

## Introduction

Angiogenesis and cell proliferation are fundamental biological processes, and their pathologic states represent two hallmarks of cancer. Central to controlling both processes is the TGF- $\beta$ /BMP signaling pathway, for which the SMAD proteins serve as major downstream effector molecules. Upon ligand binding, TGF- $\beta$  or BMP type I/II receptor heterotetramers phosphorylate receptor-regulated SMADs (R-SMADs). TGF- $\beta$ -receptor binding leads to SMAD2/3 phosphorylation, whereas BMP-receptor binding induces phosphorylation of SMADs 1/5/8. Phosphorylated SMADs (pSMADs) then form a complex with the co-SMAD SMAD4, at which point they undergo nuclear translocation, where they operate as transcriptional modulators of target genes. The magnitude of SMAD phosphorylation and functional output (e.g., stimulated angiogenesis or attenuated cell proliferation) depends on both the ligand concentration and the duration of exposure. In this study, we quantitatively evaluate the phosphorylation states of R-SMADs, as well as their downstream biological effects at three different biological scales: 1) biochemical; 2) cellular; and 3) tissue/3D models. At the biochemical level, the temporal and dose-dependent phosphorylation of R-SMADs in response to growth factor stimulation can be quantified using a quantitative ELISA approach. Growth factor stimulation can be quantified at the cellular level by high-content imaging of SMAD phosphorylation and nuclear translocation, while simultaneously monitoring cell proliferation rates. Lastly, high-resolution confocal imaging can be employed to observe developmental consequences of phospho-SMAD signaling using a HUVEC-based 3D angiogenesis model system.

## Experimental

### Cell lines

Immortalized human fibroblasts (CI-huFIB, part number INS-CI-1010) were a kind gift from InSCREENex GmbH (Braunschweig, Germany). Human Umbilical Vein Endothelial Cells (HUVECs) were purchased from Lonza Bioscience (part number C2519A; Basel, Switzerland). A549 lung carcinoma cells were purchased from ATCC (part number CCL-185; Manassas, VA, USA). CI-huFIBs and A549s cells were cultured in Advanced DMEM (part number 12491; Gibco Thermo Fisher Scientific; Waltham, MA, USA) containing 10% FBS and 1x penicillin/streptomycin/glutamine. HUVECs were maintained in Endothelial Cell Grown Medium-2 BulletKit (EGM-2; part numbers CC-3162; Lonza).

### Growth factor treatment

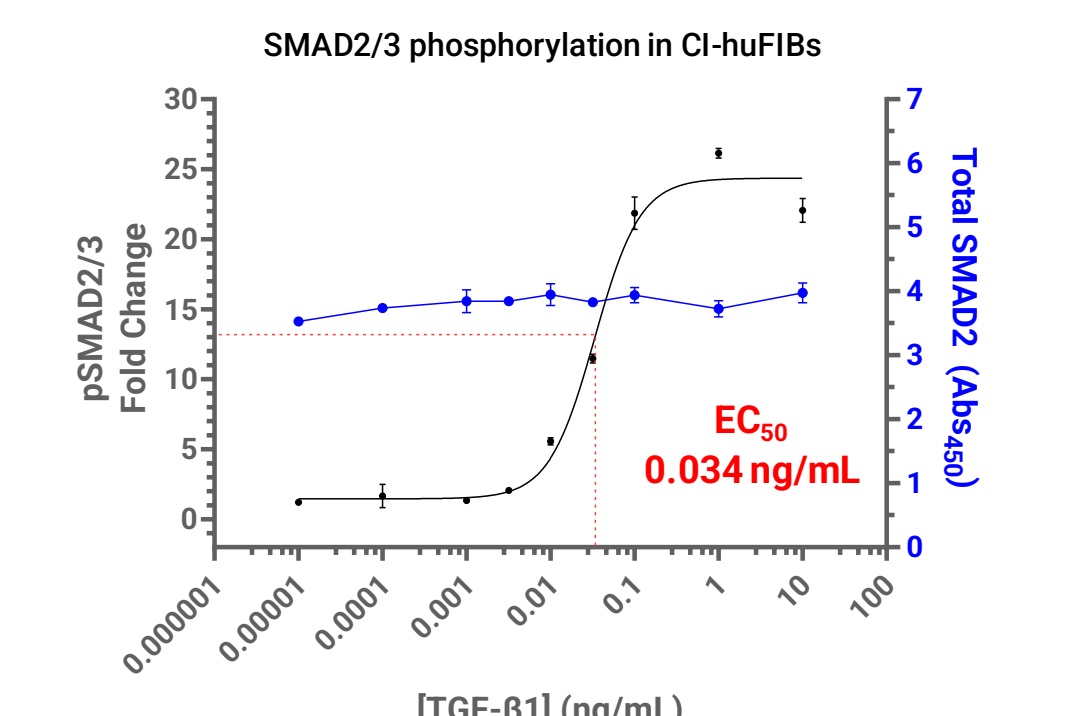
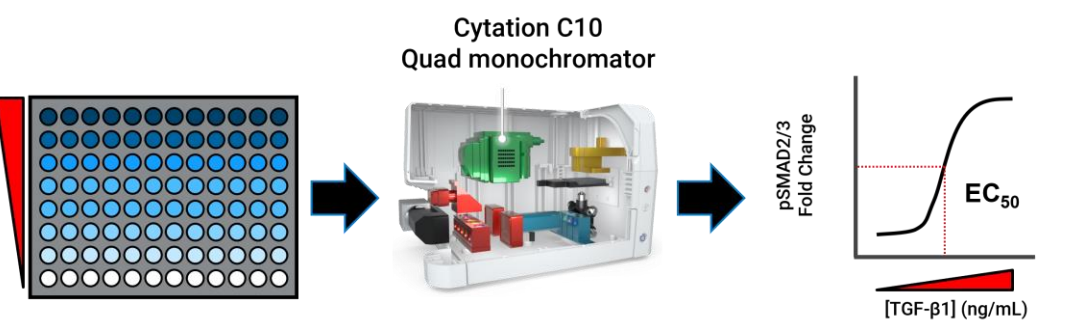
Before exposure with defined concentrations of growth factors, cells seeded in culture plates (12-well for PathScan sandwich ELISA and 96-well for 2D cultures) were subjected to an overnight, 18-hour serum-starvation period using the same basal culture medium as listed above, lacking FBS (A-DMEM) or BulletKit supplements (EBM). Growth factor treatment was conducted by adding hTGF- $\beta$ 1, hBMP-2 (part numbers 75362 and 4697, respectively; Cell Signaling Technology, Danvers, MA, USA), or rhBMP-6 (part number 507-BP; R&D Systems, Minneapolis, MN, USA) diluted in serum-free media to adherent cells at the indicated concentrations and allowed to incubate for 1 hour, followed by 4% paraformaldehyde fixation and immunostaining.

## Results and Discussion

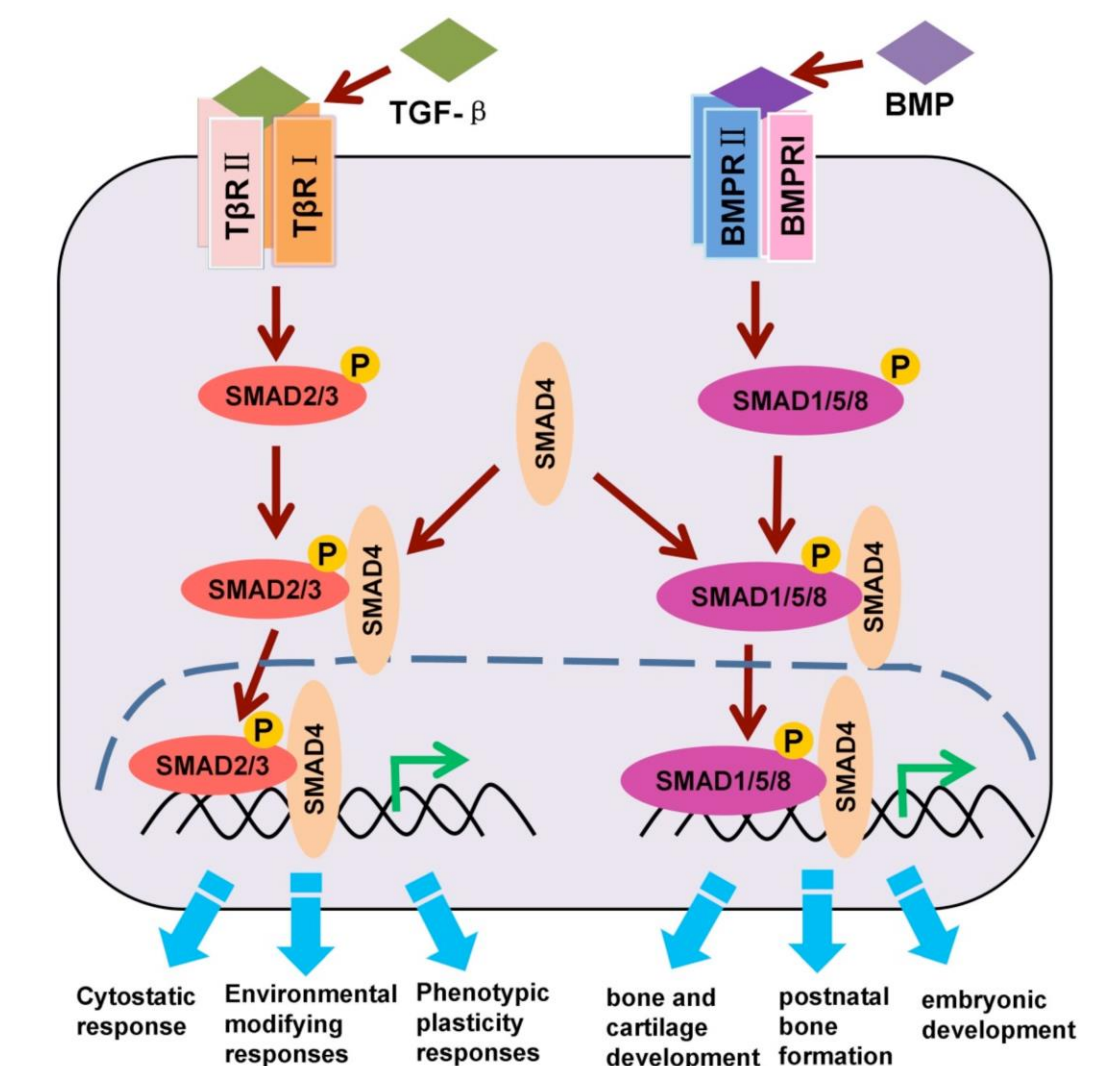
### Biochemical analysis: high-throughput ELISA assay

Cell Signaling Technology PathScan sandwich ELISA kits:

- Phospho-SMAD2 (S465/467)/SMAD3 (S423/425), part number #12001
- Total SMAD2/3, part number #12000



### Canonical TGF- $\beta$ /SMAD and BMP/SMAD signaling pathway



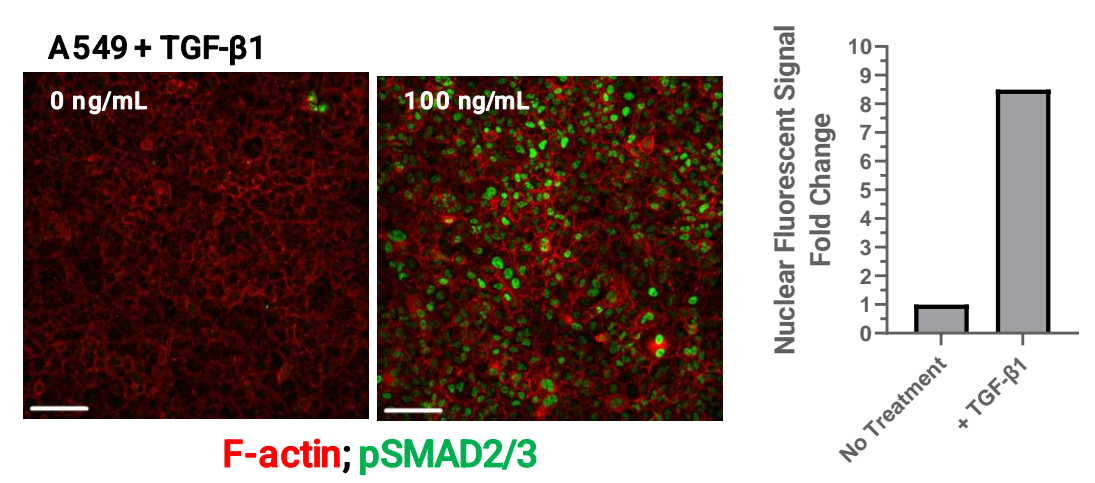
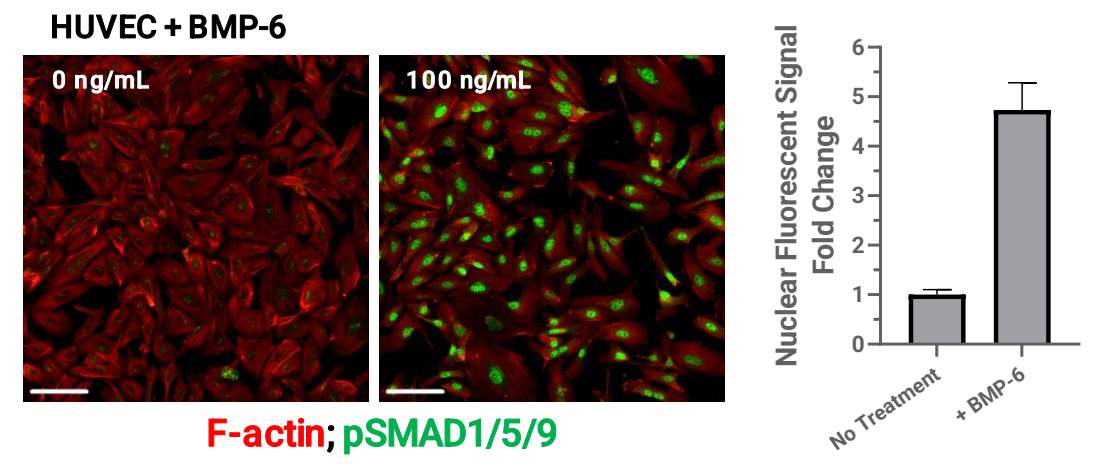
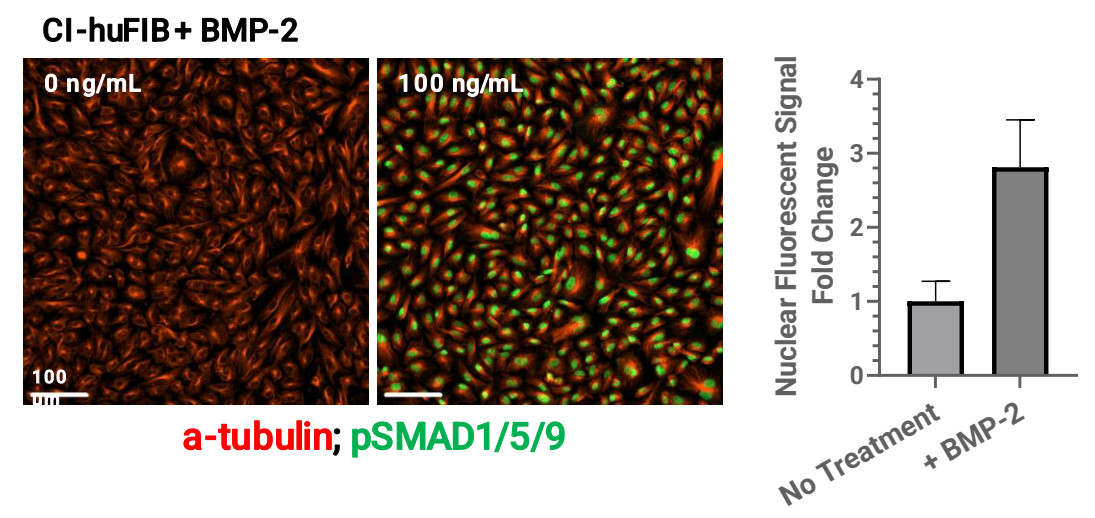
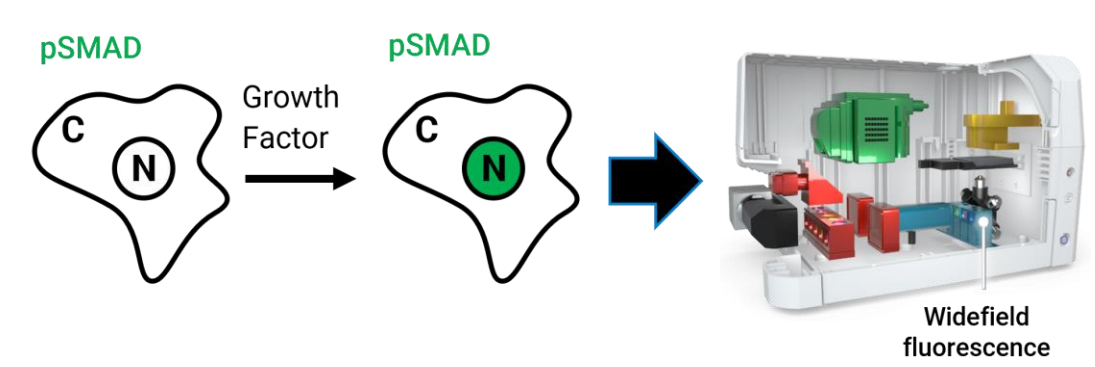
Cells. 2014, 3 (4), 981–993. doi: [10.3390/cells3040981](https://doi.org/10.3390/cells3040981)  
Attribution 4.0 International (CC BY 4.0)

## Results and Discussion

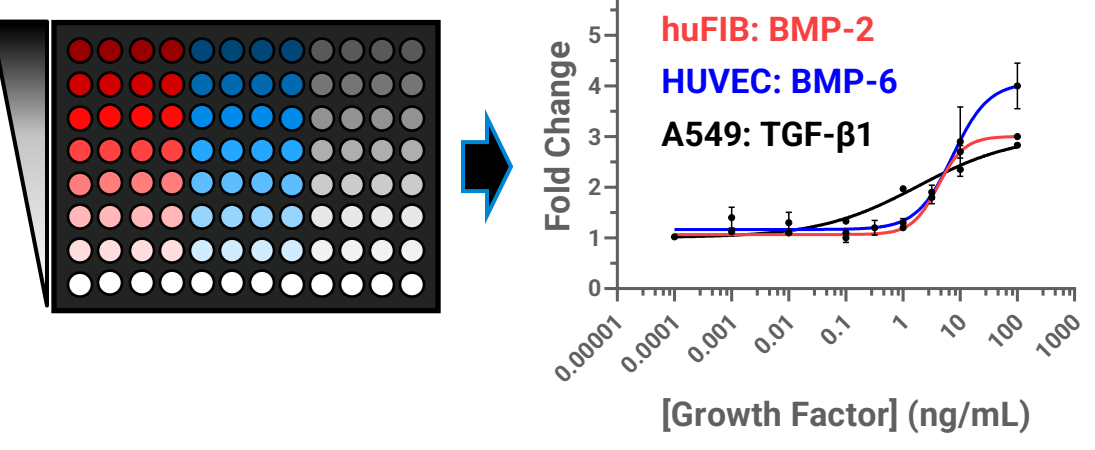
### Cellular assay: nuclear translocation

Cell Signaling Technology antibodies:

- Phospho-SMAD2 (S465/467) (E8R3R), part number #56532
- Phospho-SMAD1 (S463/465)/SMAD5 (S463/465)/SMAD9 (S465/467) (D3B10), part number 55605



### The nuclear translocation assay can be scaled to a high-throughput 96-well microplate array

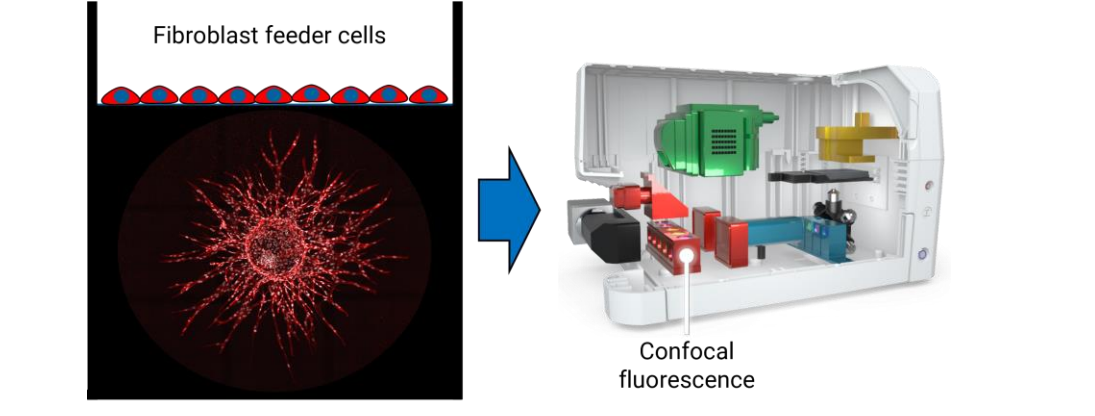


## Results and Discussion

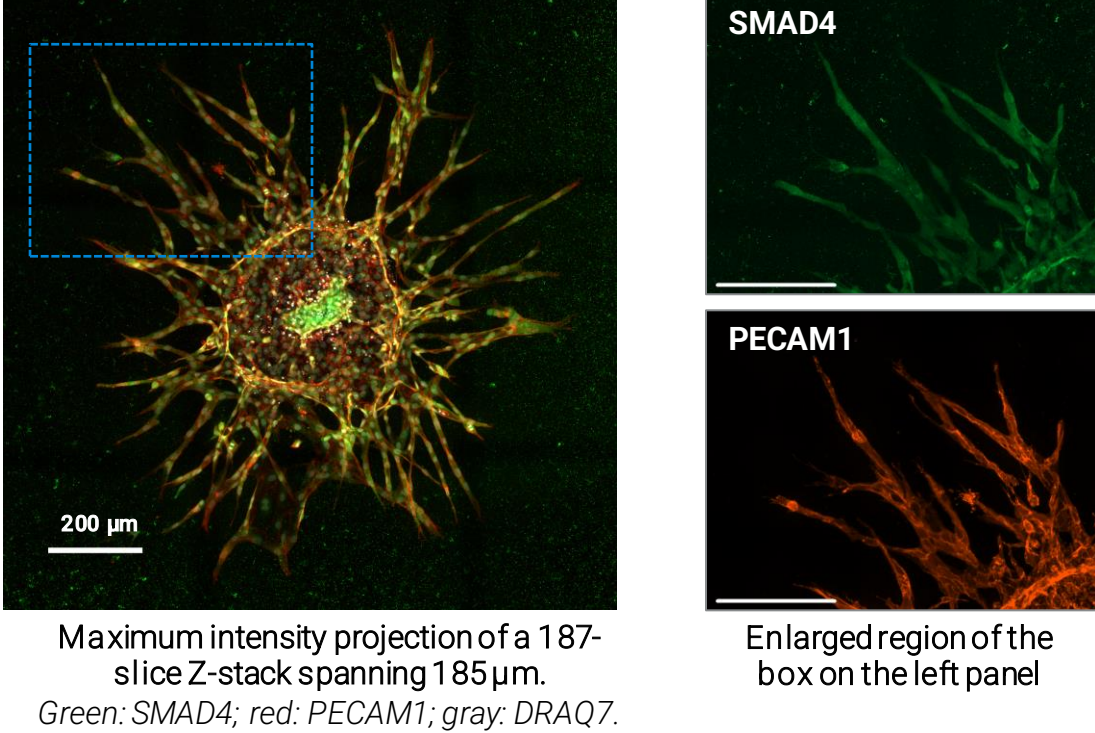
### 3D assay: spheroid sprouting angiogenesis, and EMT model

Cell Signaling Technology antibodies:

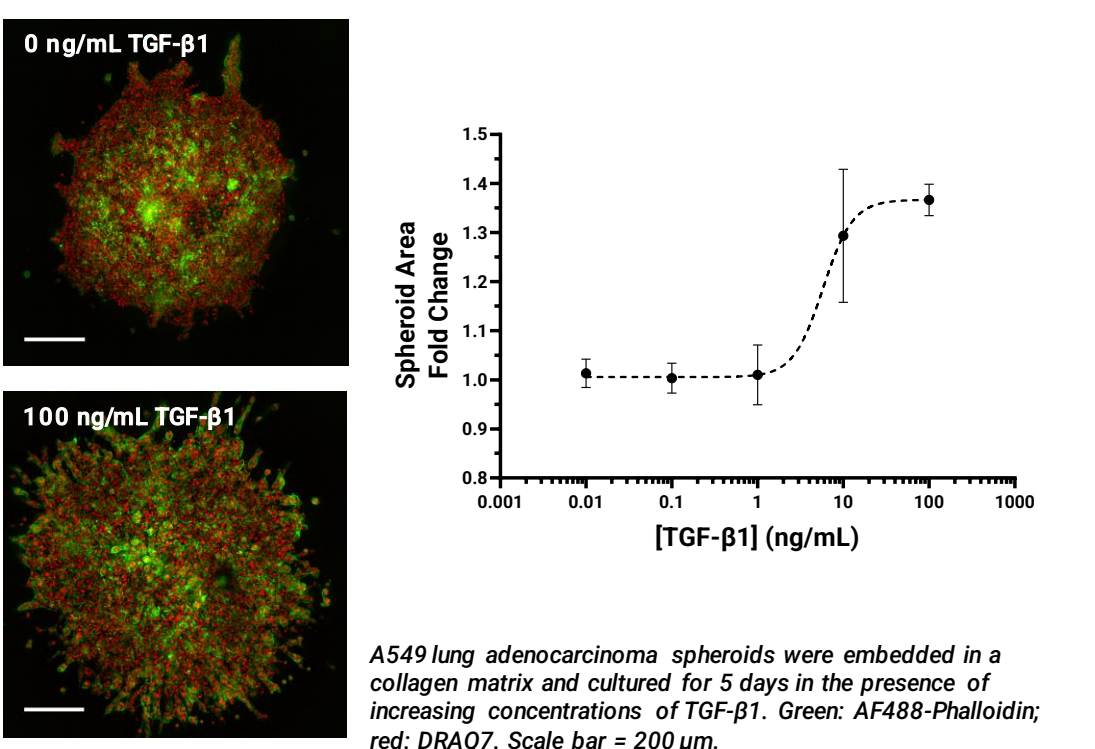
- SMAD4 (D3R4N), part number #24483



### CST reagents can be used for antibody-based detection of SMAD4 expression during in vitro angiogenesis development



### A549 lung adenocarcinoma spheroids serve as a 3D model for EMT when cultured in the presence of TGF- $\beta$ 1



## Conclusion

The high-quality reagents from Cell Signaling Technology combined with the multifunctional capabilities of the Agilent BioTek Cytation C10 confocal imaging reader enable quantitative evaluation of the TGF- $\beta$ /BMP pathway activation at different biological scales.

For Research Use Only. Not for use in diagnostic procedures. RA45026.4146412037