

Identification of drug candidates in MASH fast and cost-effectively with MASH CALL



Join the MASH CALL Initiative

Deadline September 6th

insphero

The Model of Excellence™

InSphero MASH CALL Initiative

Fast and Reliable **Identification** and **Comparison** of Clinical Candidates



Why engaging in a MASH CALL?

Fast and translational MASH phenotypic candidate selection process

Cost-effective human pre-clinical model

On demand combinatorial drug approach

Possibility to a comprehensive phenotypic analysis follow-up

Benchmarking with clinical candidates



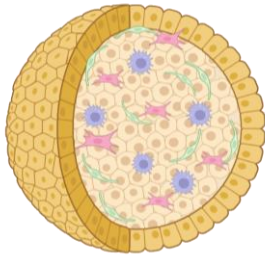
**PATH TO A SUCCESSFUL
MASH CLINICAL CANDIDATE**

How do you benefit? MASH CALL content

MASH CALL – 40 days since start until report delivery



Drug testing in Human Liver
Microtissue 3D InSight™
MASH model



- 6-concentration dose-response
- 3 replicates

ENDPOINTS FROM SAME ASSAY

Toxicity

Extracellular LDH Time course

Steatosis

Triglycerides

Fibrosis

Secreted Pro-Collagen 1 (2 time points)

CLINICAL AND TRANSLATIONAL BENCHMARKING

- Comparison against up to 20 clinical candidates
- Comparison against reference controls and healthy model

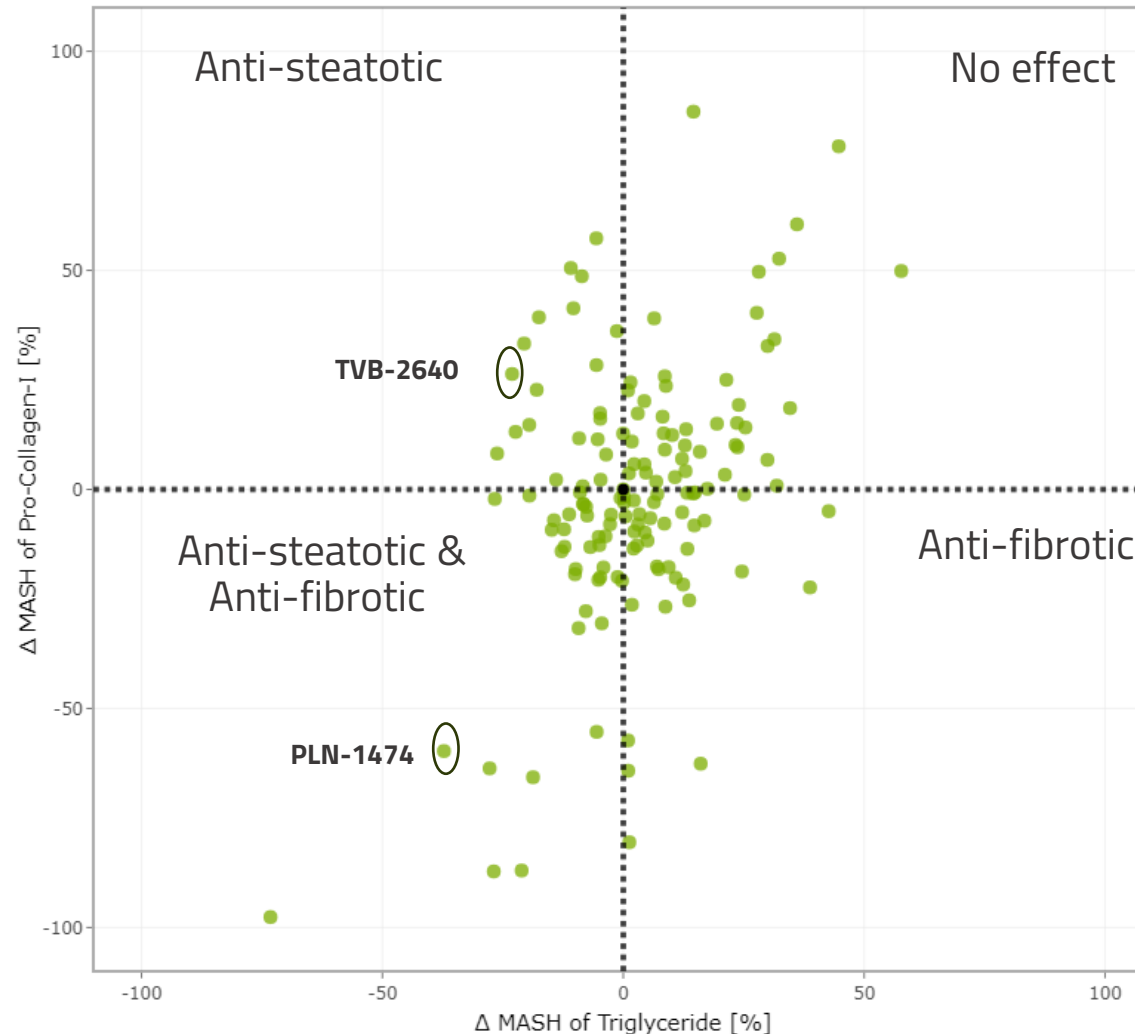
Anti-steatotic and anti-fibrotic comparison of 20 clinical candidates

Set up call for a deeper dive into the data!

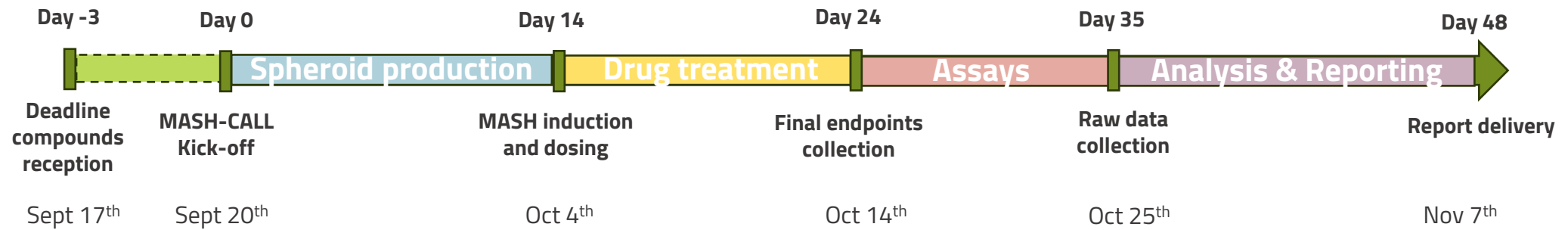
Examples of effects of clinical candidates matching MOA:

TVB-2640, a fatty-acid synthase enzyme inhibitor, has **anti-steatotic** effects

PLN-1474, a small-molecule selective inhibitor of the integrin $\alpha v \beta 1$, shows both **anti-steatotic and anti-fibrotic** effects



Timelines MASH CALL



Phase 1 MASH CALL and Phase 2 follow-up

Phase 1

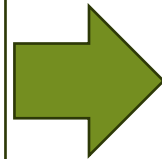
MASH CALL

Drug candidate screening and selection

With a 6-concentration dose-response range in 3 replicates

ENDPOINTS

- LDH
- Triglycerides
- Pro-Col1



Phase 2: Validation study and extended phenotypic analysis

- Increased sample size (n=6)
- Optimized concentrations (selected from phase 1)
- Validation of results
- Additional confirmation with comprehensive phenotypic analysis
- Flexibility of selection of endpoints (modularity) and time of execution

Pathophysiological Hallmarks	Offered Modular Assays
Steatosis	Total intracellular triglycerides
	Bodipy staining
Inflammation	Cytokine multiplex 6 markers (TNF- α , IL-6, IL-8, MCP-1, MIP-1 α , IP-10)
Fibrosis	Pro-collagen I (ELISA) – 2 time points
	Pro-collagen III (ELISA) – 2 time points
	MMP1, MMP2, MMP3, MMP9, TIMP1 (Luminex)
	Histology Sirius Red Staining
Samples collection	Collection of media supernatants for follow-up assays
	Pre-collection of samples for fibrosis quantification (Fibronest, Pharmanest)
	Pre-collection of samples for transcriptomics
Additional modules	Fibrosis quantification digital pathology (Fibronest, Pharmanest)
	Transcriptomics (TempO-Seq, DRUG-Seq, low-input RNA-Seq)



THANK YOU

Francisco Verdeguer, PhD
Liver Discovery

Radina Kostadinova, PhD
Senior Scientist, Deputy VP Liver Discovery

Jesús Glaus, PhD
Senior Scientist Liver Discovery

