

# C-DILI™ Assay: Integrating BSEP Inhibition and FXR Antagonism to Improve Prediction of Cholestatic Drug Induced Liver Injury

Jonathan P. Jackson, Kimberly M. Freeman, Matthew K. Palmer, Robert L. St. Claire III, Christopher B. Black, and Kenneth R. Brouwer  
BioIVT, ADME-TOX, 2810 Meridian Parkway, Suite 100, Durham, NC 27713

## INTRODUCTION

Cholestatic Drug Induced Liver Injury (DILI) in humans has been associated with bile salt export pump (BSEP) inhibition ( $IC_{50}s < 25 \mu M$ ); however, *in vitro* BSEP  $IC_{50}$  concentrations do not correlate with *in vivo* cholestatic DILI severity. Sandwich-cultured human hepatocytes (SCHH), when treated with BSEP inhibitors, respond to the resulting increased intracellular concentrations of bile acids (BA), via activation of FXR (adaptive response). This results in decreased synthesis of BA and increased expression of basolateral and canalicular efflux of BA via OST $\alpha/\beta$ , and BSEP which prevents cholestatic hepatotoxicity. We propose that BSEP inhibition alone may not be sufficient to induce toxicity due to this adaptive response. In addition to BSEP inhibition, inhibition of basolateral efflux and/or interference with the adaptive response (FXR antagonism) may lead to increases in drug-induced cholestatic bile acid hepatotoxicity. Such mechanisms must be incorporated to accurately predict *in-vivo* cholestatic drug induced liver injury (DILI).

We have determined the time dependence of this adaptive response, linking inhibition of BSEP, increases in the intracellular concentration of BA, activation of FXR, and evaluated the effects of FXR antagonists. A predictive model for cholestatic hepatotoxicity was developed (C-DILI™ Assay) which integrates the effects of BSEP inhibition, basolateral efflux inhibition, and FXR antagonism using a human hepatocyte system.

## METHODS

**Human Hepatocytes** Cryopreserved, Transporter Certified™ human hepatocytes in a sandwich configuration were cultured using QualGro™ Media for 5 days.

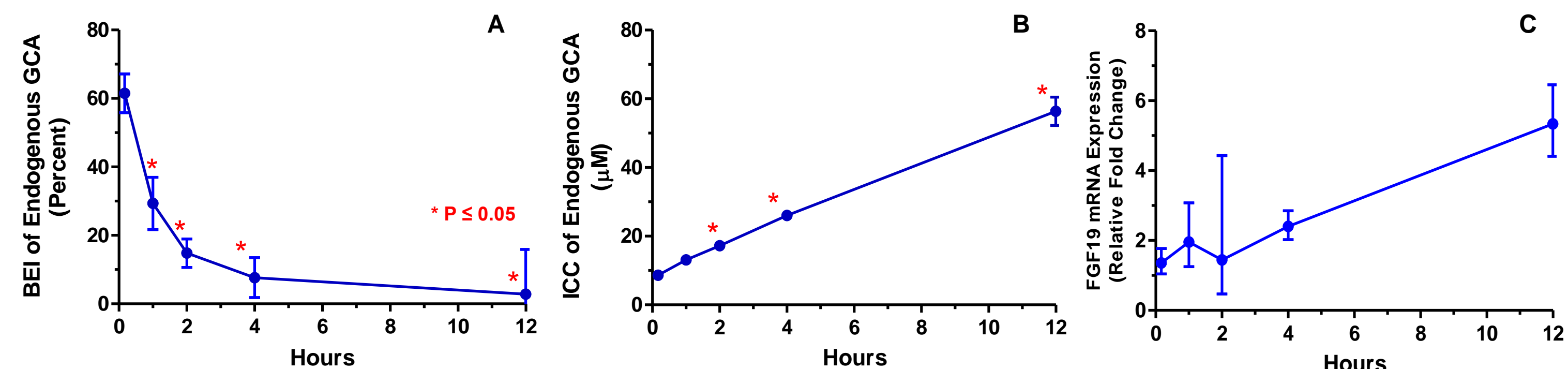
**Treatments** On Day 4 of culture, hepatocyte cultures were exposed to test compounds (Table 1) at 20X (or limit of solubility) of their systemic  $C_{max}$  in Qual-Gro™ Sensitization Media (physiologically relevant concentrations of lipids and bile acids) for 24 hours. Approximately 49 compounds with a range of BSEP  $IC_{50}s$  and associated with clinical DILI were tested. Treatment with vehicle control and standard QualGro™ Media was used as a control.

**Gene Expression** mRNA content of various transporters, synthetic enzymes, and regulatory factors from SCHH was determined from each RT reaction using gene-specific TaqMan® primer/probe sets. All reactions were normalized to the endogenous control GAPDH. Amplifications were performed on an ABI ViiA7 Real-Time PCR System in relative quantification mode. Relative-fold mRNA content was determined for each treatment group relative to the vehicle control.

**Endogenous Bile Acids** LC-MS/MS which employed reversed-phase HPLC and electrospray ionization was used to quantitate endogenously generated cholic acid (CA), CDCA, and their taurine (TCA, TCDCA) and glycine (GCA, GCDCA) conjugates in cells, bile and cell culture media.

## RESULTS AND DISCUSSION

Figure 1



Exposure to Cyclosporine A (10  $\mu M$ ), a potent BSEP inhibitor leads to a **rapid, time dependent decrease** in biliary excretion of endogenous bile acids.

Inhibition of biliary excretion leads to an increase in the intracellular concentration of endogenous bile acids.

Increased intracellular concentrations of bile acids activate FXR (increased FGF19).

Figure 2 Increased OST $\beta$  expression through activation of FXR (adaptive response) observed:

- in the presence of a BSEP inhibitor (CsA)
- Increased intracellular concentration of bile acids by addition of CDCA

Synergistic effect observed in the presence of CsA and CDCA due to greatly increased intracellular concentrations of bile acids

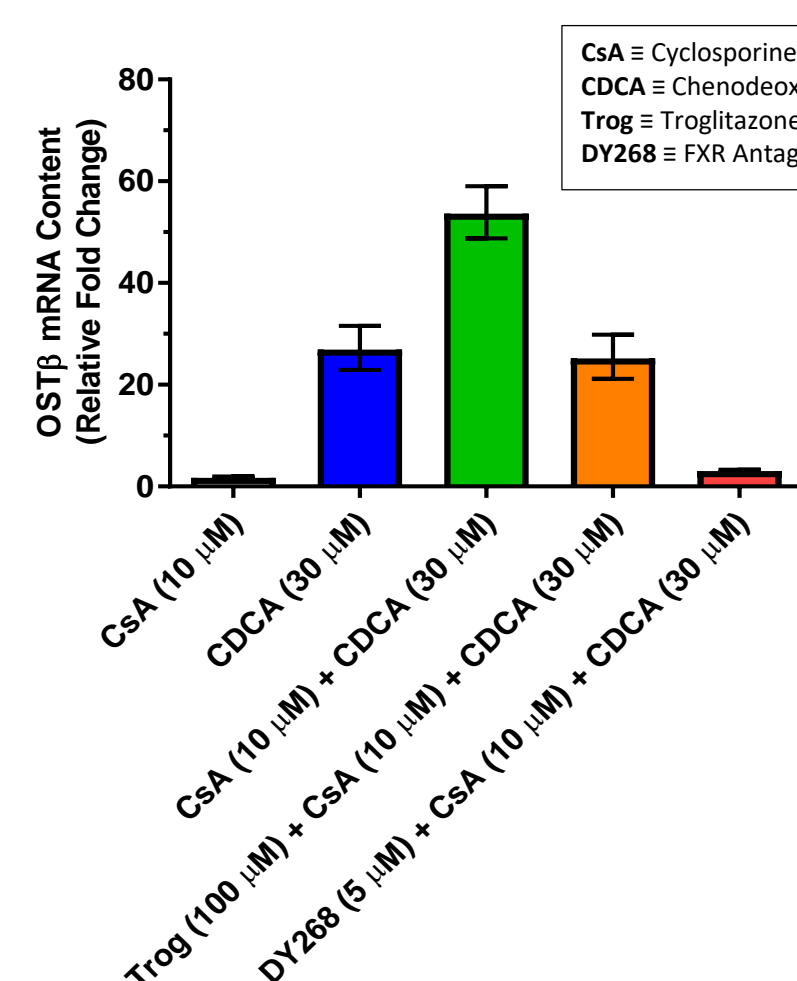


Figure 3 Decrease in OST $\beta$  mRNA expression when troglitazone (weak FXR antagonist) or DY268 (strong FXR antagonist) are co-administered with CsA and CDCA

FXR antagonism can prevent the hepatocyte from responding to high intracellular concentrations of bile acids and increasing the potential for cholestatic hepatotoxicity

Increase in Canalicular and Basolateral Efflux Clearance Following Exposure to CDCA

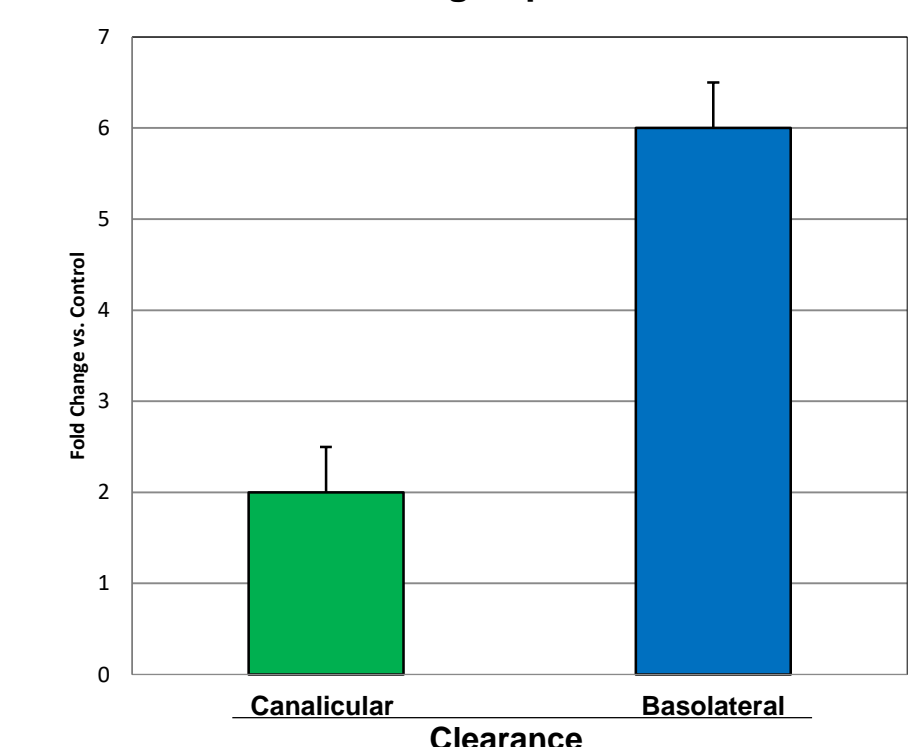


Figure 4 Chronic exposure to CDCA increases mRNA expression of BSEP (canalicular efflux transporter) and OST $\alpha/\beta$  (basolateral efflux transporters) by approximately 8X and 80X, respectively<sup>2</sup>

Molecular modeling was utilized to estimate changes in the canalicular and basolateral efflux clearances following 72 hours of exposure to CDCA (100  $\mu M$ ). Consistent with the changes in mRNA, the canalicular clearance increased by approximately 2X while the basolateral clearance increased by 6X.

In the induced state basolateral efflux clearance mediated by OST $\alpha/\beta$  represents an important elimination pathway.

**C-DILI™ Assay** The cultures were measured for ATP content (CellTiter-Glo® Promega) and LDH secretion (CytoTox-ONE™ Promega) following 24 hours of treatment. The LiverTox data base was used to identify compounds that were consistent with hepatocellular injury. LDH readout was used as a surrogate marker of cholestatic bile acid toxicity.

Table 1 Compounds were from selected publications<sup>1,3</sup> based on BSEP  $IC_{50}$  values, and hepatotoxicity information from the NIH LiverTox<sup>4</sup> database.

Drug	DILI Category	C-DILI™ Assay Potential	Correct Call	Clinical DILI
Acitretin	Idiosyncratic	Low	True negative	Yes
Amiodarone	Steatotic	High	False positive	Yes
Atorvastatin	Idiosyncratic	Low	True negative	No
Bosentan	Idiosyncratic	Low	True negative	Yes
Calcitriol	Idiosyncratic	Low	True negative	NID <sup>3</sup>
CsA	Lack of causal evidence	Low	True negative	No
Clofibrate	Cholelithiasis	Low	True negative	Yes
Deferasirox	DILI	High	True positive	Yes
Dicloxacillin	Idiosyncratic	Low	True negative	Yes
Dipyridamole	Lack of causal evidence	Low	True negative	No
Donepezil	Lack of causal evidence	Low	True negative	No
Entacapone	Lack of causal evidence	Low	True negative	No
Eryth. Estolate	Idiosyncratic	Low	True negative	Yes
Everolimus	Lack of causal evidence	Low	True negative	No
Felbitimbe	Lack of causal evidence	Low	True negative	No
Felxostat	Lack of causal evidence	Low	True negative	No
Flupirtine	NID	Low	---	NID <sup>3</sup>
Fluvastatin	DILI	High	True positive	Yes
Glyburide	Lack of causal evidence	Low	True negative	No
Iloperidone	Lack of causal evidence	Low	True negative	No
Imatinib	DILI	High	True positive	Yes
Indomethacin	Idiosyncratic	Low	True negative	Yes
Iribesant	Idiosyncratic	Low	True negative	Yes
Ketoconazole	DILI	High	True positive	Yes
Lapatinib	DILI	Low	False negative	Yes
Losartan	Idiosyncratic	Low	True negative	Yes
Megestrol Acetate	NID	Low	---	NID <sup>3</sup>
Mifepristone	NID	Low	---	NID <sup>3</sup>
Nicardipine	Lack of causal evidence	Low	True negative	No
Nifedipine	Idiosyncratic	Low	True negative	Yes
Pazopanib	DILI	Low	False negative	Yes
Pioglitazone	Lack of causal evidence	Low	True negative	No
Posaconazole	DILI	High	True positive	Yes
Pranlukast	NID	Low	---	NID <sup>3</sup>
Primaquine	No clinical evidence	Low	True negative	No
Repaglinide	Idiosyncratic	Low	True negative	Yes
Risperone	Lack of causal evidence	Low	True negative	No
Rifabutin	Lack of causal evidence	Low	True negative	No
Ritonavir	Difficult to separated from viral infection	High	True positive	Yes
Rosiglitazone	Lack of causal evidence	Low	True negative	No
Simvastatin	Idiosyncratic	Low	True negative	Yes
Sitaxsentan	Idiosyncratic	Low	True negative	Yes
Tacrolimus	Lack of causal evidence	Low	True negative	No
Telithromycin	Idiosyncratic	Low	True negative	Yes
Telmisartan	Lack of causal evidence	Low	True negative	No
Tolcapone	DILI	High	True positive	Yes
Tolvaptan	Observed in cirrhosis and ADPKD patients	Low	True negative	Yes
Troglitazone	DILI	High	True positive	Yes
Zafirlukast	Idiosyncratic	Low	True negative	Yes

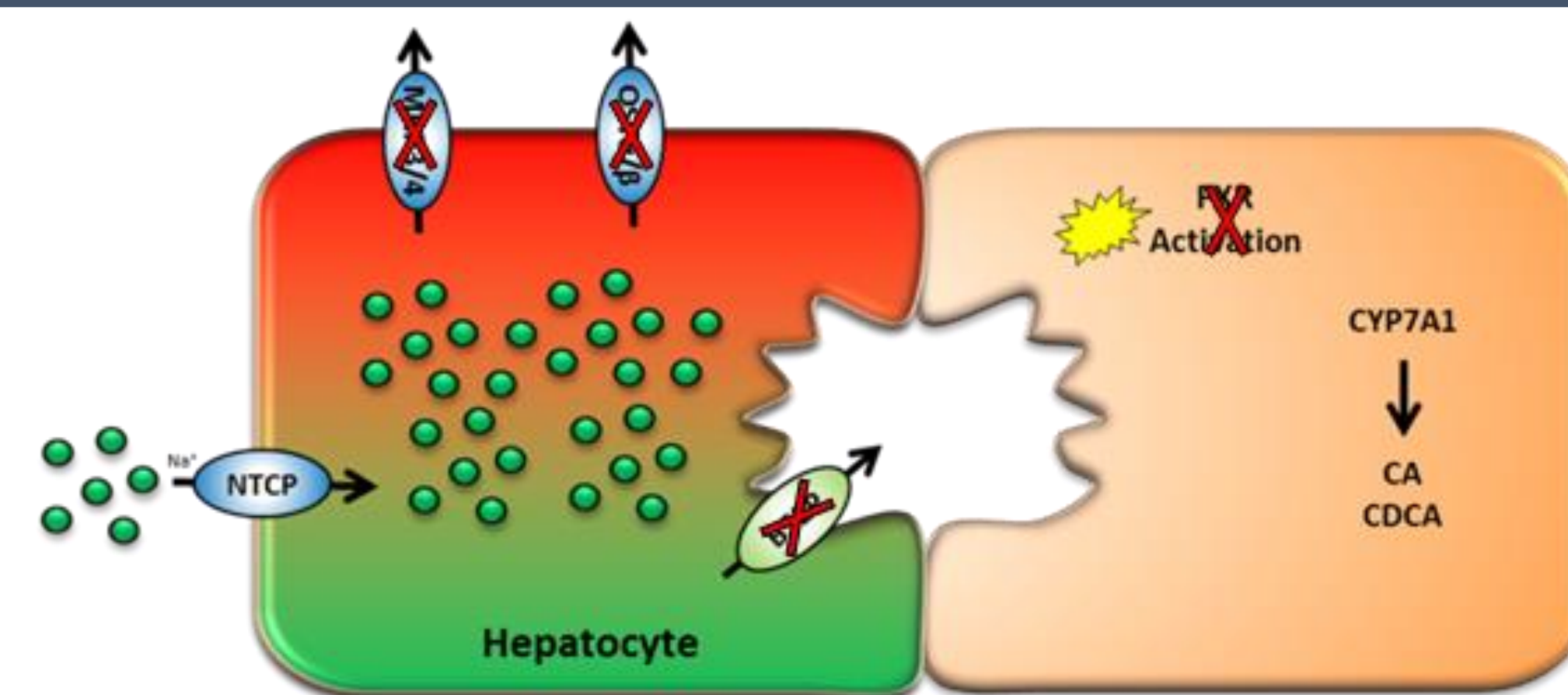


Figure 5 Inhibition of BSEP along with antagonism of FXR activation in response to increased intracellular concentrations of bile acids, and/or inhibition of bile acid basolateral efflux transporters can lead to increased cholestatic hepatotoxicity.

Drugs that impact multiple pathways or single pathways when given in combination with other drugs that affect alternate pathways could also result in hepatotoxicity. E.g. drug interaction with a BSEP inhibitor and FXR antagonist or basolateral efflux transporter inhibitor.

Table 2 The C-DILI™ Assay correctly predicted hBSEP False Positives (Pioglitazone, Rosiglitazone) as having low potential for hepatocellular cholestasis, and hBSEP False Negative (Deferasirox) as having a high potential. Troglitazone and Ketoconazole, both True Positives in the hBSEP analysis were also correctly predicted in the C-DILI™ Assay. Cyclosporine A is a potent BSEP inhibitor, however the hepatocyte can adapt through FXR activation of basolateral efflux transporters, and is used as a negative control in the C-DILI Assay. Troglitazone is a potent BSEP inhibitor (sulfate metabolite) and a weak FXR antagonist, and recently the sulfate metabolite has been shown to inhibit OST $\alpha/\beta$ <sup>5</sup>. Troglitazone is used as a positive control in the C-DILI assay.

Compound	hBSEP $IC_{50}$ ( $\mu M$ )	hBSEP Prediction	Correct Prediction <sup>4</sup>	Reports of Liver Injury	C-DILI™ Prediction
Cyclosporine A	0.5	Positive	Negative	None or rare	Low Potential
Pioglitazone	0.5	Positive	Negative	None or rare	Low Potential
Rosiglitazone	3	Positive	Negative	None or rare	Low Potential
Troglitazone	3	Positive	Positive	Hepatocellular	High Potential
Ketoconazole	3	Positive	Positive	Hepatocellular	High Potential
Simvastatin	25	Positive	Negative	None or rare	Low Potential
Imatinib	25	Positive	Negative	Hepatocellular	Direct Toxicity
Fluvastatin	36	Negative	Positive	Mixed or Rare	Moderate
Deferasirox	58	Negative	Positive	Mixed	High Potential

Compared to use of hBSEP data alone, analysis with C-DILI™ increased the accuracy from 22% to 96%.

## Applications:

### Discovery Stage

- No information on clinical concentrations
- Screen at high concentrations (50 – 100  $\mu M$ ), and then follow up hits with a dose ranging study at lower concentrations

### Pre-Clinical Stage

- Projected clinical concentrations
- Screen at concentrations that cover clinical  $C_{max}$  or  $C_{ss}$  and up to 20X to 50X to account for higher portal vein concentrations

### Clinical Stage

- Known clinical concentrations
- Screen for potential drug interactions at 20X clinical  $C_{max}$  or  $C_{ss}$  for test compound and anticipated concentration range for co-administered compound

## CONCLUSIONS

The C-DILI™ Assay correctly predicted compounds with significant clinical hepatocellular cholestatic toxicity

- Integrates effects on BSEP, OST $\alpha/\beta$ , MRP3/4, and FXR to delineate hepatotoxicity resulting from a build up of intracellular bile acids.

### Key Assay Features:

- Transporter Certified™ human hepatocytes in a model that maintains transporter expression, localization and function
- Optimized culture media
- Convenient 24 hour incubation in 96-well format
- Assess effects of parent and metabolites simultaneously
- Integrated effects of transporter inhibition (acute effects) and regulation (chronic effects) in a single toxicity readout
- Evaluate cholestatic vs. direct toxicity
- Available as a service or kit

<sup>1</sup> Dawson et.al., Drug Metab Dispos 40,130, 2012, <sup>2</sup>Jackson et.al., App In Vitro Tox 2, 4, 2016, <sup>3</sup>Morgan et.al. Tox Sci 136,1, 2013, <sup>4</sup><https://livertox.nih.gov/>

<sup>5</sup>Malinen et.al. Am J Physio – GI & Liver Physio. Feb 2018