

Immunofluorescence of Collagen type IV (green) and nuclei (blue) after TGF-β stimulation

3D InSight™ Human Liver Fibrosis Model

The InSphero 3D InSight™ Human Liver Fibrosis Model is a primary human liver microtissue for *in vitro* screening of anti-fibrotic drug efficacy, and for studying mechanisms of fibrosis induction (Figure 1). The model includes the critical liver cells needed to recapitulate the fibrotic disease state upon induction, and serves as a powerful model for fibrosis drug discovery and development.

- **Rely on a physiologically relevant model** composed of primary human hepatocytes (PHHs), hepatic stellate cells (HSCs), Kupffer cells, and liver endothelial cells (LECs)
- **Expand your assay window** by leveraging a pre-qualified model that exhibits pro-fibrotic marker induction reflecting the *in vivo* disease state: ACTA2 (>5x) and COL1A1 (>3x)
- **Perform reproducible, throughput-compatible anti-fibrotic drug screening** using a standardized, assay-ready 96-well model system, amenable to a host of certified application endpoints

Certified Applications	
Designed for assessment of:	Options:
Efficacy screening of anti-fibrotic drugs	● ● ●
Testing pro-fibrotic potential of drugs	● ●
Anti-fibrotic drug mechanism of action	● ●
Combined efficacy and toxicity testing of anti-fibrotic drugs	● ●
Mechanism of disease progression	● ●
Contribution of cell types to disease progression	● ●
● Microtissue ● Service ● Protocol	

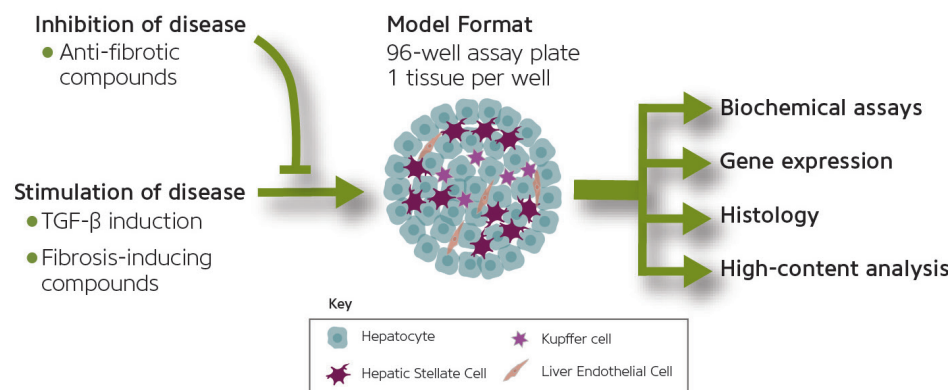


Figure 1: Disease Model Concept. The 3D InSight™ Human Liver Fibrosis Model enables the study of liver fibrosis induction and inhibition of disease progression. The screening-compatible 3D microtissue format provides maximal endpoint compatibility.

Liver Fibrosis Model Characterization Data

Induction of liver fibrosis requires the presence of HSCs, which are activated by pro-fibrotic stimuli, such as TGF-β. Kupffer cells (specialized macrophages located in the liver) and LECs influence liver fibrosis and are thus required for a physiologically, mechanistically relevant *in vitro* model system. The 3D InSight™ Human Liver Fibrosis model includes these relevant primary liver cell types, with robust induction of liver fibrosis biomarkers at the mRNA level observed upon TGF-β treatment (Figure 2). On histological examination, fibrotic biomarkers were also highly increased on the protein level (Figure 3). Concomitant treatment of TGF-β with an ALK5 inhibitor completely abolished progression of liver fibrosis (Figure 3), demonstrating suitability of the model system for testing pro- and anti-fibrotic compound efficacy.

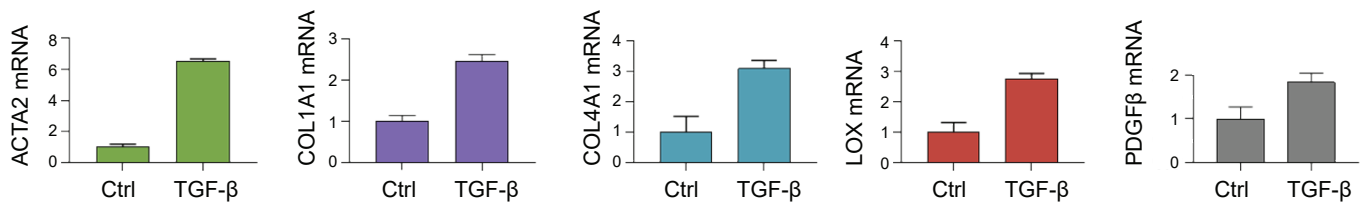


Figure 2: Robust induction of pro-fibrotic marker gene expression upon TGF-β stimulation. After TGF-β treatment, a 5-fold induction of ACTA2 (α-SMA, stellate-cell activation marker), 4-fold induction of COL1A1 and COL4A1 (ECM markers), 3-fold induction of LOX (early fibrosis marker), and 2-fold induction of PDGFβ (stellate-cell activation marker) is observed.

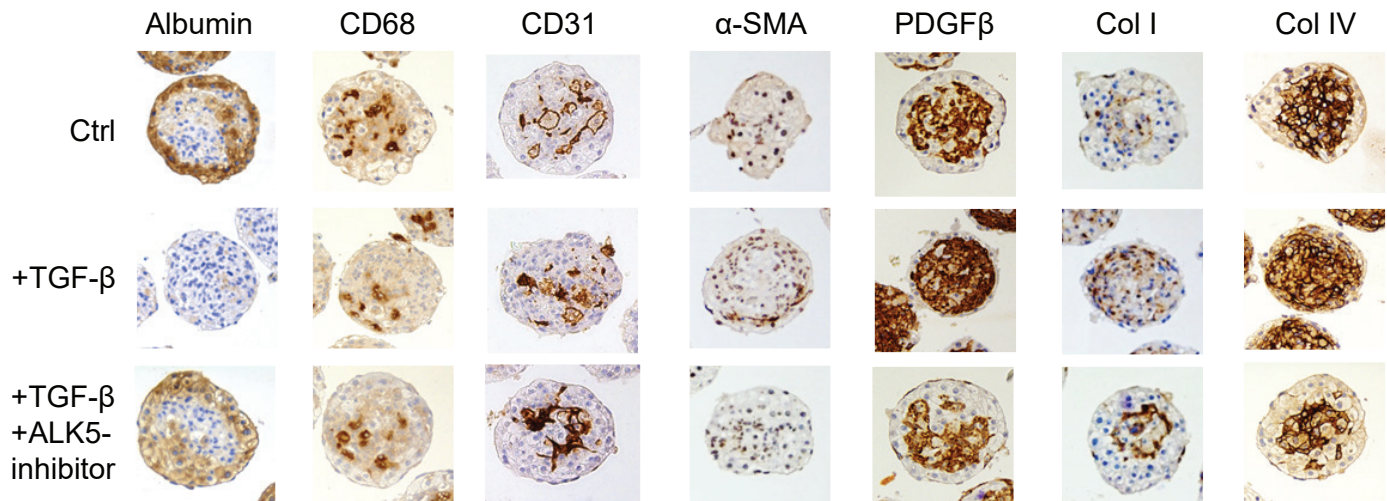


Figure 3: Histological characterization after TGF-β and disease inhibition. Control tissues exhibit presence of hepatocytes (albumin), Kupffer cells (CD68), LECs (CD31), and stellate cells (PDGFβ). TGF-β treatment induces elevated expression of α-SMA, PDGFβ, Col I, Col IV, accompanied by decreased hepatocyte function as detected by decreased albumin expression. Simultaneous treatment of TGF-β and ALK5 inhibitor halts the induction of fibrosis biomarkers and rescues hepatocyte function.

3D InSight™ Human Liver Fibrosis Model

Liver Fibrosis Disease model

Model Specifications

- **Microtissue**
3D InSight™ Human Liver Co-culture (PHHs, Kupffer cells, HSCs, and LECs)
- **Media**
InSight™ Human Liver Maintenance Medium
- **Format**
Akura™ 96 plate
96 microtissues, 1 tissue/plate

Related Models and Platforms

- **3D InSight™ Liver Disease Discovery Platform**
 - 3D InSight™ Human Liver NASH Model
 - 3D InSight™ Human Liver Steatosis Model
- **3D InSight™ Diabetes Discovery Platform**
 - 3D InSight™ Diabetes Type II Model
 - 3D InSight™ Diabetes Type I Model
- **3D InSight™ Liver Toxicology Platform**

IS-500-0029-01-Ad

InSphero AG
Schlieren, Switzerland
© +41 44 515 04 90

Email and Website
info@insphero.com
www.insphero.com

InSphero Inc.
Brunswick, ME, USA
© +1 800-779-7558

InSphero is ISO 9001:2015 certified
All rights reserved, © 2018 InSphero AG.
3D InSight and Akura are trademarks of InSphero AG. All other trademarks cited are the property of their respective owners. For life science research only. Not for use in diagnostic procedures.