

Introduction

Organic anion transporting polypeptides 1B1 and 1B3 (OATP1B1, *SLCO1B1* and OATP1B3, *SLCO1B3*) are transmembrane proteins expressed in the sinusoidal membrane of human hepatocytes and are capable of transporting a wide range of hydrophilic molecules from the blood into the hepatocyte for subsequent metabolism and/or elimination. Inhibition of OATP1B1 and/or OATP1B3 by xenobiotics can lead to pharmacokinetic drug-drug interactions whereby the exposure (AUC) of a co-administered victim drug (substrate) is increased in the presence of the inhibitory drug. For example, orally administered cyclosporin inhibits OATP-mediated uptake of pravastatin, a prototypical *in vivo* probe substrate of OATP, which causes a 9.9-fold increase in AUC.¹ The increase in systemic exposure to narrow safety margin drugs (e.g. statins) results in undesired side-effects and dose-limiting toxicities. Therefore, the FDA and EMEA recommend the evaluation of new drug candidates for the potential to inhibit OATP1B1 and OATP1B3 which is often first evaluated with *in vitro* test systems, typically IC_{50} experiments.

The recent EMEA Guideline on the Investigation of Drug Interactions (2012) recommends the calculation of K_i values when evaluating the inhibition of a transporter by investigational drugs.² Although K_i values have been reported for inhibitors of various uptake transporters, the difficulty of experimental design coupled with vigorous mathematical extrapolation can make the determination of intrinsic K_i values difficult and oftentimes reverts to the determination of extrinsic IC_{50} values. Additionally, it is important to consider clinically relevant substrates rather than prototypical probe substrates when evaluating potential DDIs of investigational drugs at the transporter level. In this study, K_i and IC_{50} values were determined for one prototypical (estradiol-17 β -glucuronide (E2G)) and one clinically-relevant (pravastatin) substrate of the hepatic uptake transporters OATP1B1 and OATP1B3 with an experimental design incorporating multiple substrate and inhibitor concentrations.

Methods

HEK293 cells over-expressing either OATP1B1 or OATP1B3 were obtained from Sekisui Medical Division, LTD (Tokai, Japan). Pravastatin and rifampin were obtained from Sigma (St. Louis, MO, USA). [³H]Estradiol-17 β -glucuronide (E2G) was obtained from Perkin Elmer Life and Analytical Sciences (Waltham, MA, USA).

In brief, HEK293 cells over-expressing OATP1B1 or OATP1B3 were cultured in 24-well plates. Cells were pre-incubated (15 min at 37°C) with four concentrations of the inhibitor, rifampin, at concentrations bracketing the experimentally determined IC_{50} value (typically 1/6 to 6-fold the experimentally determined IC_{50}). The pre-incubation was followed by incubation (1 min) with six concentrations of substrate (typically 1/3 to 3-fold the experimentally determined K_m) and inhibitor. Uptake was terminated by washing the cells three times with cold 1x PBS. For analysis by scintillation counting (E2G), cells were lysed with 0.1 M NaOH. For analysis by LC-MS/MS (pravastatin), cells were extracted with 50:50 v/v methanol:water. IC_{50} values were determined with GraphPad Prism software (version 6.01). K_m and K_i values were determined with GraFit software (version 7.0.2).

Results

Figure 1. Experimentally determined inhibition profiles (IC_{50} and K_i) of OATP1B1-mediated uptake of E2G and pravastatin into HEK-OATP1B1 cells by rifampin

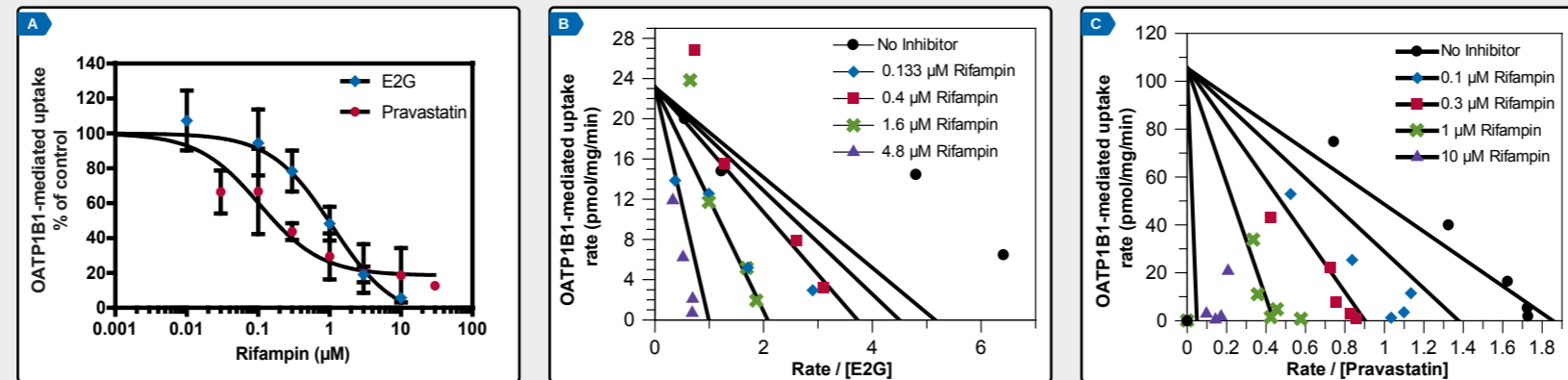


Figure 2. Experimentally determined inhibition profiles (IC_{50} and K_i) of OATP1B3-mediated uptake of E2G and pravastatin into HEK-OATP1B3 cells by rifampin

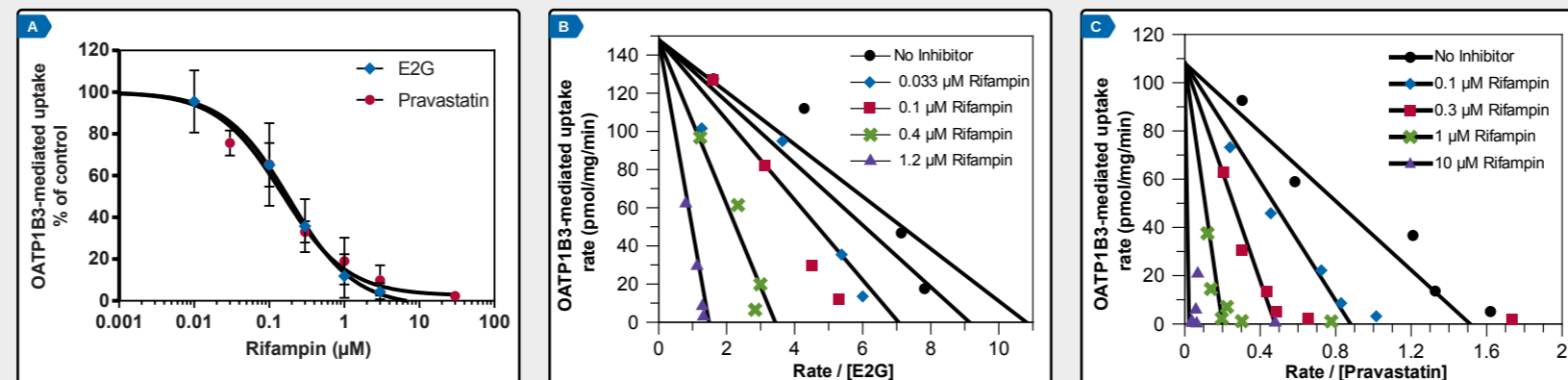


Table 1. Summary of experimentally-determined IC_{50} and K_i values for the inhibition of OATP1B1- and OATP1B3-mediated uptake by rifampin

Transporter	Substrate	Inhibitor	K_m (μ M)	IC_{50} (μ M)	K_i (μ M)	Mechanism (best fit)
OATP1B1	E2G	Rifampin	4.5 \pm 1.9	0.79 \pm 0.13	1.1 \pm 0.6	Competitive
OATP1B3	E2G	Rifampin	13.7 \pm 7.7	0.19 \pm 0.09	0.19 \pm 0.04	Competitive
OATP1B1	Pravastatin	Rifampin	56.8 \pm 15.7	0.16 \pm 0.02	0.28 \pm 0.08	Competitive
OATP1B3	Pravastatin	Rifampin	71.6 \pm 10.7	0.25 \pm 0.05	0.14 \pm 0.02	Competitive

- Figure 1** illustrates the inhibition of OATP1B1-mediated uptake of E2G and pravastatin by rifampin. Rifampin inhibited E2G and pravastatin uptake into OATP1B1-expressing cells with IC_{50} values of 0.79 and 0.16 μ M, respectively (**Figure 1A and Table 1**). Rifampin inhibited the uptake of E2G and pravastatin with K_i values of 1.1 and 0.28 μ M, respectively (**Figure 1B and 1C**).
- Experiments conducted with E2G as a probe substrate for OATP1B1 yielded comparable K_i and IC_{50} values (1.1 and 0.79 μ M, respectively). Similar results were observed with the clinically-relevant probe substrate pravastatin (K_i and IC_{50} values of 0.28 and 0.16 μ M, respectively).
- Figure 2** illustrates the inhibition of OATP1B3-mediated uptake of E2G and pravastatin by rifampin. Rifampin inhibited E2G and pravastatin uptake into OATP1B3-expressing cells with IC_{50} values of 0.19 and 0.25 μ M, respectively (**Figure 2A and Table 1**). Rifampin inhibited the uptake of E2G and pravastatin with K_i values of 0.19 and 0.14 μ M, respectively (**Figure 2B and 2C**).
- Similar to the observed results for OATP1B1, experiments conducted with E2G as a probe substrate for OATP1B3 yielded comparable K_i and IC_{50} values (0.19 and 0.19 μ M, respectively). Also, similar results were observed with the clinically-relevant probe substrate pravastatin (K_i and IC_{50} values of 0.14 and 0.25 μ M, respectively).
- However, substrate-dependent inhibition was observed in OATP1B1 cells but not in OATP1B3 cells; rifampin inhibited OATP1B1-mediated pravastatin uptake with approximately four-fold more potency than observed with E2G with K_i values of 1.1 and 0.28 μ M, respectively (**Table 1**).
- In all cases, the mechanism of inhibition for OATP1B1- and OATP1B3-mediated uptake of E2G and pravastatin by rifampin was determined to be competitive according to the best fit of the experimental data (**Table 1**).

Conclusions

- Substrate concentrations used in IC_{50} experiments were much lower than K_m and, therefore, the IC_{50} values were expected to be similar to the K_i value.
- Overall, K_i values were similar to IC_{50} values (within two-fold) for inhibition of OATP1B1 and OATP1B3.
- Substrate-dependent inhibition by rifampin was observed in OATP1B1 cells but not in OATP1B3 cells; rifampin inhibited OATP1B1-mediated pravastatin uptake with approximately four-fold more potency than observed with E2G.
- In some cases, experimentally determined K_i values may be more accurate for DDI predictions than IC_{50} values, and, in some cases, should be determined with a clinically relevant substrate.

References

- Zhang L et al. (2011) Clin Pharmacol Ther. **89**: 481-484.
- [EMA] European Medicines Agency. EMA/CHMP/EWP/125211/2012. Guideline on the Investigation of Drug Interactions, Committee for Human Medicinal Products (CHMP), London, England.