

148 Inhibitory effect of the CYP typical inhibitors under long term incubation in human liver microsomes



Correspondence : Sho Nishinoaki
sho.nishinoaki@sekisui.com

Sho Nishinoaki, Ryo Fujino, Kenta Hashizume, Tsutomu Negama

Drug Development Solutions Center, Drug Development Solutions Division, Sekisui Medical Co., Ltd.

2117 Muramatsu, Tokai, Ibaraki 319-1182, Japan

Introduction

Cytochrom P450 (CYP) isoforms are responsible for the metabolism of the majority of drugs in human. In case of the metabolic enzyme identification of CYP isoforms for the low intrinsic clearance (CL_{int}) candidate compounds, it is necessary to perform metabolic reaction with long term incubation. In addition, when performing the metabolic enzyme identification with long term incubation, the specificity and sustainability of a typical inhibitor are important. In this research, we evaluated the inhibitory effect of typical inhibitor for each CYP isoform in human liver microsomes (HLM) under long term incubation (incubation time of CYP typical inhibitors was 60 minutes, final concentration of HLM was 1 mg protein/mL) to clarify an optimal concentration of the CYP typical inhibitors under long term incubation. In this presentation, we have confirmed specificity and sustainability of each typical inhibitor for CYP isoform under long term incubation in HLM.

Method

<Reaction condition and typical inhibitor for each CYP isoform>

CYP isoform	Metabolic activity	* Model substrate concentration (μM)	Typical inhibitor
CYP1A2	Phenacetin O-deethylation	30	α -Naphthoflavone
			Furafylline
CYP2B6	Bupropion hydroxylation	100	Thio-TEPA
CYP2C8	Amodiaquine N-deethylation	2.5	Montelukast
CYP2C9	Diclofenac 4'-hydroxylation	4	Sulfaphenazole
CYP2C19	S-(+)- Mephenytoin 4'-hydroxylation	30	Benzylrinivorol
CYP2D6	(\pm)- Bufuralol 1'-hydroxylation	10	Quinidine
CYP3A4	Testosterone 6 β -hydroxylation	60	Ketoconazole
CYP3A4	Midazolam 1'-hydroxylation	1	Ketoconazole

*The model substrate concentrations were set to around K_m values based on internal data of Sekisui Medical Co., Ltd.
Incubation samples contained phosphate buffer (100 mM, pH 7.4), HLM (final concentration: 1 mg protein/mL, H0610, mixed gender, pool of 50, Sekisui Xeno Tech, LLC), NADPH generating system (Corning) and each typical inhibitor were incubated at 37 $^{\circ}$ C for 60 min. The reaction was started by addition of the each model substrate solution, and the incubation samples were incubated at 37 $^{\circ}$ C for 1 min. After incubation, the incubation samples were terminated by adding acetonitrile solution containing internal standard. The metabolite concentrations were determined by LC-MS/MS (HPLC system : LC-10ADvp, Shimadzu, MS/MS system : API4000 and 4000QTRAP, SCIEX).

<Calculation of % of control>
Remaining activity (% of control) of each CYP isoform was calculated by the following equation.

$$\% \text{ of control} = \frac{\text{Peak area ratio in typical inhibitor group}}{\text{Peak area ratio in control group}} \times 100$$

Table 3 and Figure 3 Inhibitory effect of Thio-TEPA for each CYP isoform in HLM

Thio-TEPA	% of control			
	CYP1A2	CYP2B6	CYP2C8	CYP2C9
0 $\mu\text{mol/L}$	100.0	100.0	100.0	100.0
10 $\mu\text{mol/L}$	100.2	7.2	105.4	103.2
30 $\mu\text{mol/L}$	102.7	5.4	95.1	101.0
100 $\mu\text{mol/L}$	97.0	3.5	77.3	91.6

Thio-TEPA	% of control			
	CYP2C19	CYP2D6	CYP3A4 (T)	CYP3A4 (M)
0 $\mu\text{mol/L}$	100.0	100.0	100.0	100.0
10 $\mu\text{mol/L}$	101.1	95.1	81.6	81.3
30 $\mu\text{mol/L}$	99.2	97.8	58.9	66.6
100 $\mu\text{mol/L}$	85.3	96.5	42.2	42.4

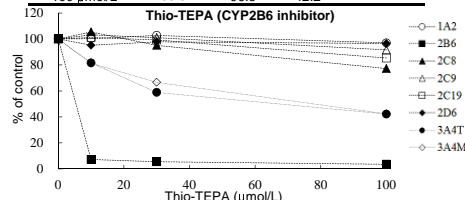


Table 6 and Figure 6 Inhibitory effect of Benzylrinivorol for each CYP isoform in HLM

Benzylrinivorol	% of control			
	CYP1A2	CYP2B6	CYP2C8	CYP2C9
0 $\mu\text{mol/L}$	100.0	100.0	100.0	100.0
1 $\mu\text{mol/L}$	92.3	82.7	101.3	91.5
5 $\mu\text{mol/L}$	86.2	74.1	84.5	93.5
25 $\mu\text{mol/L}$	93.5	65.9	67.2	100.4

Benzylrinivorol	% of control			
	CYP2C19	CYP2D6	CYP3A4 (T)	CYP3A4 (M)
0 $\mu\text{mol/L}$	100.0	100.0	100.0	100.0
1 $\mu\text{mol/L}$	78.6	94.1	93.1	94.5
5 $\mu\text{mol/L}$	13.8	93.0	77.6	77.6
25 $\mu\text{mol/L}$	2.0	114.9	49.2	37.1

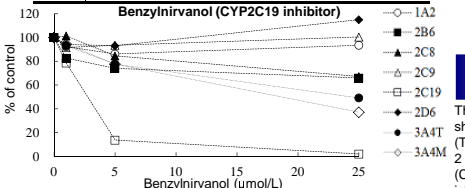


Table 9 IC_{50} of typical inhibitor for each CYP isoform in HLM

Typical inhibitor	IC_{50} ($\mu\text{mol/L}$)	
	15 min	60 min
α -Naphthoflavone (CYP1A2)	0.191	0.273
Furafylline (CYP1A2)	0.605	0.386
Thio-TEPA (CYP2B6)	1.94	1.41
Montelukast (CYP2C8)	1.16	1.69
Sulfaphenazole (CYP2C9)	0.530	0.652
Benzylrinivorol (CYP2C19)	0.441	1.88
Quinidine (CYP2D6)	0.557	1.78
Ketoconazole (CYP3A4 (T))	0.156	0.401
Ketoconazole (CYP3A4 (M))	0.185	0.467

IC_{50} values were calculated using the following equation
% of control = $(IC_{50} / (IC_{50} + \text{typical inhibitor } (\mu\text{mol/L}))) \times 100$

Results

Table 1 and Figure 1 Inhibitory effect of α -Naphthoflavone for each CYP isoform in HLM

α -Naphthoflavone	% of control			
	CYP1A2	CYP2B6	CYP2C8	CYP2C9
0 $\mu\text{mol/L}$	100.0	100.0	100.0	100.0
1 $\mu\text{mol/L}$	17.3	96.2	101.3	92.0
5 $\mu\text{mol/L}$	14.0	92.2	86.7	86.2
25 $\mu\text{mol/L}$	14.3	86.7	44.9	32.9

α -Naphthoflavone	% of control			
	CYP2C19	CYP2D6	CYP3A4 (T)	CYP3A4 (M)
0 $\mu\text{mol/L}$	100.0	100.0	100.0	100.0
1 $\mu\text{mol/L}$	89.0	100.8	92.3	98.7
5 $\mu\text{mol/L}$	68.5	106.6	79.0	119.1
25 $\mu\text{mol/L}$	27.4	98.0	16.8	36.3

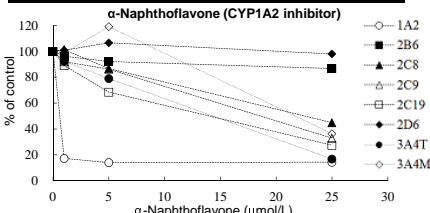


Table 4 and Figure 4 Inhibitory effect of Montelukast for each CYP isoform in HLM

Montelukast	% of control			
	CYP1A2	CYP2B6	CYP2C8	CYP2C9
0 $\mu\text{mol/L}$	100.0	100.0	100.0	100.0
0.2 $\mu\text{mol/L}$	113.1	103.3	95.5	114.3
1 $\mu\text{mol/L}$	107.6	99.7	66.9	101.4
5 $\mu\text{mol/L}$	103.9	97.2	17.1	89.4

Montelukast	% of control			
	CYP2C19	CYP2D6	CYP3A4 (T)	CYP3A4 (M)
0 $\mu\text{mol/L}$	100.0	100.0	100.0	100.0
0.2 $\mu\text{mol/L}$	99.2	95.1	97.3	110.2
1 $\mu\text{mol/L}$	99.4	97.8	93.2	111.2
5 $\mu\text{mol/L}$	89.5	96.5	92.3	103.9

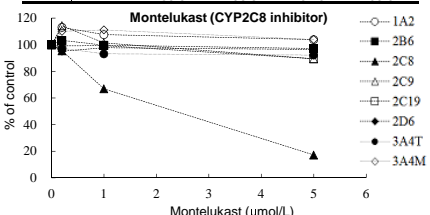
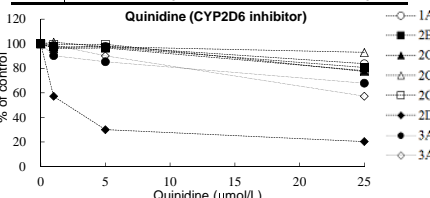


Table 7 and Figure 7 Inhibitory effect of Quinidine for each CYP isoform in HLM

Quinidine	% of control			
	CYP1A2	CYP2B6	CYP2C8	CYP2C9
0 $\mu\text{mol/L}$	100.0	100.0	100.0	100.0
1 $\mu\text{mol/L}$	96.7	97.6	96.2	101.3
5 $\mu\text{mol/L}$	97.9	97.3	96.5	97.6
25 $\mu\text{mol/L}$	83.7	80.7	78.0	93.0

Quinidine	% of control			
	CYP2C19	CYP2D6	CYP3A4 (T)	CYP3A4 (M)
0 $\mu\text{mol/L}$	100.0	100.0	100.0	100.0
1 $\mu\text{mol/L}$	99.5	57.2	90.3	99.0
5 $\mu\text{mol/L}$	99.5	30.0	85.4	90.1
25 $\mu\text{mol/L}$	77.3	20.3	67.9	57.2



Discussion

This study characterized the specificity of each CYP typical inhibitor under long term incubation in HLM. At 1 $\mu\text{mol/L}$ α -naphthoflavone (CYP1A2 inhibitor) showed specific inhibition to CYP1A2. However, at 25 $\mu\text{mol/L}$, it inhibited other CYP isoform activities (CYP2C8, CYP2C9, CYP2C19 and CYP3A4) (Table 1 and Figure 1). By contrast, furafylline (CYP1A2 inhibitor) showed specific inhibition to CYP1A2 at the maximum concentration (25 $\mu\text{mol/L}$) (Table 2 and Figure 2). At 10 $\mu\text{mol/L}$, thio-TEPA (CYP2B6 inhibitor) showed specific inhibition to CYP2B6 (Table 3 and Figure 3). At 5 $\mu\text{mol/L}$, montelukast (CYP2C8 inhibitor) showed specific inhibition to CYP2C8 (Table 4 and Figure 4). At 5 to 25 $\mu\text{mol/L}$, sulfaphenazole (CYP2C9 inhibitor) showed specific inhibition to CYP2C9 (Table 5 and Figure 5). At 5 $\mu\text{mol/L}$, benzylrinivorol (CYP2C19 inhibitor) showed specific inhibition to CYP2C19, but at 25 $\mu\text{mol/L}$, it inhibited other CYP isoform activities (CYP2B6, CYP2C8 and CYP3A4) (Table 6 and Figure 6). At 5 $\mu\text{mol/L}$, quinidine (CYP2D6 inhibitor) showed specific inhibition to CYP2D6 (Table 7 and Figure 7). At 1 $\mu\text{mol/L}$, ketoconazole (CYP3A4 inhibitor) showed specific inhibition to CYP3A4 (Table 8 and Figure 8). We also compared the IC_{50} values of the CYP typical inhibitors under the different incubation times (15 min and 60 min) (Table 9). As for the typical inhibitors of which the IC_{50} values after 60 min incubation were higher (α -naphthoflavone, montelukast, sulfaphenazole, benzylrinivorol, quinidine and ketoconazole), the inhibitory effects might be weakened by the metabolism of the typical inhibitors themselves. As for the typical inhibitors of which the IC_{50} values after 60 min incubation were lower (furafylline and thio-TEPA), time-dependent inhibition was indicated. In this presentation, we have confirmed the specificity and sustainability of CYP typical inhibitors under the tough condition (incubation time of inhibitor was 60 min, final concentration of HLM was 1 mg protein/mL). Therefore, these results will help us to identify metabolic enzymes of the low intrinsic clearance (CL_{int}) compounds metabolized by CYP isoforms.

References

- 1) Drug Metab. 2005;6(5):413-54.
- 2) Drug Metab Dispos. 2000;28(10):1176-83.
- 3) Br J Clin Pharmacol. 2000;49(4):343-51.
- 4) Drug Metab Dispos. 1995;23(1):154-8.
- 5) U.S. Food and Drug Administration. Guidance for Industry (DRAFT GUIDANCE): Drug Interaction Studies-Study Design, Data Analysis, and Implications for Dosing and Labeling (2006).