

Join InSphero at SOT 2022

March 27-31, San Diego Convention Center

Tuesday, March 29

Exhibitor-hosted session | 12 – 1 PM | Room 23A

The X-Species DILI Validation Consortium: Progress and First Results

In partnership with: Pfizer, Sanofi, Merck, Genentech, AbbVie, Takeda, Philips Morris International, MicroMatrices



[Add to your conference agenda](#)

Speakers:

Prof. Dr. Armin Wolf, CSO, InSphero

Dr. Bruno Filippi, VP Liver Safety, InSphero

Dr. Björn Titz, Manager Proteomics and Lipidomics, Philip Morris International

Dr. Simon Plummer, Managing Director, MicroMatrices

Dr. Sue Grepper, Scientific Ambassador, InSphero – *Session Moderator*

Using animal models to predict human DILI has known issues and high failure rates. The [X-Species DILI Validation consortium](#) brings together representatives from biopharma to compare drug effects in regulatory animal models and human 3D *in vitro* liver models and recapitulate *in vivo* findings in animals by specific DILI mechanisms to improve the translations. Join the special scientific session to learn about the latest advancement within the pre-competitive consortium and the first findings. A light assortment of snacks and drinks will be provided, and we look forward to seeing you in the audience!



Poster Sessions

Wednesday, March 30



10:45 AM – 12:30 PM | Exhibit Hall | Session: **Liver II: In Vitro** | **4435/P435**

A Rapid 3D *In Vitro* Screening-Based Discovery Approach for Selecting and Prioritizing NASH Drug Candidates

Presenting Author: Dr. Sue Grepper, InSphero



2:30 PM – 4:15 PM | Exhibit Hall | Session: **Safety Assessment: Pharmaceutical-Drug Development II** | **4809/P874**

3D InSight™ Liver Microtissues as Promising Preclinical Safety Assessment *In Vitro* Tools: Mechanistic Investigation of Aflatoxin B1 Toxicity in Dog, Rat, and Human 3D InSight™ Liver Microtissues

Presenting Author: Dr. Bruno Filippi, InSphero



Top 10 abstracts at SOT 2022



2:30 PM – 4:15 PM | Exhibit Hall | Session: **Safety Assessment: Pharmaceutical-Drug Development II** | **4813/P878**

Mechanistic Investigation of Drug-Induced Liver Toxicity Using Human 3D InSight™ Liver Model

Presenting Author: Dr. Bruno Filippi, InSphero



2:30 PM – 4:15 PM | Exhibit Hall | Session: **Oxidative Injury and Redox Biology** | **4308/P242**

Relation between Oxidative Stress and Antioxidant Defense: Revisited in 3-Dimensional Primary Human Liver Microtissues

Presenting Author: Prof. Dr. Armin Wolf, InSphero

ToxExpo Booth 1126

Monday, March 28 – Wednesday, March 30

Meet InSphero Toxicology and 3D cell-based assay experts at SOT 2022 to learn how you can leverage tailored 3D *in vitro* solutions. Stop by our booth and get an exclusive look at our new all-in-one kit.

[Book a meeting](#) with an InSphero representative at SOT 2022.

We look forward to seeing you in San Diego!

**NEW
PRODUCT
LAUNCH**