

Introduction

The identification of the metabolic pathways (reaction phenotyping) for new drug candidates is an essential component of drug development, as these studies increase the understanding of how a drug is cleared and the potential for the drug to be subject to drug-drug interactions with co-administered medications. Typically *in vitro* reaction phenotyping is conducted in early drug development by measuring disappearance of parent drug, as metabolite standards are typically unavailable. In the industry, these assays have been conducted with various approaches, including the use of recombinant CYP enzymes, antibody inhibition of CYP enzymes and correlation analysis.

Another approach used to identify which cytochrome P450 (CYP) enzymes are responsible for the metabolism of a drug is the use of specific chemical inhibitors in test systems such as pooled human liver microsomes (HLM) or pooled cryopreserved human hepatocytes (CHH) (Ogilvie *et al.*, 2008; Khojasteh *et al.*, 2011; Nirogi *et al.*, 2014). These studies are typically more cost effective and offer logistical advantages over the other approaches. However, the design of these studies is critical because of factors such as metabolic depletion of the inhibitor, protein binding and insufficient enzyme inactivation. Additionally, the measurement of enzyme inactivation under initial rate substrate conditions is crucial to obtain an accurate assessment of the true levels of inhibition (Parkinson *et al.*, 2011).

In the present study, we examined the selectivity and duration of CYP inactivation, with a range of protein concentrations and incubation times, for a variety of commonly used CYP inhibitors all under initial rate probe substrate conditions in both HLM and CHH with the goal of optimizing assay conditions.

Materials & Methods

Chemicals

Furafylline, ketoconazole, mibefradil, paroxetine, phencyclidine and quinidine were purchased from Sigma-Aldrich (St. Louis, MO). Esomeprazole, gemfibrozil glucuronide and CYP3cide were purchased from Toronto Research Chemicals (Toronto, ON, Canada). Tienilic acid was purchased from Cypex (Dundee, Scotland, UK). Troleandomycin was purchased from US Pharmacopeia (Rockville, MD). The sources of all other reagents have been described previously (Parkinson *et al.*, 2011).

Test system

Pooled human liver microsomes (HLM, n = 200, mixed gender) and pooled cryopreserved human hepatocytes (CHH, n = 100, mixed gender) were prepared from non-transplantable livers and characterized at XenoTech, LLC (Lenexa, KS) as described previously (Pearce *et al.*, 1996; Parkinson *et al.*, 2004).

In vitro chemical inhibition under initial rate conditions in HLM and CHH

Briefly as described in Figure 1, NADPH-fortified pooled HLM (n = 200) at 0.1, 0.5 and 1 mg/mL or pooled CHH (n = 100) at 1 million cells/mL plated in 48-well plates, were pre-incubated for 30 min at 37°C with various CYP inhibitors (as shown in Table 1), namely furafylline (10 µM), phencyclidine (up to 30 µM), gemfibrozil glucuronide (100 µM), tienilic acid (20 µM), esomeprazole (10 µM), quinidine (5 µM), paroxetine (1 or 5 µM), ketoconazole (1 or 4 µM), CYP3cide (2.5 µM), mibefradil (1 µM) and troleandomycin (50 µM). Following the pre-incubation step, marker substrate ($\approx K_m$) incubations were performed for up to 5 min (10 min for CHH) at three time points (0, 30, 120 min; simulating the time course of an unknown drug) following the pre-incubation step to determine CYP1A2 (phenacetin), CYP2B6 (bupropion), CYP2C8 (amodiaquine), CYP2C9 (diclofenac), CYP2C19 (S-mephenytoin), CYP2D6 (dextromethorphan), CYP2E1 (chlorzoxazone) and CYP3A4/5 (midazolam) residual activities under initial rate conditions. Reactions were terminated with the addition of one volume of acetonitrile containing deuterated internal standard, followed by protein precipitation (10 min, 920 RCF). Metabolite formation was determined by LC-MS/MS analysis as described previously (Parkinson *et al.*, 2011).

Figure 1. Incubation process for the assessment of chemical inhibition at initial rate conditions

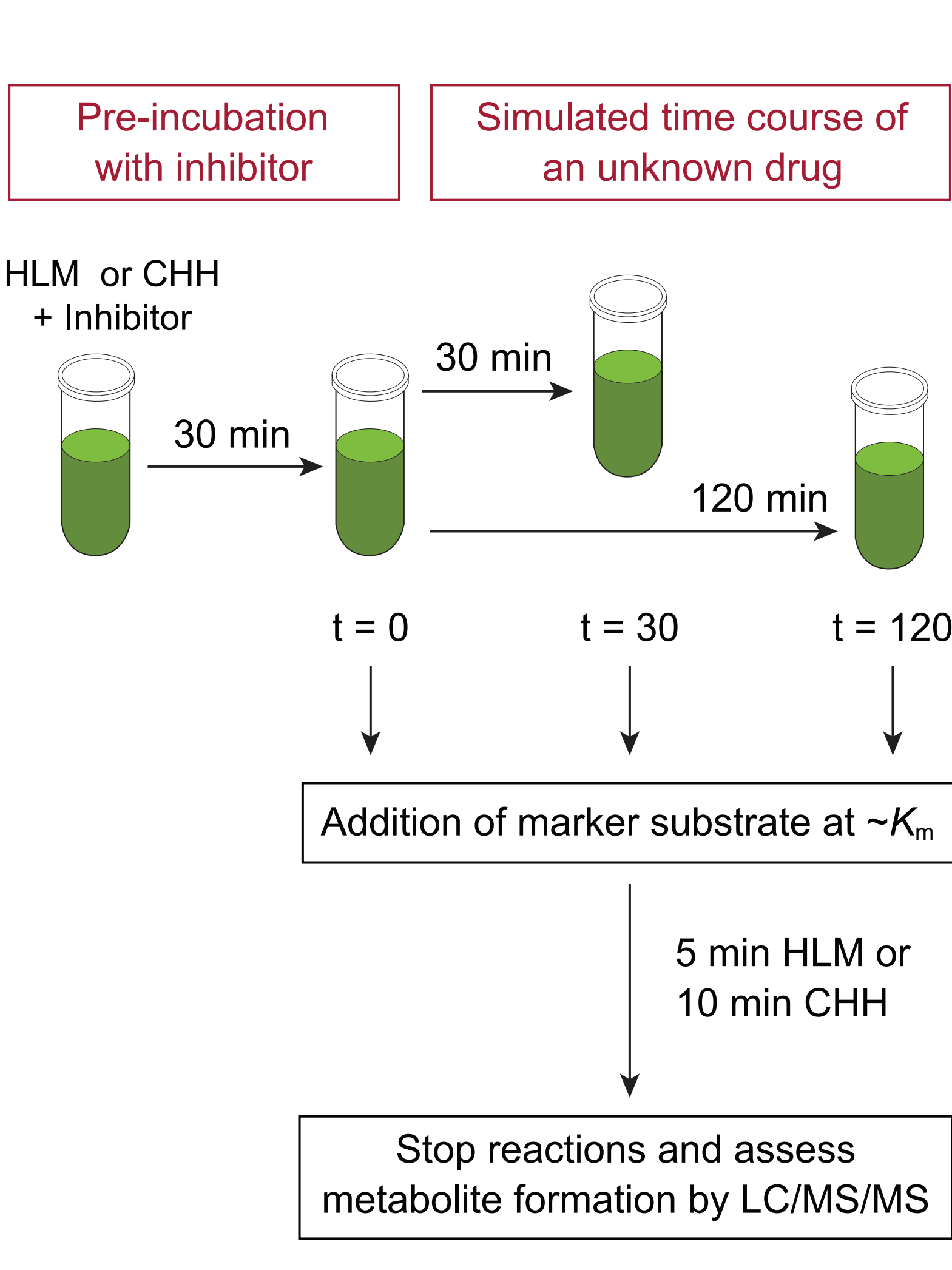


Table 1. Summary of inhibitor concentrations in each test system

Inhibitor	HLM	CHH
Furafylline		10 µM
Phencyclidine	30 µM	10 µM
Gemfibrozil gluc		100 µM
Tienilic acid		20 µM
Esomeprazole	10 µM	
Paroxetine	5 µM	1 µM
Quinidine		5 µM
Mibefradil		1 µM
CYP3cide		2.5 µM
Troleandomycin		50 µM
Ketoconazole	1 µM	4 µM

Results

Results indicated that, in both HLM and CHH, inhibition of specific CYP enzymes can be accomplished over the duration of 120 minutes after pre-incubation with the inhibitors used. As summarized in Figure 2, in HLM at 0.1, 0.5 and 1 mg/mL over the time course of the incubation, furafylline selectively inactivated CYP1A2 activity by 63-80%; phencyclidine selectively inactivated CYP2B6 activity by 48-91%; gemfibrozil glucuronide selectively inactivated CYP2C8 activity by 96-98%; tienilic acid selectively inactivated CYP2C9 activity by 96-100%; paroxetine inactivated CYP2D6 activity by 55-90% with moderate inhibition of CYP2B6; quinidine selectively inactivated CYP2D6 by 55-90%; and ketoconazole, CYP3cide, mibefradil and troleandomycin inactivated CYP3A4/5 activity by 78-97%, 81-88%, 72-91% and 92-94% respectively. Esomeprazole was found to potentially inhibit CYP2C19 activity by 85-100% with some moderate inhibition of CYP2E1 (53-66%) and CYP3A4/5 (31-68%).

Figure 3 summarizes the results from CHH (1 million cells/mL) over the incubation time course. Briefly, furafylline selectively inactivated CYP1A2 activity by 83-87%; phencyclidine inactivated CYP2B6 activity by 63-90% with moderate inhibition of CYP2D6 (37-56%); gemfibrozil glucuronide selectively inactivated CYP2C8 activity by 32-97%; tienilic acid selectively inactivated CYP2C9 activity by 100%; paroxetine inactivated CYP2D6 activity by 72-85% with moderate inhibition of CYP2B6; quinidine selectively inactivated CYP2D6 activity by 87-94%; and ketoconazole, CYP3cide, mibefradil, and troleandomycin inactivated CYP3A4/5 activity by 96-97%, 83-91%, 63-89% and 92-95% respectively. Esomeprazole was found to potentially inhibit CYP2C19 activity by 89-95% with some moderate inhibition of CYP2E1 (68-69%), consistent with results from HLM.

Figure 2. Time course of inactivation of CYP enzymes in HLM at 0.1, 0.5 and 1 mg/mL by various chemical inhibitors under initial rate conditions

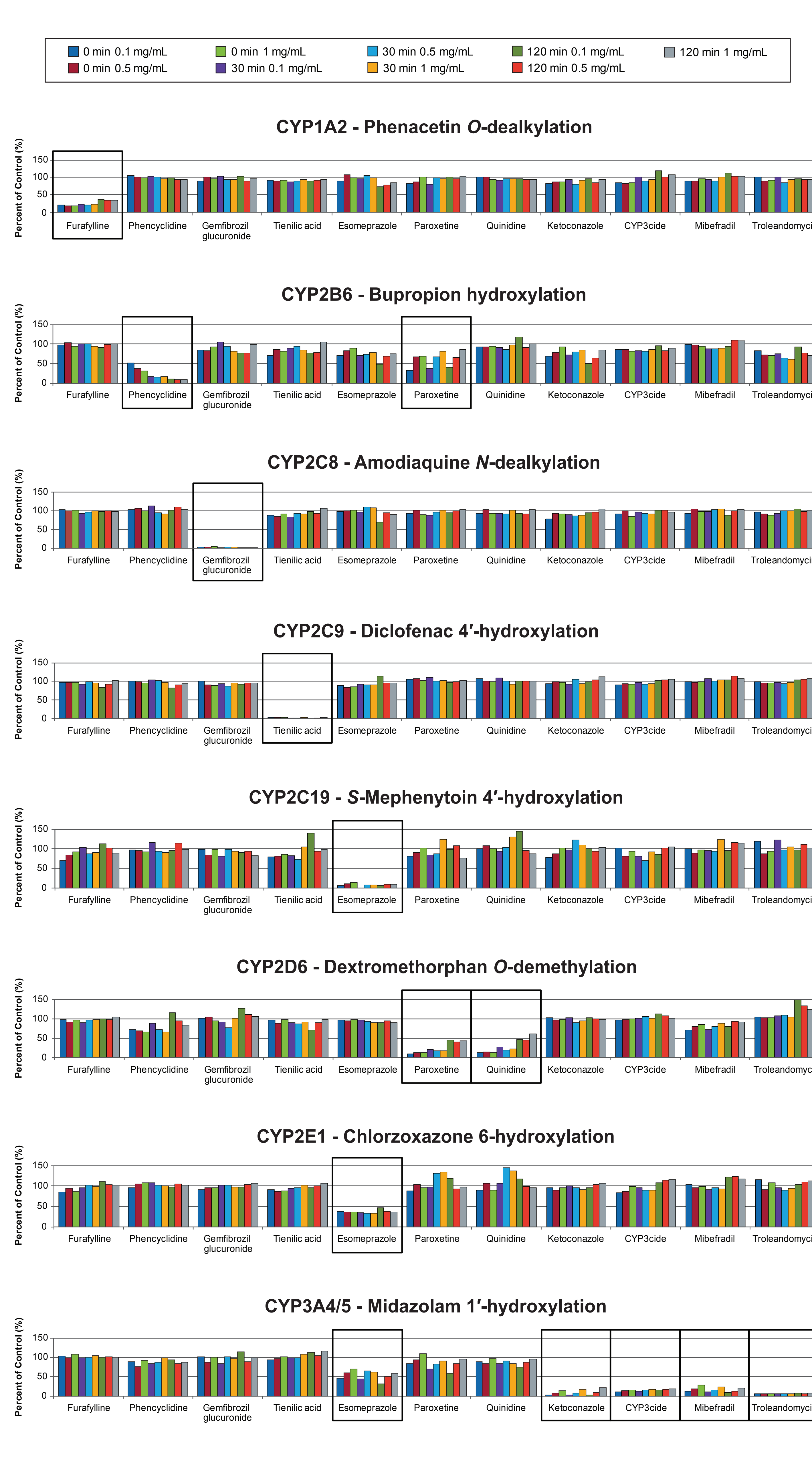
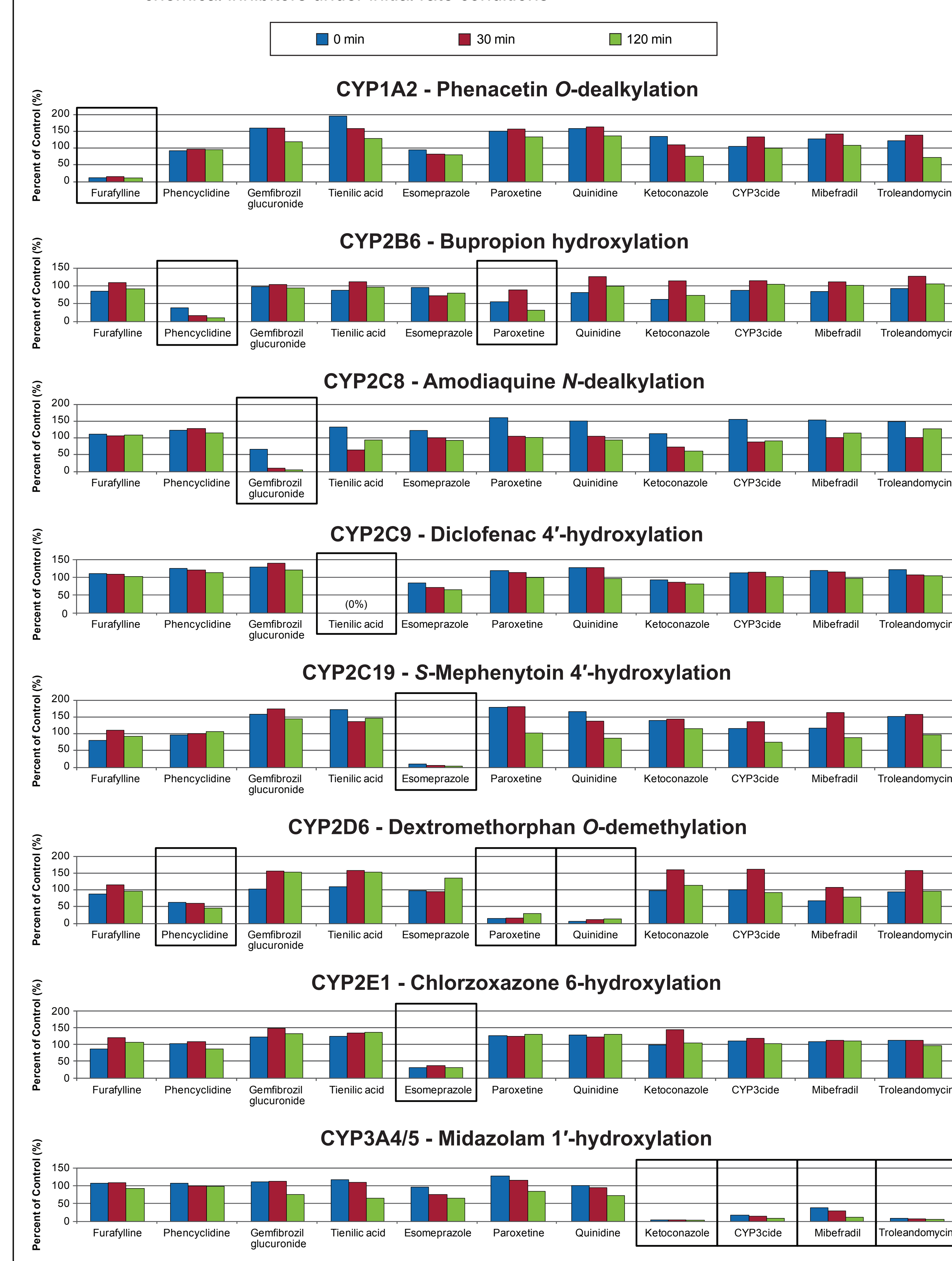


Figure 3. Time course of inactivation of CYP enzymes in CHH at 1 million cells/mL by various chemical inhibitors under initial rate conditions



Conclusions

- 11 chemical inhibitors were found to be selective for various CYP enzymes over 120 min after pre-incubation in both HLM at multiple protein concentrations and CHH.
- In summary, these findings demonstrate the suitability and optimized conditions for the use of common chemical inhibitors in CYP reaction phenotyping studies, with implications for the design of such studies.

References

- Khojasteh SC, Prabhu S, Kenny JR, Halladay JS, and Lu AY (2011) Chemical inhibitors of cytochrome P450 isoforms in human liver microsomes: a re-evaluation of P450 isoform selectivity. *Eur J Drug Metab Pharmacokin* 36:1-16.
- Nirogi R, Palacharla RC, Uthukam V, Manoharan A, Srikalopu SR, Kalaikadhiban I, Boggavarapu RK, Ponnamaneni RK, Ajjala DR, and Bhyrapuneni G (2014) Chemical inhibitors of CYP450 enzymes in liver microsomes: combining selectivity and unbound fractions to guide selection of appropriate concentration in phenotyping assays. *Xenobiotica* doi: 10.3109/00498254.2014.945196.
- Ogilvie BW, Usuki E, Yerino P, and Parkinson A (2008) In vitro approaches for studying the inhibition of drug-metabolizing enzymes and identifying the drug-metabolizing enzymes responsible for the metabolism of drugs (Reaction Phenotyping) with emphasis on cytochrome P450, in: *Drug-Drug Interactions* (Rodrigues AD ed), pp 231-358, Informa Healthcare USA Inc., New York, NY.
- Parkinson A, Kazmi F, Buckley DB, Yerino P, Paris BL, Holsapple J, Toren P, Otradovec SM, and Ogilvie BW (2011) An Evaluation of the Dilution Method for Identifying Metabolism-dependent Inhibitors (MDIs) of Cytochrome P450 (CYP) Enzymes. *Drug Metab Dispos* 39:1370-1387.
- Parkinson A, Mudra DR, Johnson C, Dwyer A, and Carroll KM (2004) The effects of gender, age, ethnicity, and liver cirrhosis on cytochrome P450 enzyme activity in human liver microsomes and inducibility in cultured human hepatocytes. *Toxicol Appl Pharmacol* 199:193-209.
- Pearce RE, Rodrigues AD, Goldstein JA, and Parkinson A (1996) Identification of the human P450 enzymes involved in lansoprazole metabolism. *J Pharmacol Exp Ther* 277:805-816.