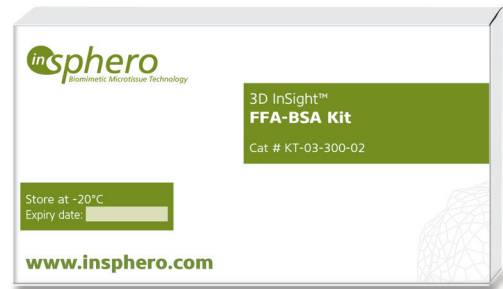


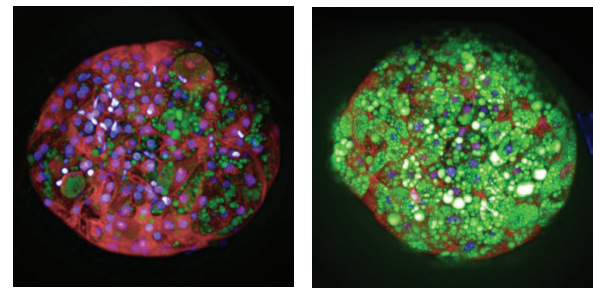
3D InSight™ FFA-BSA Kit



for liver steatosis induction with free-fatty acids

The 3D InSight™ FFA-BSA Kit contains conjugated, quality controlled, free fatty acids (FFA) bound to bovine serum albumin (BSA). Designed for *in vitro* studies with 3D InSight™ Human Liver Disease Discovery Platform models, this kit ensures efficient induction of steatosis to recapitulate the fatty liver disease state. The bioavailable, conjugated FFA substrates are readily taken up, metabolized and stored in primary human hepatocytes in our human liver microtissues models, enabling the study of non-alcoholic fatty liver disease (NAFLD) and other dietary-induced liver diseases.

- **Start running lipid loading immediately with ready-to-use substrates** that eliminate the need for the tedious, time-consuming variable conjugation process
- **Receive a comprehensive, 3D optimized protocol for steatosis induction** in Human Liver Microtissues and support from our expert application scientists
- **Effectively model non-alcoholic fatty liver disease (NAFLD), non-alcoholic steatohepatitis (NASH)** and other liver diseases using lipid-loaded human liver models.



Control

Fatty-acid-loaded

Induction of NAFLD in 3D InSight™ Human Liver Microtissues using the FFA-BSA kit. Immunofluorescence of lipids (green), nuclei (blue), membranes (red) in Human Liver Microtissues before and after treatment with free fatty acids. Stack imaging was performed on the PerkinElmer Opera Phenix™ High-Content Screening system.

3D InSight™ FFA-BSA Kit*

Kit Contents (enough to treat 4x96 microtissues)

- BSA (FFA-free) vehicle control
- BSA-coupled Oleate
- BSA-coupled Palmitate

Delivered at -20° C

*Available only as part of a 3D InSight™ Liver Disease Discovery Platform Bundle.

Related Models and Platforms

3D InSight™ Liver Disease Discovery Platform

- 3D InSight™ Human Liver NASH Model
- 3D InSight™ Human Liver Fibrosis Model
- 3D InSight™ Human Liver Steatosis Model

3D InSight™ Diabetes Discovery Platform

- 3D InSight™ Diabetes Type II Model
- 3D InSight™ Diabetes Type I Model