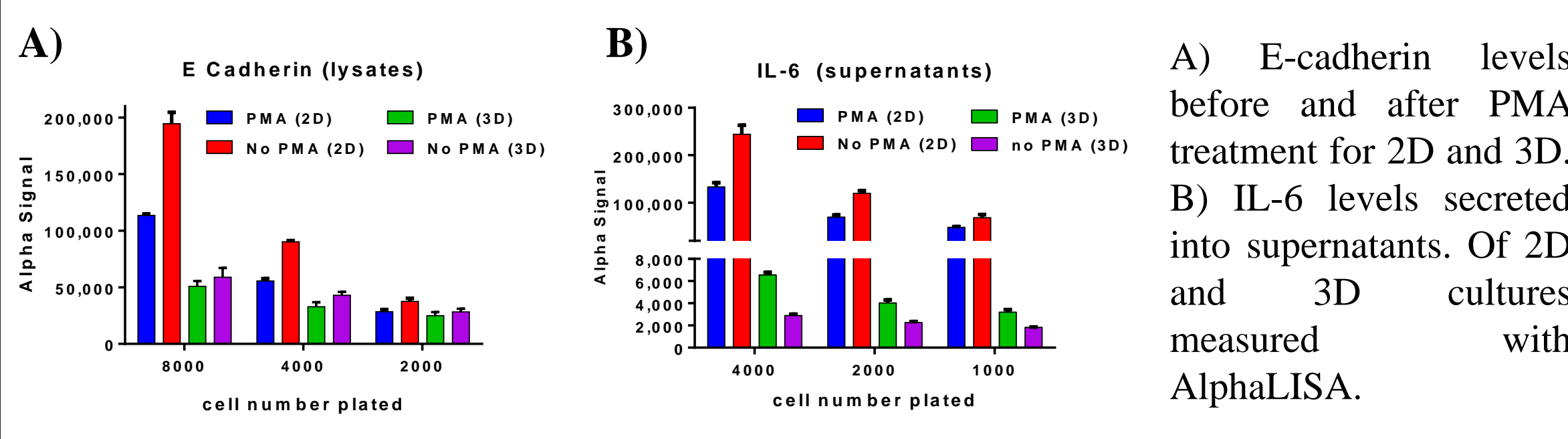


## 7 Examining TGF- $\beta$ effects on Fibronectin expression with LANCE TR-FRET

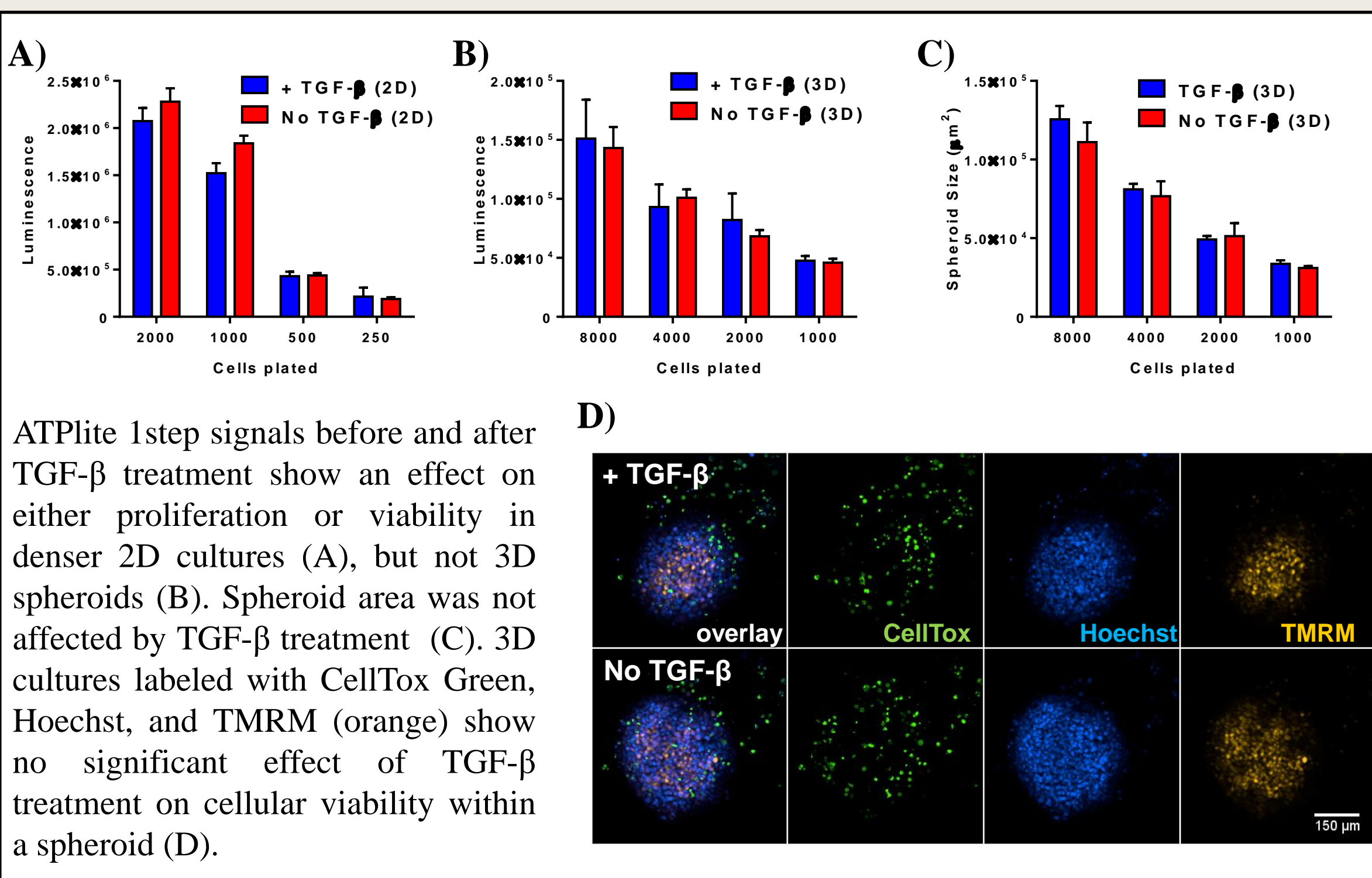


## 8 PMA -induced changes in E-cadherin and IL-6 expression

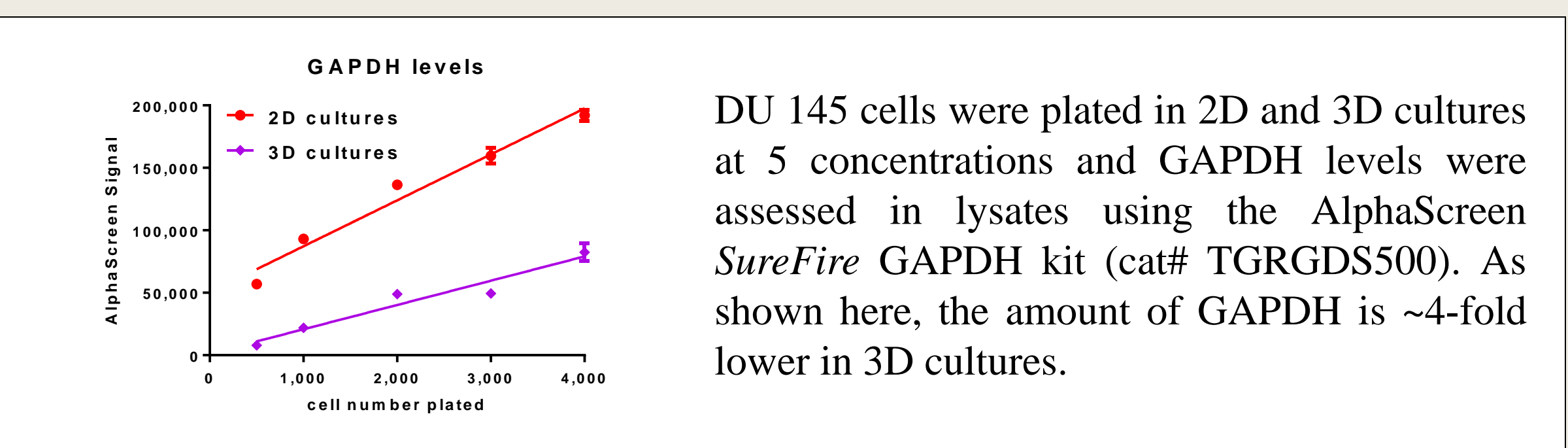
PMA has been shown to induce EMT also. Here, we show E-cadherin levels are decreased in 2D and less so in 3D after treatment with 150 nM PMA. Interestingly, the IL-6 secretion levels increase for 2D and decrease for 3D after PMA treatment.



## 9 Affects of TGF- $\beta$ on cell viability

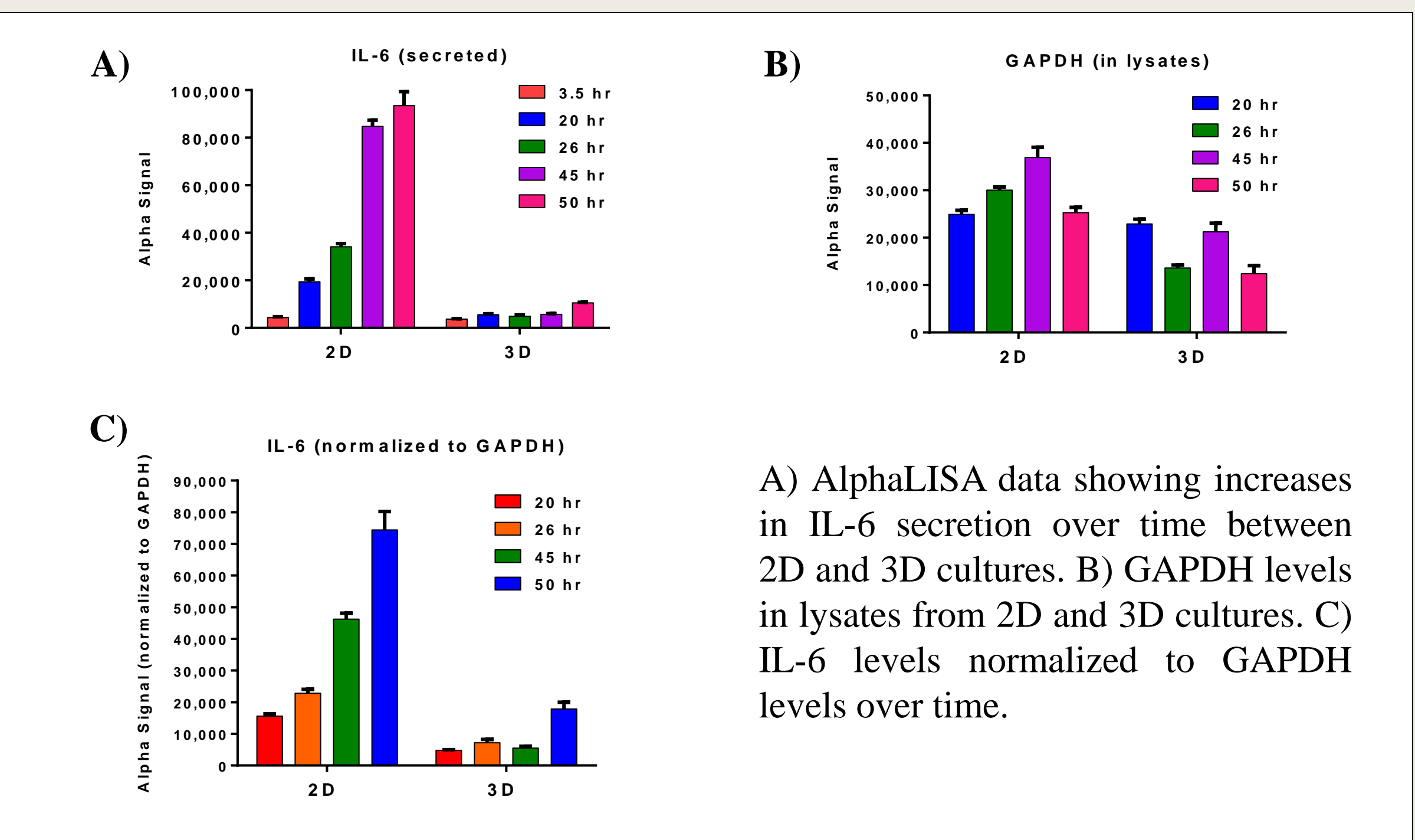


## 10 GAPDH levels may be used to correlate cell number differences between 2D and 3D over time



## 11 Changes in IL-6 levels over time between 2D and 3D

To assess the changes in IL-6 secretion over time, 4000 cells were plated in multiple wells in 2D and 3D. At each time point, supernatants were removed and cells lysed and aliquots were frozen at -20°C. AlphaLISA was used to measure IL-6 levels in the samples and GAPDH was used to normalize to potential differences in cell number due to differences in growth rates. The dramatic difference in IL-6 secretion over time in 2D compared to 3D cannot simply be explained by differences in cell numbers.



## 12 Summary

- We confirmed that treatment of DU 145 cells with TGF- $\beta$  is sufficient for inducing changes in both EMT biomarker expression and cellular morphology in monolayer cultures.
- AlphaLISA and LANCE (TR-FRET) biomarker assays can be used to measure ECM-associated protein modulation caused by human transforming growth factor-beta (TGF- $\beta$ ) induction of EMT in a 3D Spheroid model of human prostate carcinoma.
- E-Cadherin is downregulated by TGF- $\beta$  in both 2D and 3D cultures, whereas Fibronectin is increased significantly only in 2D monolayer cultures.
- Monolayer cultures proliferate considerably more than cells in spheroid cultures.
- PMA and TGF- $\beta$  treatment induced significant differences in IL-6 secretion levels between 2D and 3D cultures.
- TGF- $\beta$  affects proliferation or viability of cells in 2D cultures differently than in 3D cultures.
- These data illustrate the differences in protein expression levels and in cellular tolerance for treatment between a human prostate cell line grown in monolayers versus 3D spheroids.