

Accumulation of taurocholic acid in B-CLEAR® human sandwich-cultured hepatocytes in the presence of clinically relevant cholestatic agents

Caleb D. Istringhausen,¹ Lydia M. Vermeer,¹ Kimberly Freeman,² Chris Black,² Kenneth D. Brouwer,² and Greg J. Loewen¹

¹XenoTech, LLC, 16825 W. 116th St., Lenexa, KS, USA ²Qualyst Transporter Solutions, LLC, 2810 Meridian Parkway, Durham, NC, USA

Abstract

Purpose: Drug induced liver injury (DILI) has led to the withdrawal of drugs from the market. One postulated mechanism of DILI is cholestasis caused by inhibition of the bile salt efflux pump (BSEP). Inhibition of BSEP can be measured *in vitro* with BSEP expressing vesicles. However, not all drugs identified as BSEP inhibitors cause DILI as there are many pathways involved in cholestatic toxicity. An *in vitro* system which models *in vivo* transporter expression and functionality more completely may better predict the potential of drugs to cause cholestasis or DILI. We evaluated the hepatobiliary disposition of taurocholic acid (TCA) in the presence of cholestatic agents using B-CLEAR® human sandwich-cultured hepatocytes.

Methods: Compounds were screened for effect on uptake, efflux, and intracellular concentrations of [³H]-TCA using B-CLEAR® human sandwich-cultured hepatocytes in the presence of 4% BSA. Cells were incubated with inhibitor for 10 min and then co-incubated with inhibitor and TCA for 30 min. Follow up work examined a range of concentrations for compounds that modified the intracellular concentration of TCA, namely: of fluvastatin, ezetimibe, cyclosporine A, nicardipine, and cisapride monohydrate. In all assays, cells were lysed with 0.1% NaOH and analyzed using LSC.

Results: Results of screening assays are described in the table. The biliary excretion index (BEI), intracellular concentration (ICC), and Kp (ratio of ICC to extracellular concentration) of TCA are important factors in determining how exposure to a compound effects hepatobiliary disposition. A higher ICC and Kp value may indicate inhibition of TCA efflux, and a higher potential for toxicity, while a lower ICC and Kp value may indicate inhibition of uptake.

Conclusion: Transporter based interactions with bile acids are complex and an integrated system is required to evaluate uptake and efflux inhibition of bile acids. Studying the effect of a compound on the ICC and Kp ratio for TCA may provide information on the integrated effects of inhibition of uptake and efflux by a xenobiotic. Overall, the use of an *in vitro* system which more closely models *in vivo* conditions can lead to better predictions of hepatic transporter inhibition and toxicity.

Background & Purpose

Drug-induced liver injury (DILI) is characterized as liver injury due to intake of medications or xenobiotics, which leads to liver abnormalities or dysfunction.¹ There are a large number of drugs that have been withdrawn from the market due to acute DILI; including, troglitazone (antidiabetic and anti-inflammatory), benzbramarone (gout), and sitaxsentan (pulmonary arterial hypertension).² The mechanism behind DILI is multifaceted and can be exacerbated by unpredictable metabolism and bodily response, as well as the complex relationship between an individual's genetic makeup and environmental risk factors.³ Furthermore, several factors may contribute to hepatocellular injury and cholestasis; including, acute hepatocyte necrosis, the production of reactive metabolites (and oxidative stress) during cytochrome P450 metabolism of the parent drug, activation of stress signaling, and mitochondrial dysfunction.⁴

Bile acids are water soluble end products of cholesterol metabolism which are highly regulated by metabolism, excretion, absorption, and feedback mechanisms, in order to limit their intracellular accumulation. It is hypothesized that dysfunction to the bile salt efflux pump (BSEP) may contribute to the mechanism of action behind DILI.⁵ Current preclinical, *in vitro* models of BSEP inhibition may not accurately predict the potential for DILI. Therefore, an *in vitro* model which better reflects *in vivo* transporter expression and activity may better predict the potential for novel xenobiotics to cause DILI in the clinic. The purpose of this work was to evaluate the use of B-CLEAR® human sandwich-cultured hepatocytes to investigate the hepatobiliary disposition of taurocholic acid (TCA) in the presence of cholestatic agents.

Materials & Methods

Materials: Human hepatocytes were isolated and purified at XenoTech, LLC and characterized for use at Qualyst Transporter Solutions, LLC. Seeding, overlay, and culturing media, and Plus (+) and Minus (-) Assay Buffers were obtained from Qualyst Transporter Solutions. [³H]-taurocholic acid (TCA) was obtained from Perkin Elmer Life and Analytical Sciences (Waltham, MA, USA). Bovine serum albumin, carbamazepine, cisapride monohydrate, cyclosporine A, fluphenazine, glyburide, nicardipine, omeprazole, sulpiride, and troglitazone were obtained from Sigma (St. Louis, MO, USA). Ezetimibe and olmesartan medoxomil were obtained from Cayman Chemicals (Ann Arbor, MI, USA).

Cell Culture: Human hepatocytes were seeded at a density of 1.3x10⁶ cells/mL in 24 well cell culture plates. Approximately 3 hours after seeding, matrigel overlay (0.25 mg/mL) was added to the wells and allowed to polymerize overnight. Cells were fed everyday for 5 days and kept at 37°C, 95% relative humidity, and 5% CO₂.

B-CLEAR® Assay: On day 6, cells were removed from the incubator and rinsed 3 times with Plus (+) Buffer. Cells were then pre-incubated in 300 µL Plus (+) or Minus (-) Buffer for 10 min at 37°C with one of the following inhibitors: carbamazepine, cyclosporine A (CspA), cisapride monohydrate, ezetimibe, fluphenazine hydrochloride, fluvastatin, glyburide, Nicardipine, olmesartan, omeprazole, sulpiride, troglitazone. Pre-incubation buffer was then removed and replaced with 300 µL Plus (+) buffer containing 2.5 µM taurocholic acid (TCA) and the inhibitors previously listed and incubated in the presence of 4% BSA for an additional 30 min at 37°C. At the end of the incubation, cells were washed three times in Plus (+) buffer, lysed with 0.1N NaOH and the lysate was analyzed using liquid scintillation counting. In follow up studies, a dose-response was generated with seven concentrations of selected compounds (nicardipine, ezetimibe, fluvastatin, cyclosporine A, and cisapride monohydrate) under the same conditions described above.

Results

Table 1. Summary of screening results for clinically relevant cholestatic agents

Inhibitor	Predicted Cholestasis Concern from BSEP IC50	Inhibitor Dose (µM)	ICC of TCA (µM)	BEI (%)	Kp	% Change of Control Kp
Solvent Control	NA	NA	17.5	31.4	6.83	100%
Carbamazepine	Unknown	100	17.9	42.9	7.18	105%
Cisapride Monohydrate	Yes	20	24.3	14.6	9.72	142%
Cyclosporine A	Yes	1	9.91	11.8	3.96	42%
Ezetimibe	No	100	9.26	43.3	3.70	46%
Fluphenazine	Yes	100	19.7	46.1	7.89	116%
Fluvastatin	Maybe	100	14.1	43.4	5.65	83%
Glyburide	Yes	1	19.4	31.5	7.78	114%
Nicardipine	Yes	5	22.0	27.6	8.78	129%
Olmesartan medoxomil	No	1	16.4	43.5	6.57	96%
Omeprazole	Yes	30	16.4	45.5	6.55	96%
Sulpiride	Yes	500	19.2	36.9	7.68	112%
Troglitazone	Yes	1	15.8	34.9	6.31	92%

Table 1 lists the results of screening a panel of clinically relevant cholestatic agents for their effect on intracellular concentration (ICC), biliary excretion index (BEI), and Kp (ratio of ICC to extracellular concentration). Results indicate that compared to solvent control (0.1% DMSO), there was a dramatic shift in the ICC of TCA in the hepatocytes for cisapride monohydrate, cyclosporine A, ezetimibe, fluvastatin, and nicardipine. Furthermore, BEI and Kp were also greatly impacted by exposure to these inhibitors, leading to changes in the hepatobiliary disposition. Follow up studies investigated a range of concentrations for the bolded inhibitors below.

Figure 1. Nicardipine total accumulation (pmol/mg) and average BEI (%) (n=3)

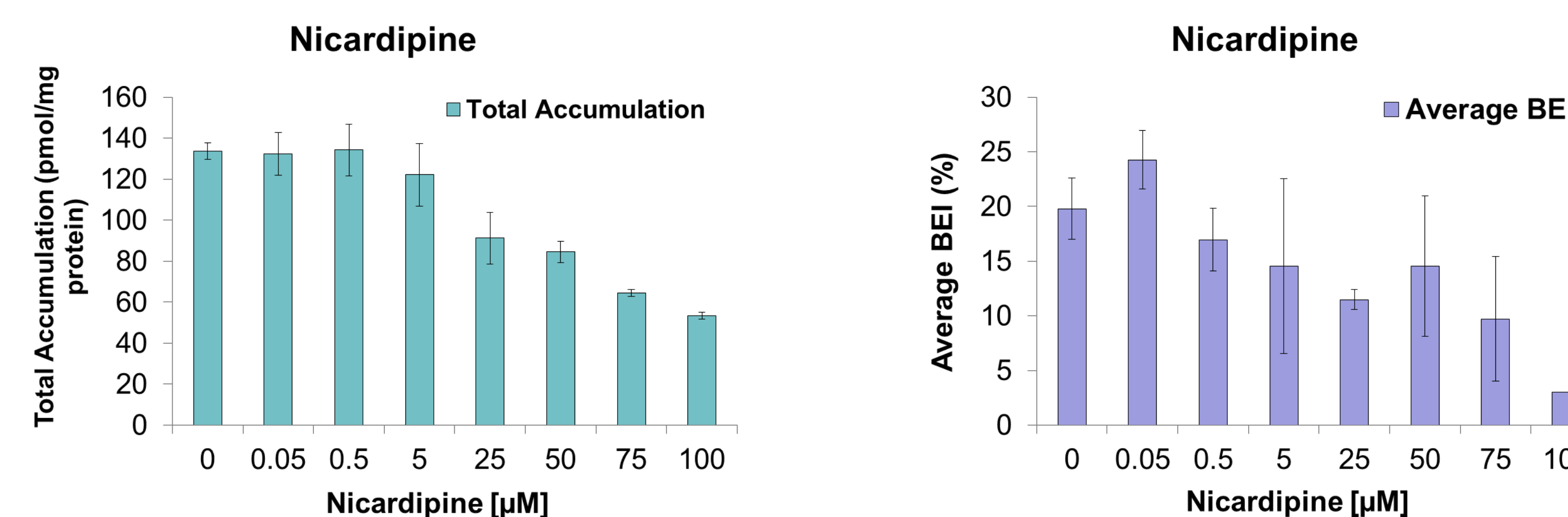


Figure 2. Ezetimibe total accumulation (pmol/mg) and average BEI (%) (n=3)

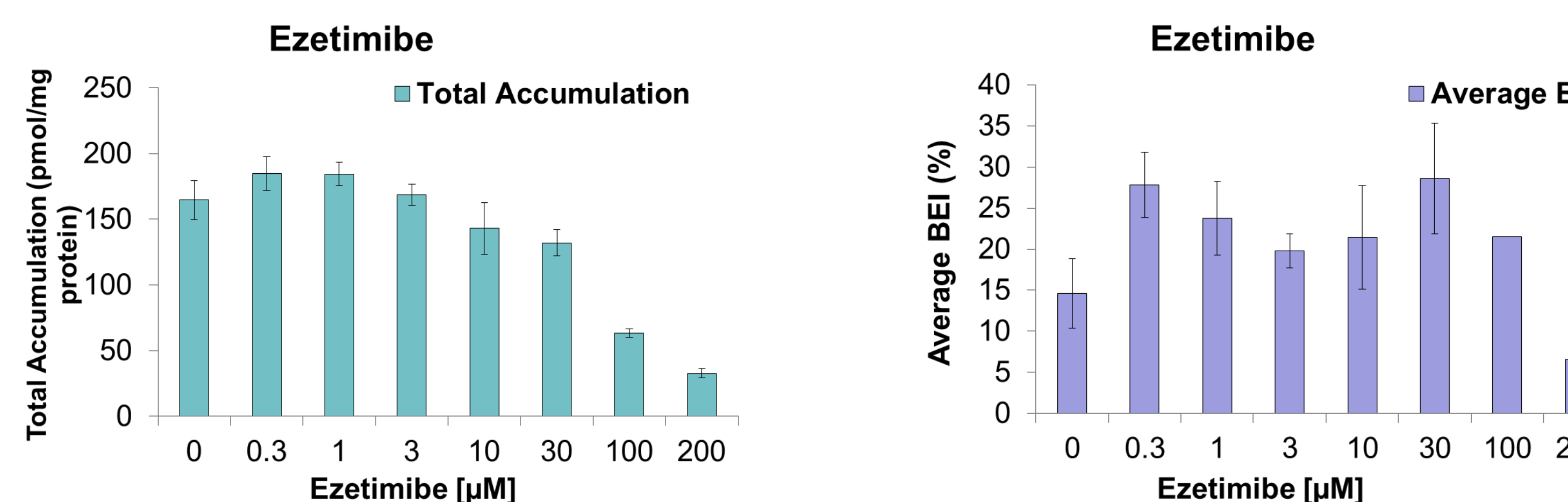


Figure 3. Fluvastatin total accumulation (pmol/mg) and average BEI (%) (n=3)

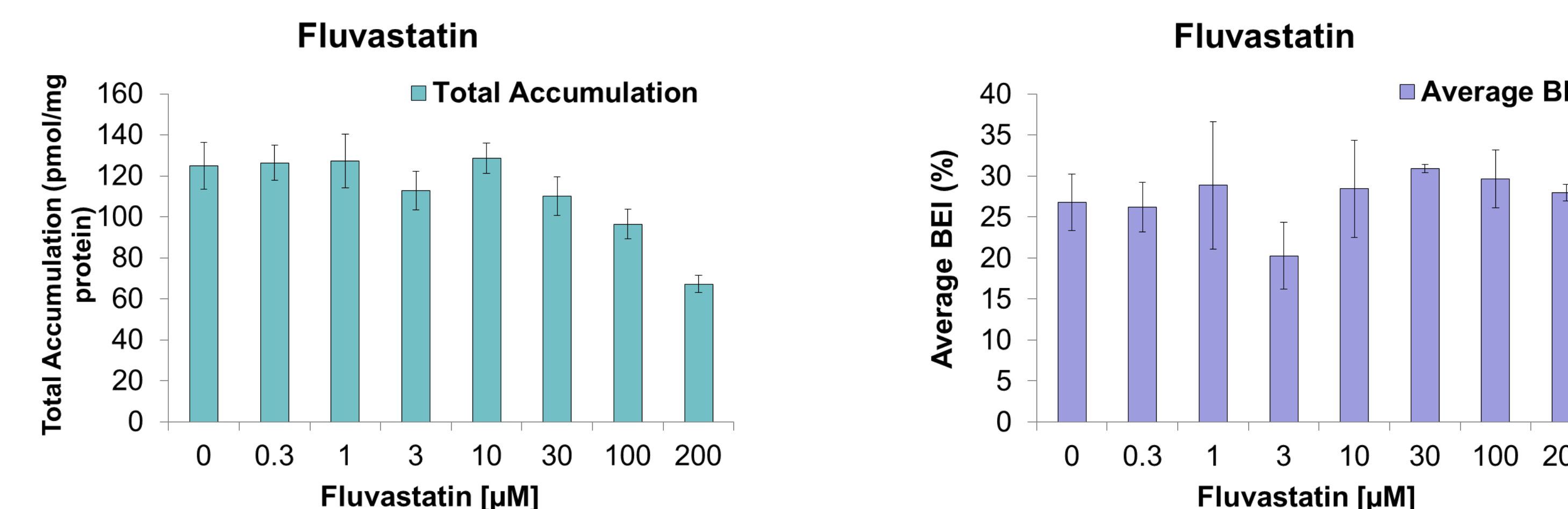


Figure 4. Cyclosporine A total accumulation (pmol/mg) and average BEI (%) (n=2)

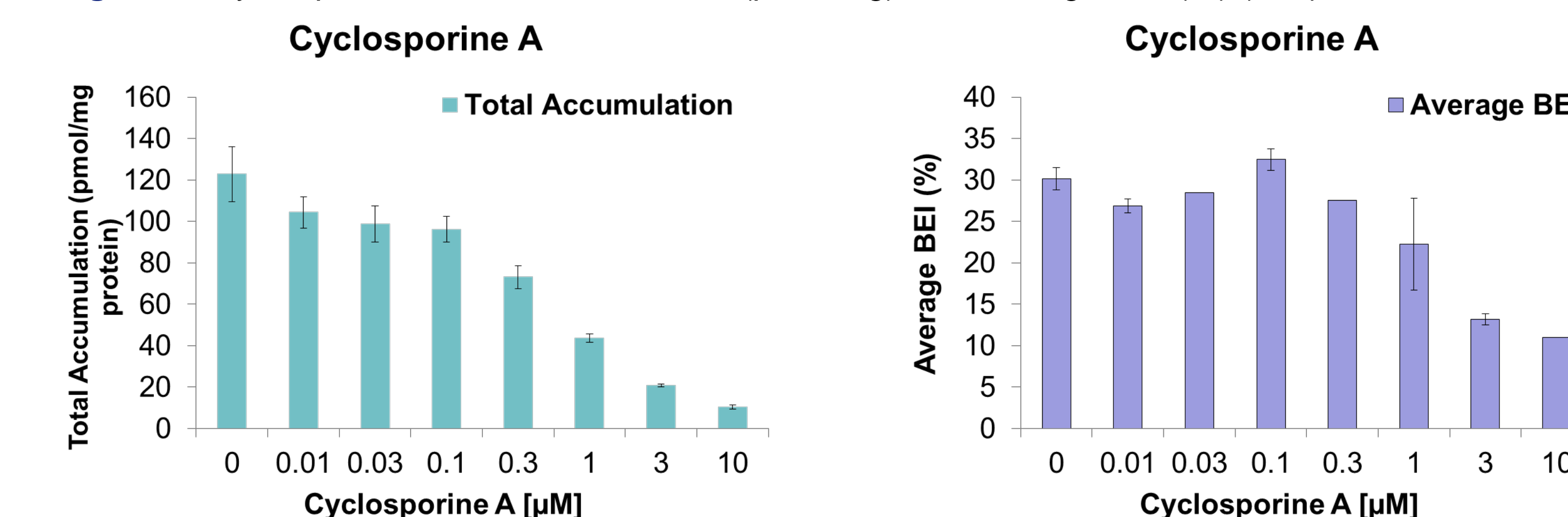
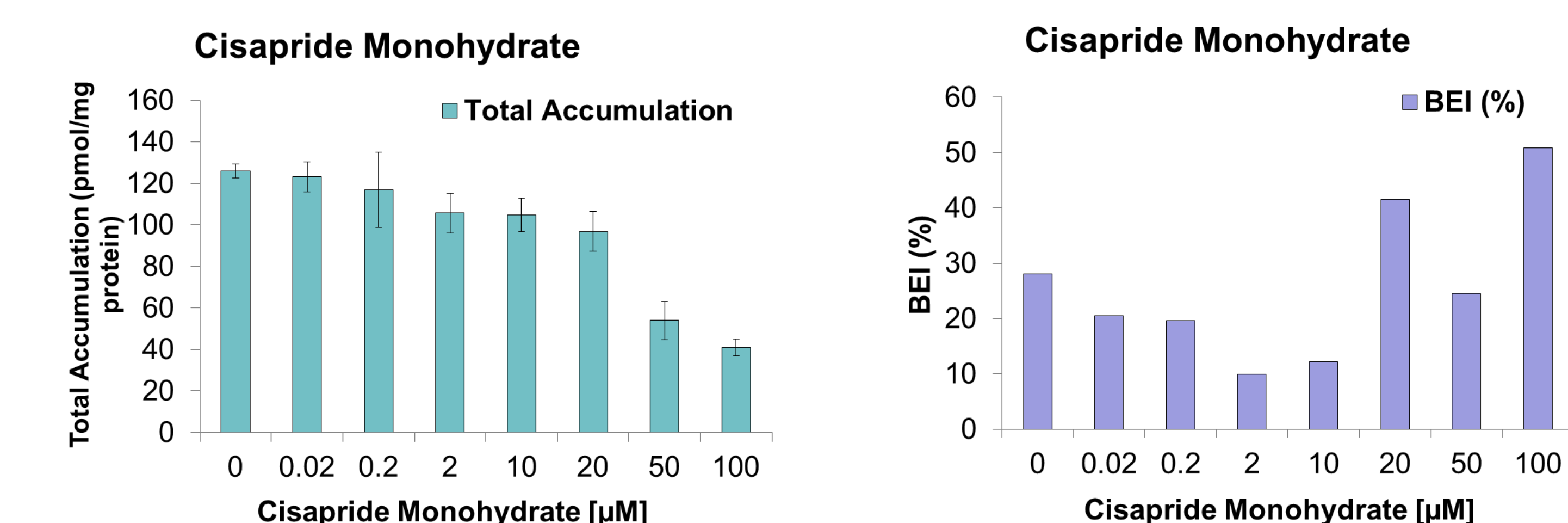


Figure 5. Cisapride monohydrate total accumulation (pmol/mg) and BEI (%) (n=1)



Figures 1-5 show the total accumulation of TCA (2.5 µM) and the average biliary excretion index (BEI, %) following a 30 min incubation with the probe substrate. Nicardipine (Figure 1), demonstrated a concentration-dependent decrease in total accumulation of TCA at concentrations greater than 5 µM and a concentration dependent decrease in the BEI. Ezetimibe (Figure 2) demonstrated a decrease in total accumulation of TCA as the concentration of the inhibitor increased, but displayed an initial increase in BEI, followed by a subsequent decrease in the average BEI. While total accumulation of TCA in the presence of fluvastatin decreased in a concentration-dependent manner (Figure 3), the average BEI was not altered. Figure 4 displays the results for cyclosporine A, which exhibited a concentration-dependent decrease in total accumulation, with an initial increase in BEI, followed by a decrease as the level of compound in the system rose. Finally, cisapride monohydrate (Figure 5), demonstrated a concentration dependent decrease in total accumulation of TCA, and conversely, exhibited an increase in BEI as levels of cisapride increased.

Conclusions

The B-CLEAR® human hepatocyte sandwich cultured assay demonstrates the ability to utilize the accumulation of a probe substrate to assess both inhibition of uptake as well as efflux. Many xenobiotics which inhibit efflux out of hepatocytes can lead to a risk of hepatotoxicity. It is important to note though, that a compound capable of inhibiting uptake into hepatocytes may lead to a risk of systemic toxicity. Therefore, the use of an *in vitro* system which more accurately models *in vivo* conditions can help to better predict the overall effect of xenobiotics on uptake, clearance and potential toxicity. These data presented here demonstrate that hepatobiliary uptake and clearance are complicated processes which are influenced by a number of factors. The Qualyst Transporter Solutions B-CLEAR® technology paired with XenoTech-produced sandwich-cultured hepatocytes are able to better assess inhibition of both uptake and efflux in hepatocytes. Future directions include further investigation of toxicity utilizing human sandwich-cultured hepatocytes in order to better understand the effect of acute and chronic exposure to these clinically relevant cholestatic agents.

References

- Suk, Ki Tae, and Dong Joon Kim. "Drug-Induced Liver Injury: Present and Future." *Clinical and molecular hepatology* 18.3 (2012): 249–257. *PMC*. Web. 24 Mar. 2015.
- Morgan, R.E., Trauner, M., van Staden, C.J., Lee, P.H., Ramachandran, B., Eschenberg, M., Afshari, C.A., Qualls, C.W., Lightfoot-Dunn, R., Hamadeh, H.K., "Interference with Bile Salt Export Pump Function is a Susceptibility Factor for Human Liver Injury in Drug Development." *Toxicol. Sci.*, (2010) 118, (2), 485-500.
- Chen, M., Bisgin, H., Tong, L., Fang, H., Borlak, J., Tong, W., "Toward predictive models for drug-induced liver injury in humans: are we there yet?" (2014) *Biomarkers Med.*, 8 (2), 201-213.
- Yuan L, Kaplowitz N. "Mechanisms of Drug Induced Liver Injury." *Clinics in liver disease*. 2013;17(4):507-518.
- Morgan, R.E., van Staden, C.J., Chen, Y., Kalyanaraman, N., Kalanzi, J., Dunn II, R.T., Afshari, C.A., Hamadeh, H.K., "A Multifactorial Approach to Hepatociliary Transporter Assessment Enables Improved Therapeutic Compound Development." *Toxicol. Sci.* (2013), 136 (1), 216-241.