

Introduction

Epithelial-mesenchymal transition (EMT) is a cellular differentiation process whereby epithelial cells lose many epithelial features, while acquiring mesenchymal, fibroblast-like properties, leading to reduced cell-cell contacts and increased motility. While recognized as a fundamental process required for normal embryonic development, EMT is understood to be co-opted by malignant epithelial tumors to facilitate their metastatic spread. Central to stimulating EMT is the TGF- β superfamily of growth factor ligands, which elicit receptor-mediated responses in cells, primarily via TGF- β /SMAD signaling pathways. In these pathways, receptor-mediated SMAD (R-SMAD) proteins serve as the primary downstream effector molecules, the activities of which are regulated through receptor-mediated phosphorylation. The magnitude and duration of ligand-induced receptor activation influences the level of SMAD phosphorylation, which, in turn, influences the magnitude of the downstream cellular responses. Activation of the TGF- β -SMAD2/3 pathway is known to upregulate the expression levels of numerous EMT-associated genes, including *HMGA2*, *ZEB1*, *Snail1*, and *Slug*, all of which have furthermore been associated with cancer metastasis. In this study, we describe high-throughput methods to quantitatively evaluate the biochemical and cellular responses to TGF- β /SMAD pathway activation in a cellular model of TGF- β -induced EMT. Effects of pathway activation are examined at different levels of biological complexity—biochemical, cellular, and multicellular—using 2D and 3D (spheroid) models. Collectively, these approaches enable a comprehensive evaluation of TGF- β /SMAD pathway activation that is amenable to high-throughput analysis platforms.

Experimental

Cell lines

A549 lung adenocarcinoma cells were purchased from ATCC (part number CCL-185; Manassas, VA) and cultured in Advanced DMEM (A-DMEM, part number 12491; Gibco Thermo Fisher Scientific; Waltham, MA) containing 10% FBS and 1x penicillin/streptomycin/glutamine).

Growth factor treatment

Before TGF- β 1 exposure, adherent cells, or spheroids were subjected to an overnight, 18-hour serum-starvation period using A-DMEM without FBS. Human TGF- β 1 recombinant protein (part number 75362; Cell Signaling Technology, Danvers, MA) was diluted in serum-free media and added to cell and spheroid cultures at indicated concentrations. 2D cell cultures were treated with TGF- β 1 for 1 hour to measure SMAD protein phosphorylation responses, or 48 hours to measure expression changes in EMT-associated proteins. 3D spheroids were treated with TGF- β 1 for 4 days to enable spheroid outgrowth. Cells and spheroids were then fixed in 4% buffered paraformaldehyde for subsequent immunostaining.

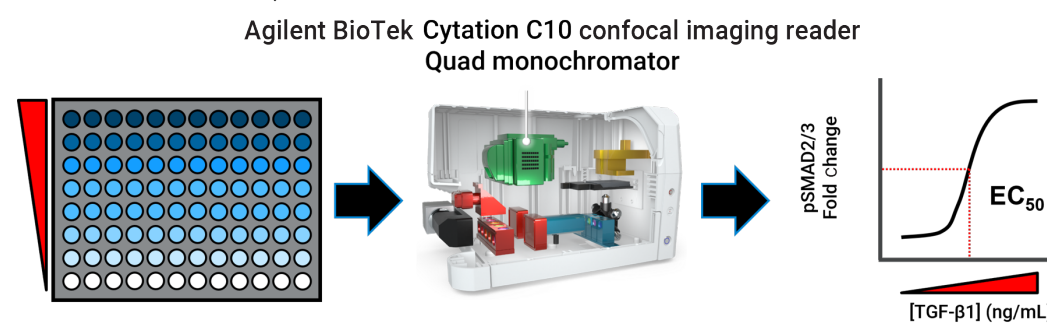
Imaging

Using the Agilent BioTek Cytation C10 confocal imaging reader equipped with a 60x 1.2 NA *water immersion* objective, a single Z-plane of adherent cells was imaged in three channels: GFP (F-actin; AF488-conjugated phalloidin), CY5 (phosphorylated SMAD2/3; CF633-conjugated goat anti-rabbit), and DAPI (nuclei; Hoechst 34580), with the DAPI channel used to set Z-focus height using laser autofocus. With the same three channels, multichannel Z-stacks of spheroids were imaged in confocal mode using the *deep-sectioning* 60 μ m disk and a 20x 0.4 NA objective. Multichannel Z-stacks underwent a background reduction transformation step before generating a maximum intensity project.

Results and Discussion

Biochemical analysis—high-throughput ELISA assay

- Cell Signaling Technology PathScan® Sandwich ELISA kits:
- Phosphorylated SMAD2 (S465/467)/SMAD3 (S423/425), #12001
 - Total SMAD2/3, #12000



SMAD2/3 phosphorylation in A549 cells

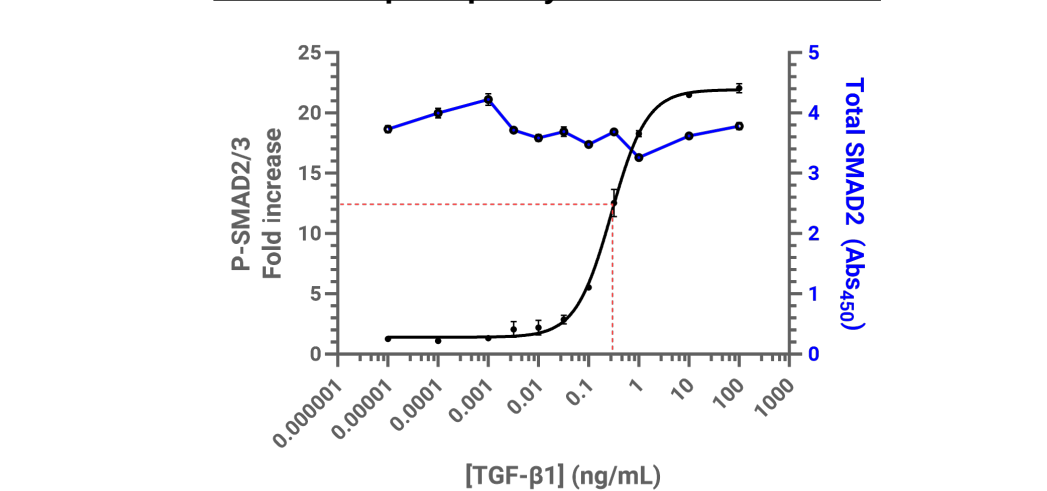


Figure 2. Dose-dependent SMAD2/3 phosphorylation (black) in response to TGF- β 1. The phosphorylated signal is normalized to total SMAD2 (blue).

Results and Discussion

Cellular assay—high-throughput nuclear translocation assay

Cell Signaling Technology antibodies:

- Phosphorylated SMAD2 (Ser465/Ser467) (E8R3R), #18338
- HMGA2 (D1A7) Rabbit mAb, #8179
- ZEB (E2G6Y) XP Rabbit mAb, #70512
- Slug (C18G7) Rabbit mAb, #9585

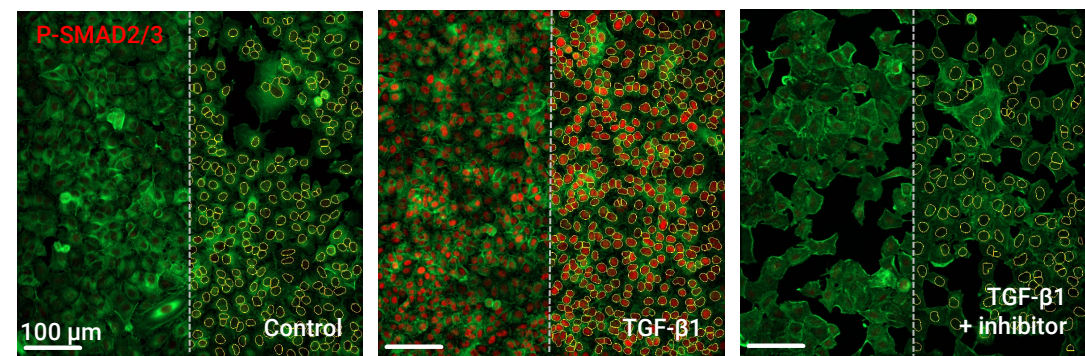
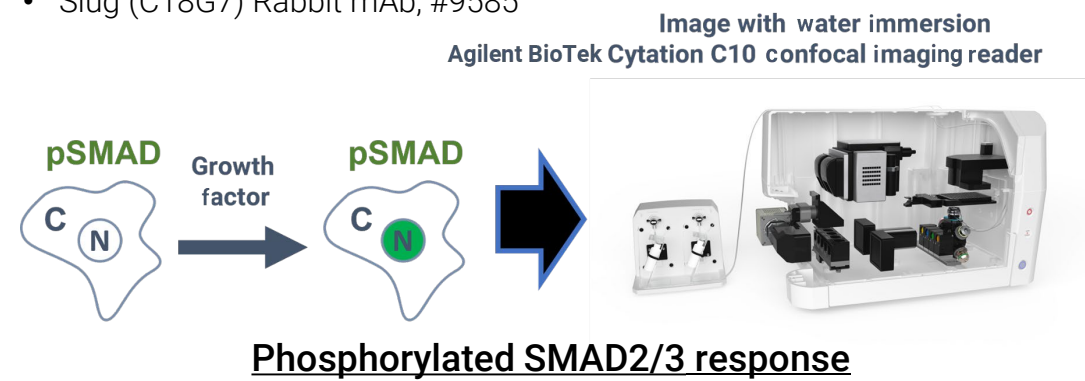


Figure 3A. TGF- β 1-induced SMAD2/3 phosphorylation and nuclear translocation, which is blocked by the TGF- β 1 receptor inhibitor SB431542. The right half of each image illustrates the image analysis carried out in Agilent BioTek Gen5 software, where phosphorylated SMAD2 fluorescence was quantified within each nuclei.

Changes in EMT-associated protein expression

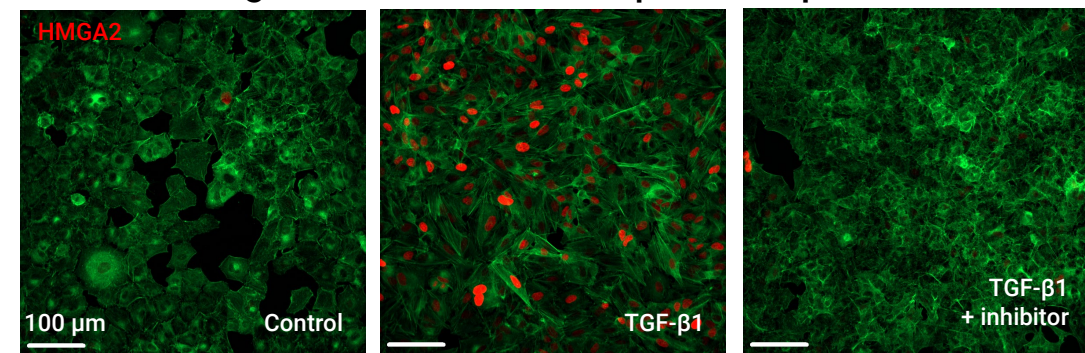


Figure 3B. Increased HMGA2 expression in response to TGF- β 1 treatment (48-h). Increased HMGA2 expression is blocked by the TGF- β 1 receptor inhibitor SB431542.

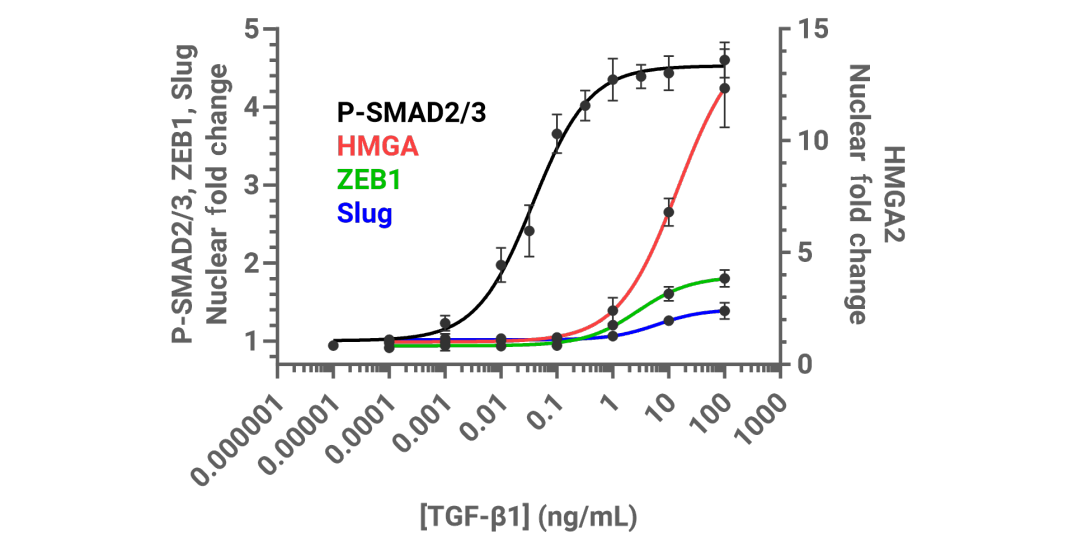


Figure 3C. Dose-response curves of TGF- β 1-induced nuclear expression of phosphorylated SMAD2/3 (P-SMAD2/3), HMGA2, ZEB1, and Slug.

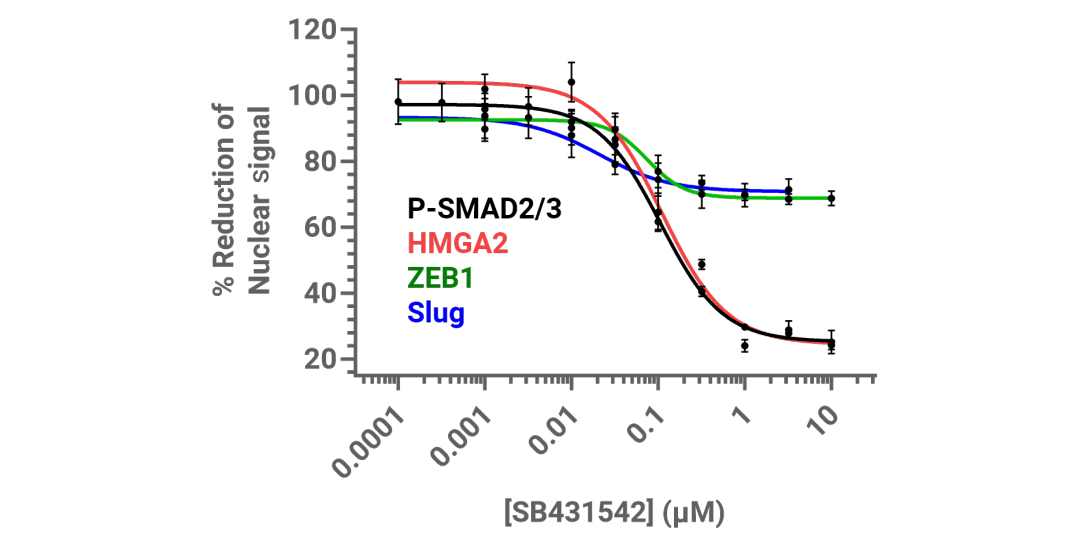


Figure 3D. Dose-dependent inhibition of TGF- β 1-induced SMAD2/3 phosphorylation, and expression of HMGA2, ZEB1, and Slug.

Results and Discussion

3D assay—spheroid model of EMT

Cell Signaling Technology antibodies:

- Phosphorylated SMAD2 (Ser465/Ser467) (E8R3R) Rabbit mAb, #18338
- SMAD4 (D3R4N) XP Rabbit mAb, #46535

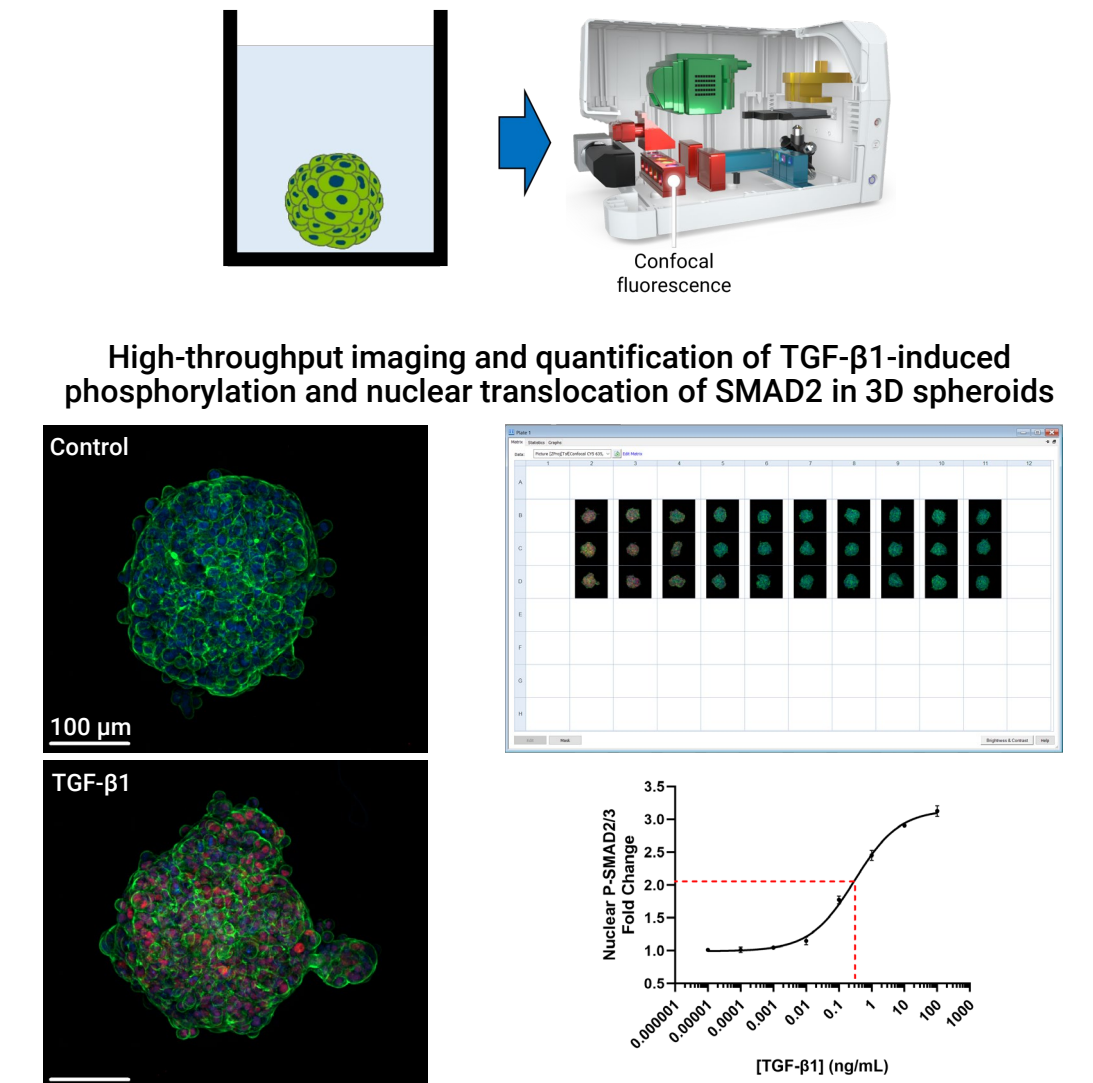


Figure 4. (Left) Maximum intensity projections of A549 spheroids treated with vehicle control, or 100 ng/mL TGF- β 1. Red: SMAD2. Green: F-actin, make unavailable. Nuclei. (Right, top) Plate layout of an experiment where triplicate spheroids were treated with decreasing amounts of TGF- β 1 (left to right). (Right, bottom) Quantified nuclear phosphorylated SMAD2 at each concentration assayed, with an EC₅₀ of 0.31 ng/mL.

Quantitative analysis of TGF- β 1-induced total SMAD4 protein nuclear translocation in 3D spheroids.

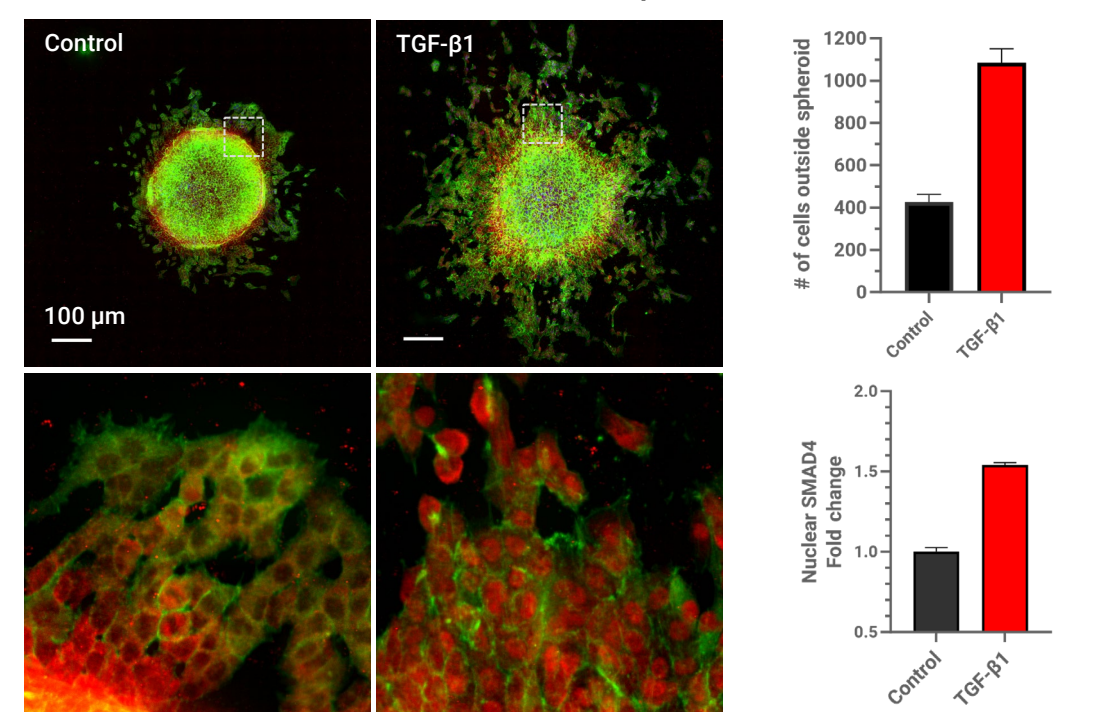


Figure 5. A549 spheroids were cultured on top of a collagen gel for 4 days in the presence/absence of TGF- β 1. Green: F-actin. Red: SMAD4. Bottom panels show an enlarged view of boxed regions in the top panels.

Conclusion

Recombinant monoclonal antibodies from Cell Signaling Technology combined with the multifunctional capabilities of the Agilent BioTek Cytation C10 confocal imaging reader enables quantitative evaluation of TGF- β 1-induced EMT pathway activation at different biological scales.

Canonical TGF- β /SMAD and BMP/SMAD signaling pathway

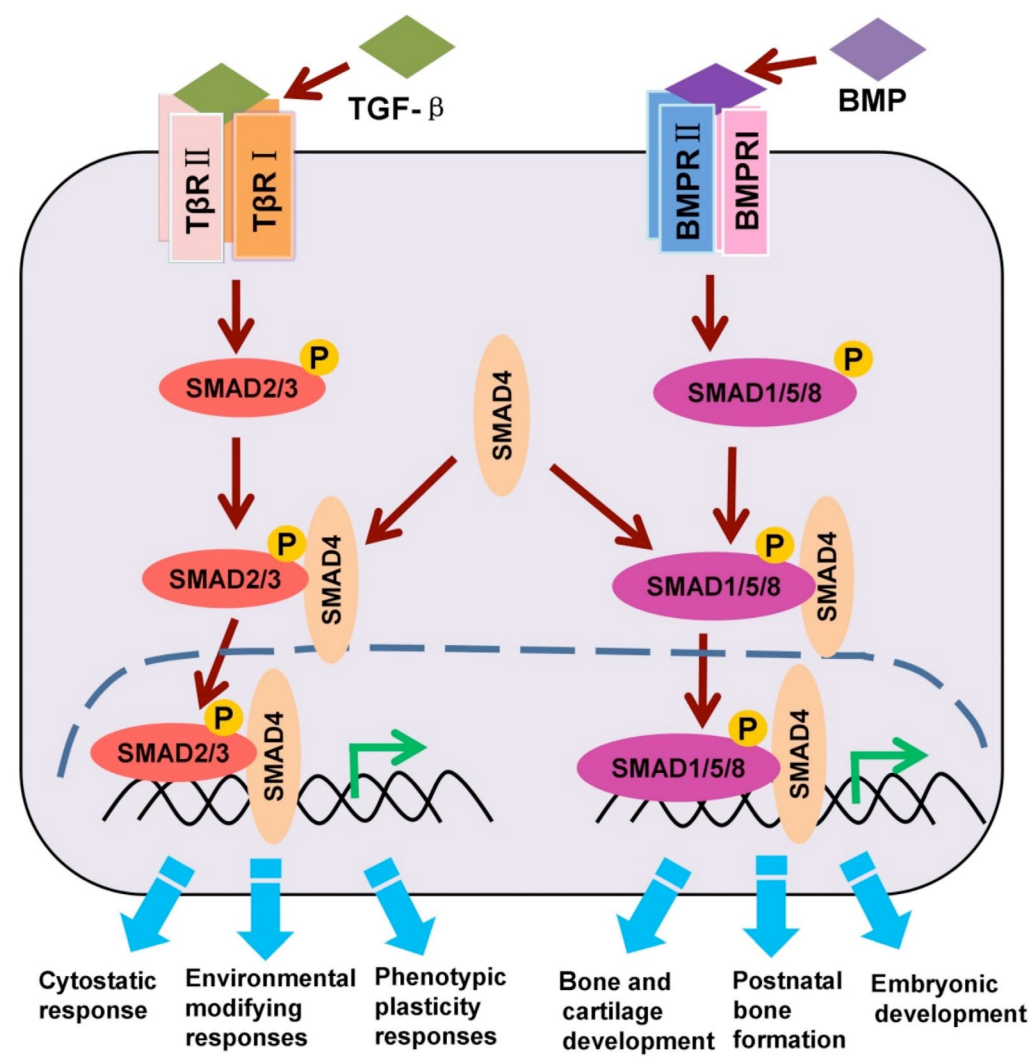


Figure 1. The canonical TGF- β /BMP/SMAD signaling pathway (adaptation). DOI: 10.3390/cells3040981