

PXB-Mice with Humanized Livers Services Provide Better Predictions of Human Pharmacokinetics Compared to CB17-SCID Mice

Chien-Chang Shen¹, Matthew Baginski², Wei-Chung Tseng¹, Yi-Kai Cheng¹, Hui-Erh Chang¹, Woan-Ing Twu¹, Ting-Jing Shen¹

- ¹ Pharmacology Discovery Services Taiwan, Ltd., Partner Lab of Eurofins Discovery I New Taipei City, Taiwan
- ² PhoenixBio USA Corporation I New York, USA

Abstract

Purpose

For drug discovery and pharmaceutical development, accurate predictions of human

PK Analysis of Naltrexone in PXB-Mice and CB17-SCID Mice

Α									
IV	t _{1/2} (h)	C ₀ (ng/mL)	AUC _{last} (h*ng/mL)	AUC _{Inf} (h*ng/mL)	AUC/D (h*kg*ng/mL/mg)	AUC Extr (%)	MRT (h)	Vss (L/kg)	CL (mL/min/kg)
CB17-SCID								_	

pharmacokinetics (PK), metabolite profiles, and drug-drug interactions (DDIs) are crucial for patient safety and drug efficacy. As the liver is the major organ participating in drug metabolism *in vivo*, several *in vitro* methods using human hepatic microsomes and hepatocytes, as well as chimeric mice with humanized livers have been employed for better preclinical predictions for drug metabolism. However, it is inconclusive as to which system provides the best predictions representing the human body using the *in vivo* platform. In this study, the aim is to investigate the translational value of PXB-mice, with humanized livers, in drug metabolism as a comparison to CB17-SCID mice by testing the PK profiles of published reference articles.

Methods

Test compounds, including amlodipine, naltrexone, and meloxicam, were administrated intravenously (IV) and orally (PO) to mice at 1 and 10 mg/kg, respectively. Blood aliquots were collected from three alternative mice at six or nine designed time points. The plasma samples were processed using acetonitrile (ACN) precipitation and analyzed by liquid chromatography-tandem mass spectrometry (LC-MS/MS). The fundamental PK parameters of each dosing compound after IV and PO administrations were obtained from the non-compartmental analysis (NCA) using WinNonlin[®]. The oral bioavailability (F%) was also calculated.

Results and Discussion

In mice administered with amlodipine, the F% in CB17-SCID mice (171%) was higher



Figure 2. The PK parameters and concentration-time profile of naltrexone after single IV (1 mg/kg) or PO (10 mg/kg) administration to PXB-mice and CB17-SCID mice. The plasma samples were obtained from these animals at 0.05, 0.5, 1, 2, 6, and 24 h after IV and 0.167, 0.5, 1, 2, 6, and 24 h after PO administration. A. The plasma PK parameters. B. Plasma concentration-time profile in PXB-mice. C. Plasma concentration-time profile in CB17-SCID mice.

PK Analysis of Meloxicam in PXB-Mice and CB17-SCID Mice

than that in PXB-mice (64%). In mice treated with naltrexone, the F% in PXB-mice (18%) was three times greater than in CB17-SCID mice (5%). Moreover, the F% of meloxicam was comparable in both strains (109% in CB17-SCID mice and 98% in PXB-mice). Taken together, the data demonstrate similarity and consistency with published oral bioavail-ability results in mouse and human PK, with respect to PXB-mice and CB17-SCID mice, suggesting PXB-mice serve as a reliable strain for use in human PK predictions.

Conclusions

PXB-mice provide high translational values with clinical relevance for predictions of human PK during the preclinical phase.

PK Analysis of Amlodipine in PXB-Mice and CB17-SCID Mice

Α													
IV	t _{1/2} (h)	C ₀ (ng/ml	AL _) (h*r	JC _{last} g/mL)	AUC _{Inf} (h*ng/mL)	ې (h*kg'	AUC/D *ng/mL/mg)	AUC Ext (%)	tr MI (ł	RT 1) (Vss _/kg)	CL (mL/min/kg)	
CB17-SCID mice	9.5	136	8	819	970		970	16	1	1	12	,	17
PXB-mice	9.4	73	3	887	455		455	15	1	1	24	3	37
PO	t _{1/2} (h)	T _{max} (h)	C _{max} (ng/mL)	AUC _{la} (h*ng/n	ast AU nL) (h*ng	C _{Inf} g/mL)	AUC/D (h*kg*ng/ mL/mg)	AUC Extr (%)	MRT (h)	Vz/F (L/kg)	C (ml	CL/F _/min/ kg)	F (%)
CB17-SCID mice	30.3	6.0	393	1357	3 165	549	1655	18	41	26		10	171

IV	t _{1/2} (h)	C ₀ (ng/mL	AU(.) (h*ng	AUC _{last} (h*ng/mL)		AUC _{Inf} (h*ng/mL)		AUC/D *ng/mL/mg)	AUC Ex (%)	tr N	MRT (h)		ss kg)	CL (mL/min/kg)		
CB17-SCID mice	2.4	2953	7532		9103		9103		17		3)	2		
PXB-mice	3.5	6135	333	888	33680		33680		1		3		0		0	
PO	t _{1/2} (h)	T _{max} (h)	C _{max} (ng/mL)	AUC (h*ng/	last mL)	AU((h*ng	C _{Inf} /mL)	AUC/D (h*kg*ng/ mL/mg)	AUC Extr (%)	MRT (h)	V. (L.	z/F /kg)	C (mL k	L/F /min/ (g)	F (%)	
CB17-SCID mice	2.5	0.5	26050	991	71 992		87	9929	0	3		0		2	109	
PXB-mice	5.6	1.0	42628	3134	-54	3302	289	33029	5	6		0		1	98	

В **PXB-Mice CB17-SCID** Mice entration (ng/mL) () 100000 g Meloxicam (10 mg/kg, PO) 100000 Meloxicam (10 mg/kg, PO) Meloxicam (1 mg/kg, IV))g∩ Meloxicam (1 mg/kg, IV) 0000 1000 100-100 Ŭ Plasma 30 20 30 20 Time (h) Time (h)

Figure 3. The PK parameters and concentration-time profile of meloxicam after single IV (1 mg/kg) or PO (10 mg/kg) administration to PXB-mice and CB17-SCID mice. The plasma samples were obtained from these animals at 0.05, 0.5, 1, 2, 6, and 24 h after IV and 0.167, 0.5, 1, 2, 6, and 24 h after PO administration. A. The plasma PK parameters. B. Plasma concentration-time profile in PXB-mice. C. Plasma concentration-time profile in CB17-SCID mice.





Figure 1. The PK parameters and concentration-time profile of amlodipine after single IV (1 mg/kg) or PO (10 mg/kg) administration to PXB-mice and CB17-SCID mice. The plasma samples were obtained from these animals at 0.05, 0.5, 1, 2, 6, and 24 hours (h) after IV and 0.167, 0.5, 1, 2, 6, 24, 36, 48, and 72 h after PO administration. A. The plasma PK parameters. B. Plasma concentration-time profile in PXB-mice. C. Plasma concentration-time profile in CB17-SCID mice.

Summary

	PXB-Mice	CB17-SCID Mice
Amlodipine	Moderate F% as expected	High F% as expected
Naltrexone	Relatively higher F% as expected	Relatively lower F% as expected
Meloxicam	High F% as expected	High F% as expected

- 1. CB17-SCID mice display higher F% for amlodipine, while PXB-mice show superior values for naltrexone.
- 2. For meloxicam, F% remains consistent across both strains.
- 3. PXB-mice provide a suitable model to study human PK, as evidenced by moderate F% in naltrexone and high F% in meloxicam.

References

- Stopher DA, Beresford AP, Macrae PV, Humphrey MJ. The metabolism and pharmacokinetics of amlodipine in humans and animals. *J Cardiovasc Pharmacol*. 1988;12 Suppl 7:S55–9. doi: 10.1097/00005344-198812007-00012. PMID: 2467130.
- 2. Licko V. Overview of human pharmacokinetics of naltrexone. NIDA Res Monogr. 1981;28:161–71. PMID: 6791002.
- Busch U, Schmid J, Heinzel G, Schmaus H, Baierl J, Huber C, Roth W. Pharmacokinetics of meloxicam in animals and the relevance to humans. *Drug Metab Dispos*. 1998 Jun;26(6):576–84.
 PMID: 9616195.