

**INTEGRATED *IN VITRO* AND *IN VIVO***  
**APPROACHES TO DRUG METABOLISM**  
**INVESTIGATION - A STUDY CASE**

---

Massimiliano Fonsi

**Meet the Experts Transporter Conference**  
**Seoul - November 14 2019**

EVERY STEP OF THE WAY

# Introduction

An ideal *in vitro* DMPK model should include all the pathways that are relevant for the PK *in vivo*. Quantitative scaling to whole organ/body is essential for the prediction of exposure of parent drug and its metabolites *in vivo*.

The liver is generally considered the organ most involved in metabolic transformations and *in vitro* screening systems are based essentially on liver tissue fractions. More frequently, metabolism in the gut is also considered. Conversely, metabolism in other organs is rarely taken into consideration.

The case study presented today will offer an example of

- 1) The experimental evidence that should trigger the investigation of extrahepatic metabolism
- 2) The differences that may exist between preclinical species and humans\*
- 3) The impact of extrahepatic metabolism on screening cascades and decisions process in a drug discovery project.

\*Extrahepatic metabolism may complicate the IMVC in rats.  
Fonsi M - Drug Metab Lett. 2014;8(1):51-66.

# Discovery of MK-4827 (Niraparib)

*Poly(ADP-ribose)polymerase (PARP) as a context specific anti-cancer target*

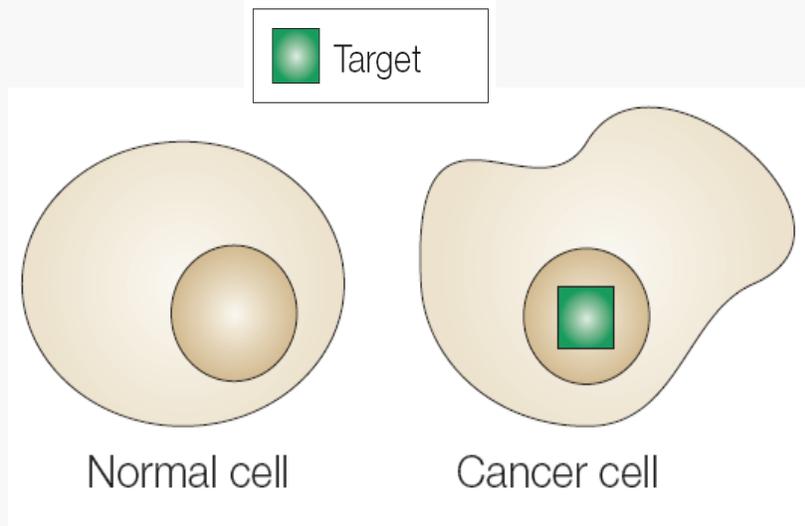


# Can we selectively kill tumor cells? The Synthetic Lethality approach

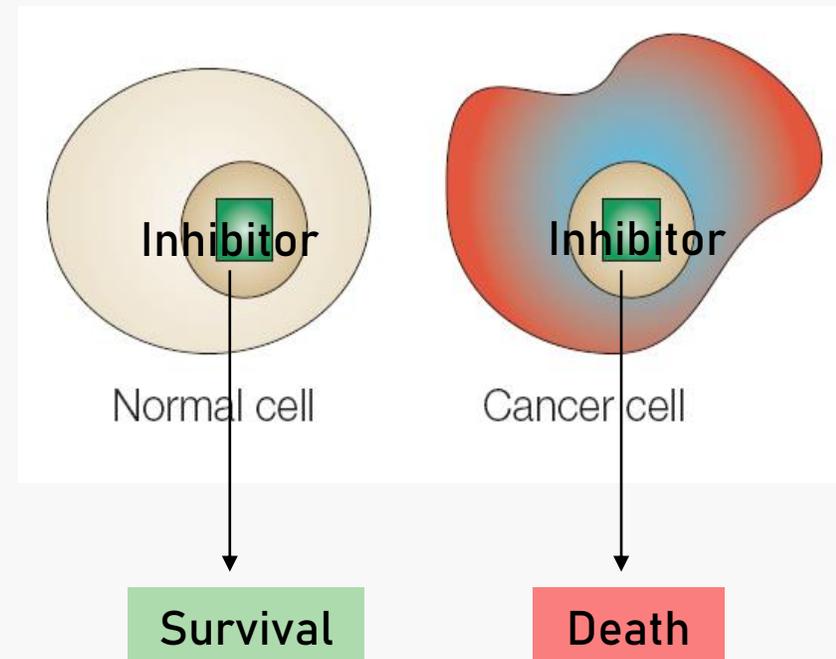
- Most anticancer drugs in use today have a low therapeutic index
- Two strategies to improve high therapeutic index

## *Target-driven*

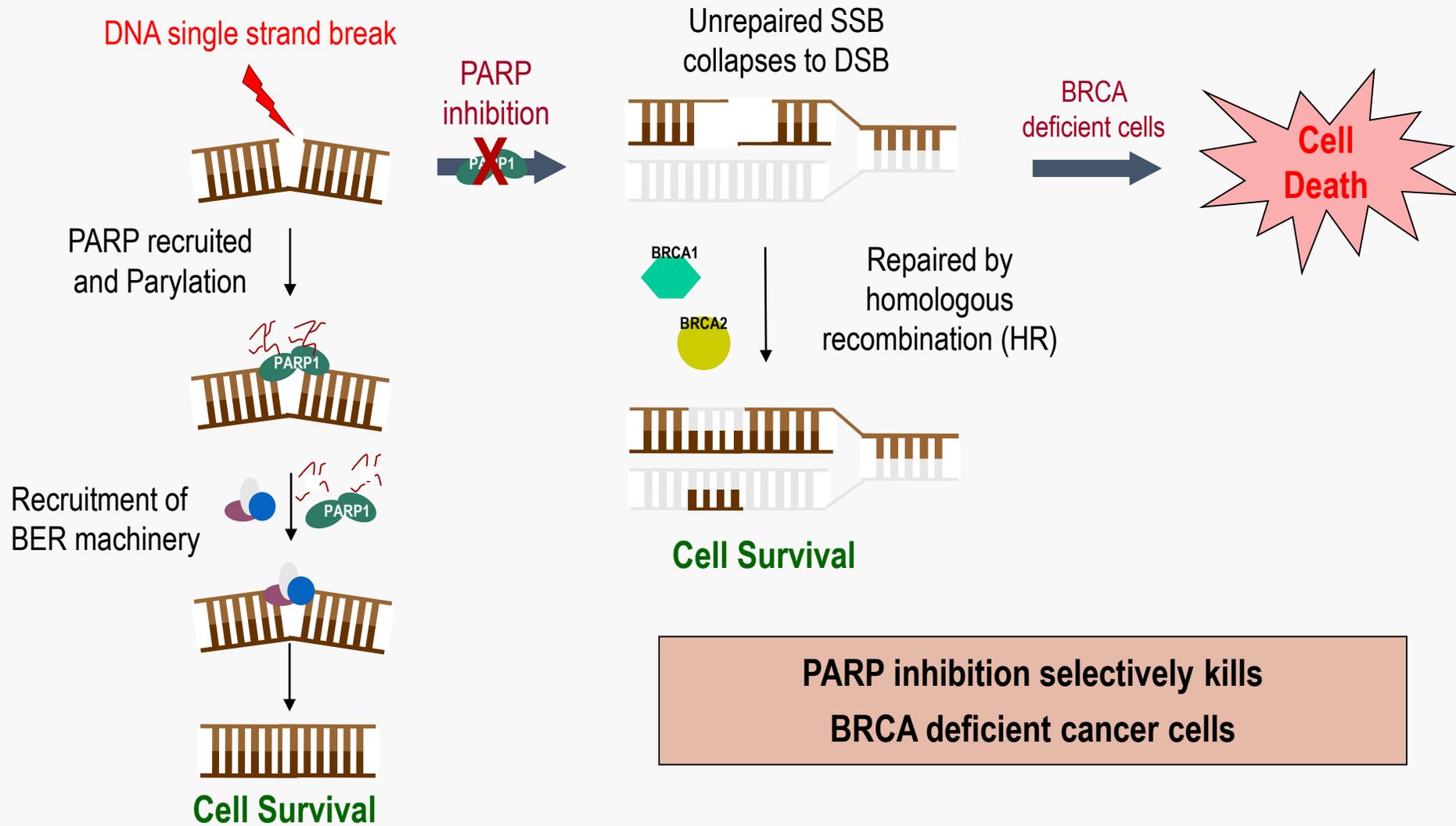
- Target only in cancer cells



## *Context-driven*

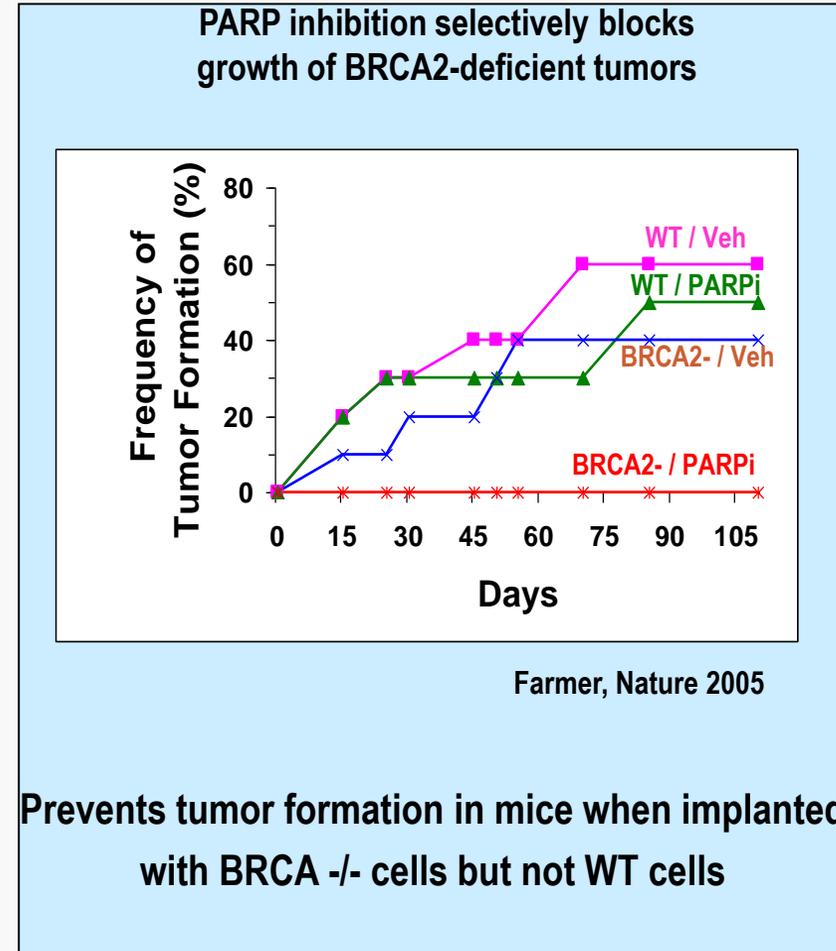
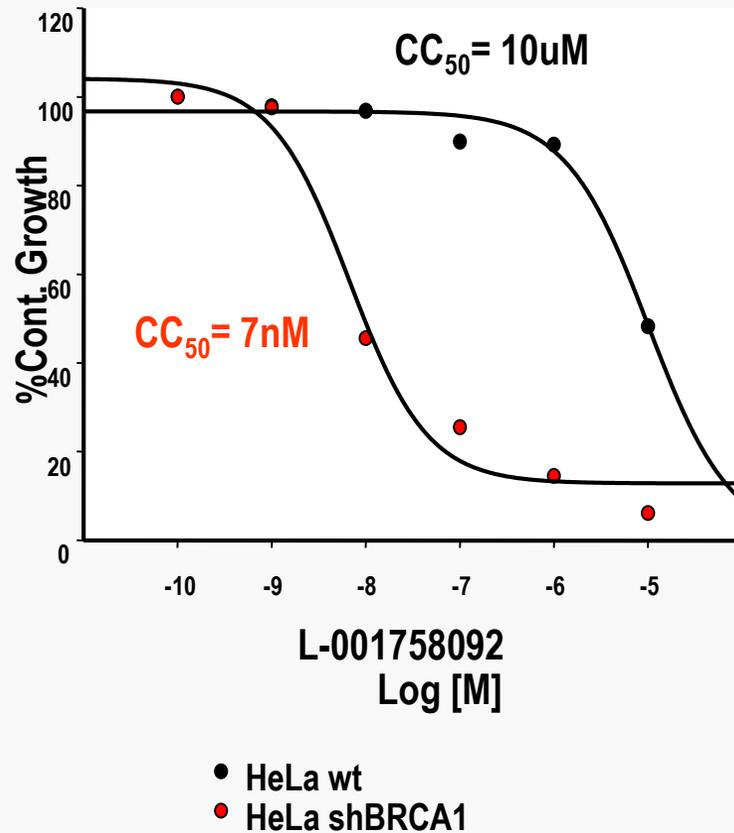


# PARPi as context-specific anti-cancer agents



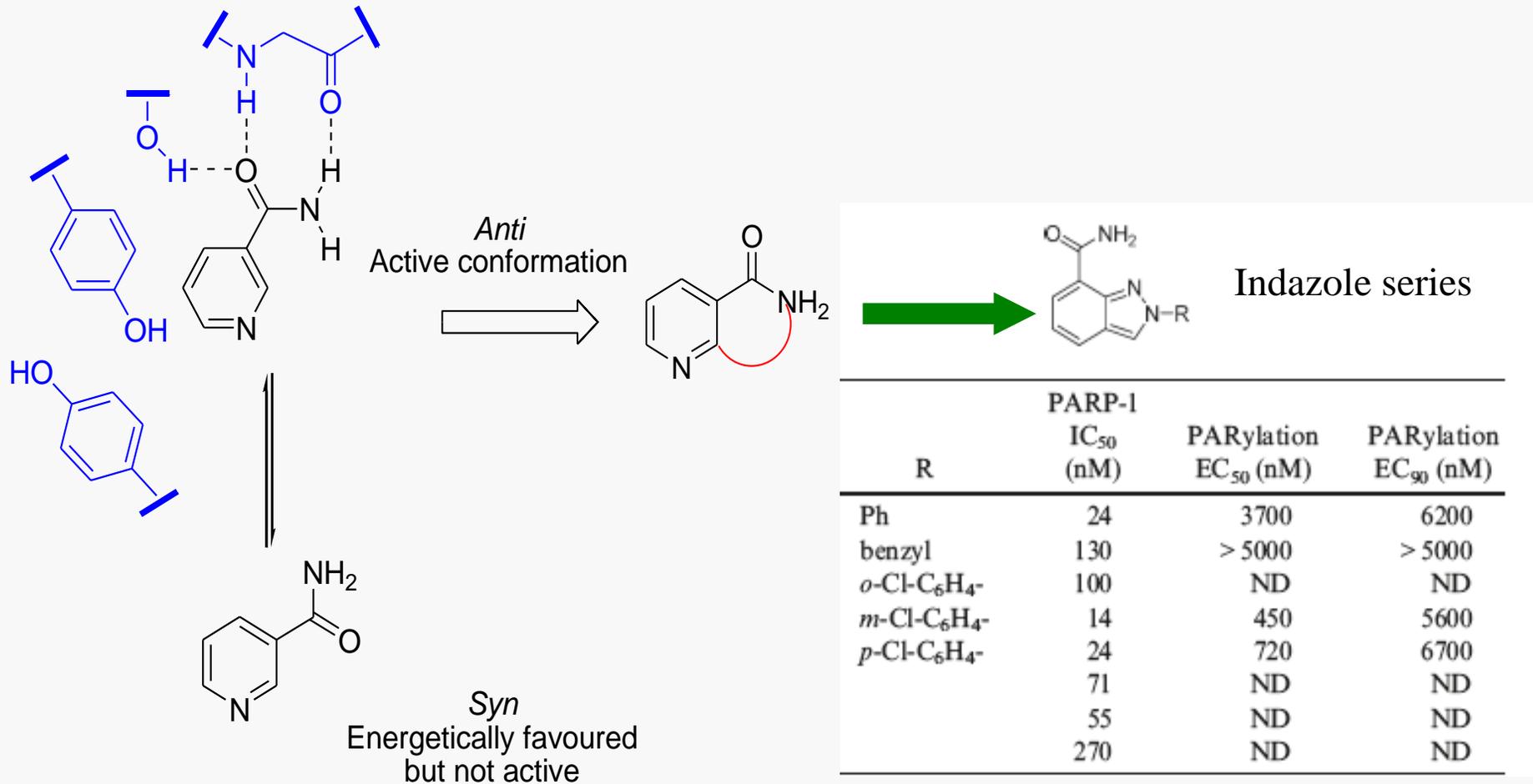
# PARPi's in BRCA-deficient cells

Selectively kill BRCA-deficient cells *in vitro*

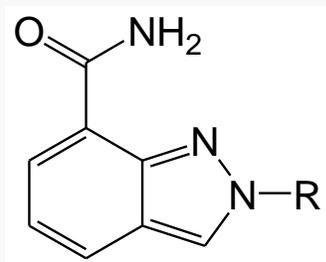


# First Generation of PARP inhibitors

- First PARPi were based on the cleavage product - nicotinamide



# Initial profiling of the prototype PARP inh: early detection of series related liabilities



Where R= Ph

**Solubility** 0.02 mg/mL (water);  
**LogP** = 2.4 – **TPSA**=61

RLM Cl <sub>int</sub> ((μL/min)/ mgP)	HLM Cl <sub>int</sub> ((μL/min)/ mgP)	rat Cl ((mL/min)/ kg)	F (%)
123	138	30	41

**Permeability (P<sub>app</sub>) = 5 (cm\*10E-6/s) (LLC-PK1);**

**ER (BA/AB in LLC-PK1 humMDR1) = 12**

**ER (BA/AB in LLC-PK1 mouseMdr1) = 17**

### Enzyme Inhibition

#### **CYP450 (IC<sub>50</sub>)**

3A4, 2D6, 2C9, 2C19, 1A2 >50 μM

### Enzyme Induction

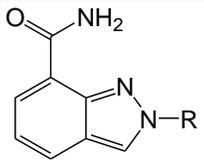
hPXR 74% rifampicin effect @ 25 μM (weak inducer)

No CYP3A4 induction in HH @ 30 μM

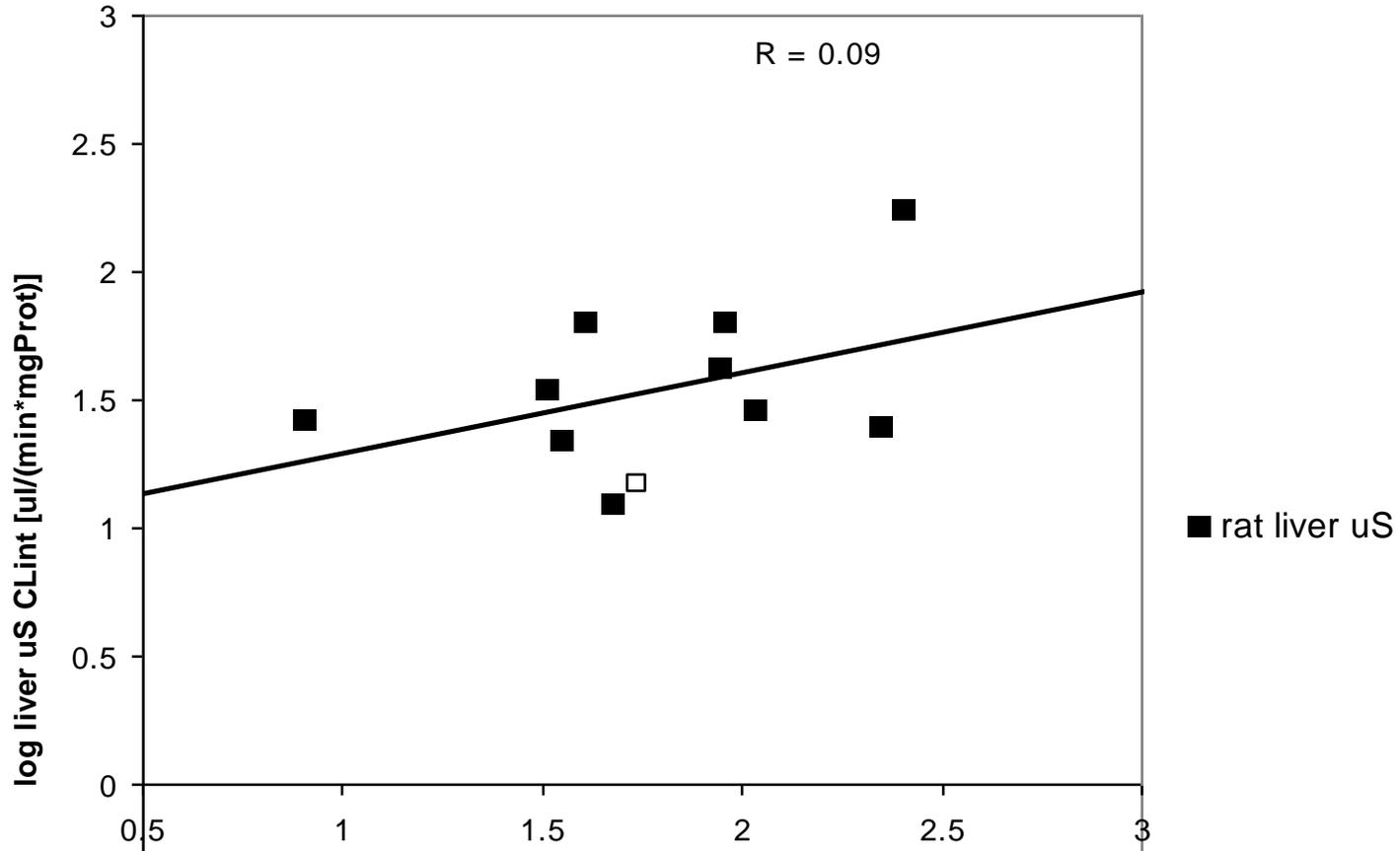
### **Off-Target**

hERG  
Na<sub>v</sub>1.5 (0.2 + 3 Hz)

9.7 μM  
19μM



### Rat liver microsomes



No better correlation using Hepatocyte  
Cpds were stable in rat plasma  
No improvements correcting CLp by B-to-P ratio or by renal CL

# Investigating mechanism of extrahepatic CL in vivo of cpd 10...

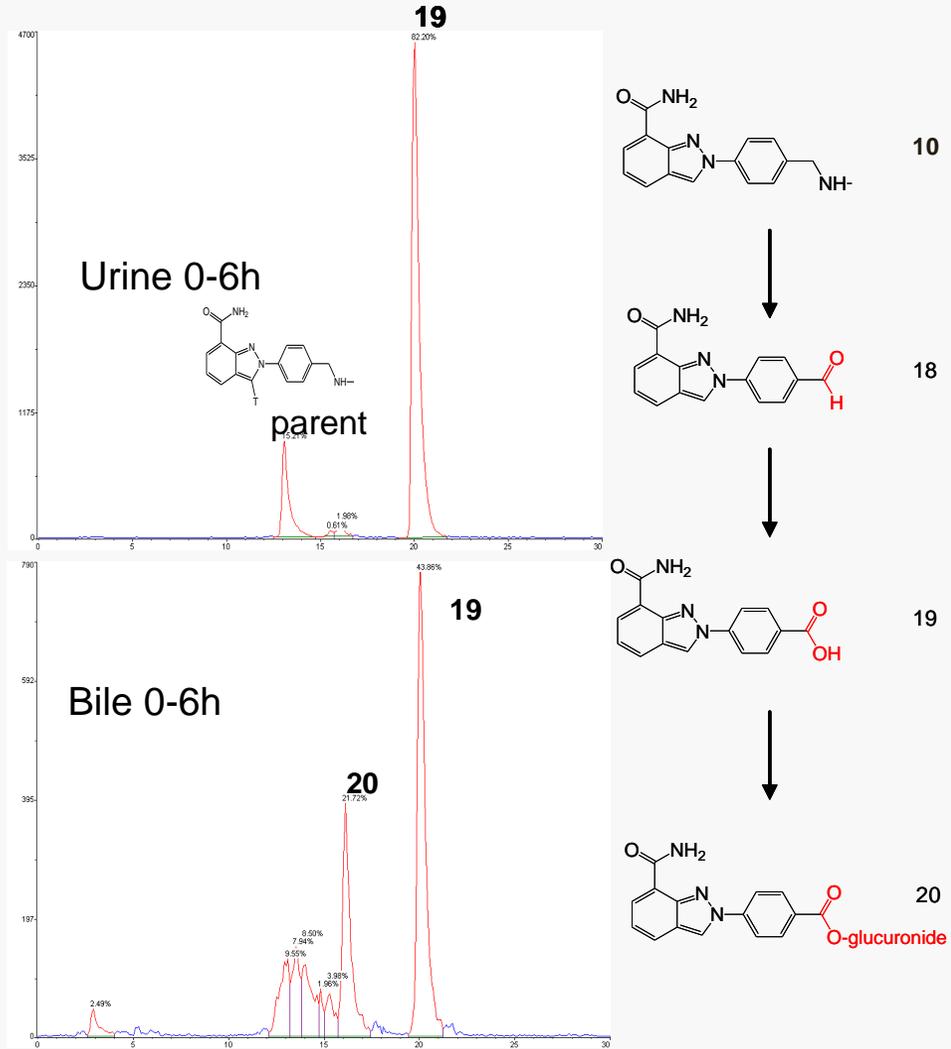
IV (treatment) Dose: 3 mg/kg  
Formulation: DMSO/PEG400/Water (20%/60%/20%)

		Mean	std. Dev.
AUC <sub>(0-∞)</sub>	μM·hr	<b>0,83</b>	0,08
Cl	mL/min/kg	<b>217</b>	20
t <sub>1/2</sub>	hr	<b>1,4</b>	0,2
Vd <sub>ss</sub>	L/kg	<b>22,1</b>	0,8

PO (treatment) Dose: 3 mg/kg  
Formulation: PEG 400 (SUSPENSION)

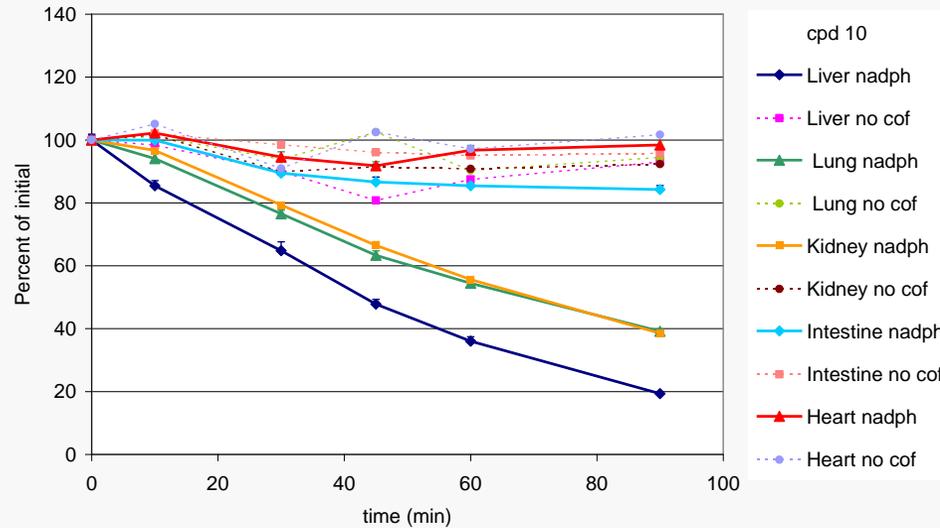
		Mean	std. Dev.
C <sub>max</sub>	μM	<b>0,07</b>	0,04
T <sub>max</sub>	hr	<b>2,2</b>	2,5
AUC <sub>(0-∞)</sub>	μM·hr	<b>0,37</b>	0,14
t <sub>1/2</sub>	hr	<b>5,4</b>	3,2
F <sub>(0-∞)</sub>	%	<b>45</b>	17,0

	% of dosed radioactivity	% of the dose recovery transformed into 19 or 20
dose	100%	
urine 0-6h	20%	19%
urine6-24h	21%	19%
bile 0-6h	30%	16%
bile 6-24h	10%	6%
<b>recovery</b>	<b>81%</b>	<b>59,4%</b>

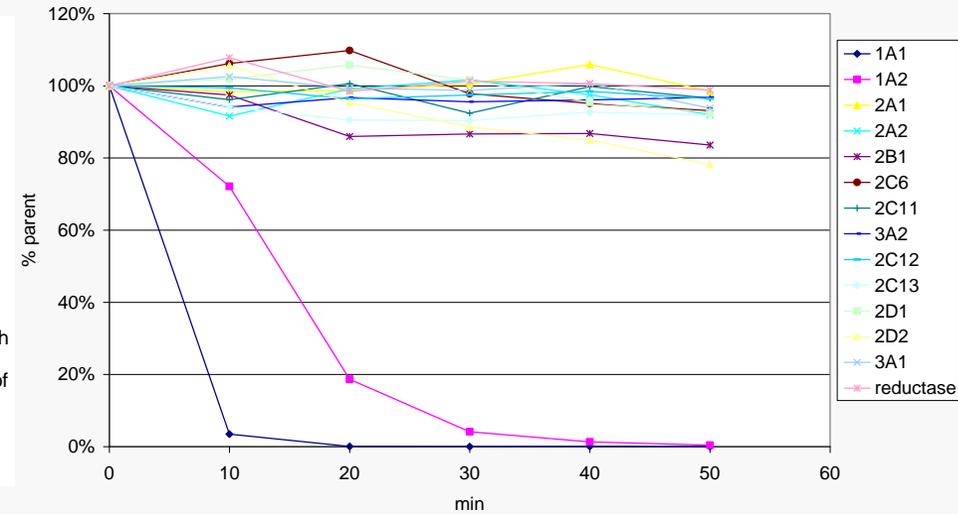


# Trying to establish an in vitro model of extrahepatic metabolism

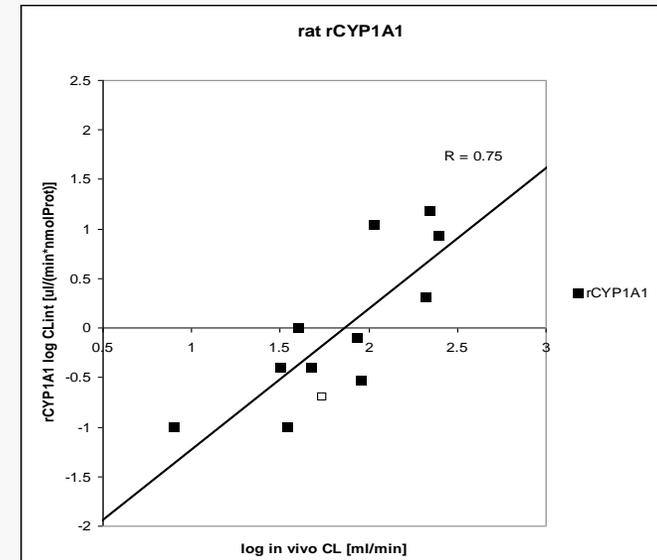
Metabolic stability in liver, lung, kidney, intestine microsomes and heart homogenate in presence of NADPH



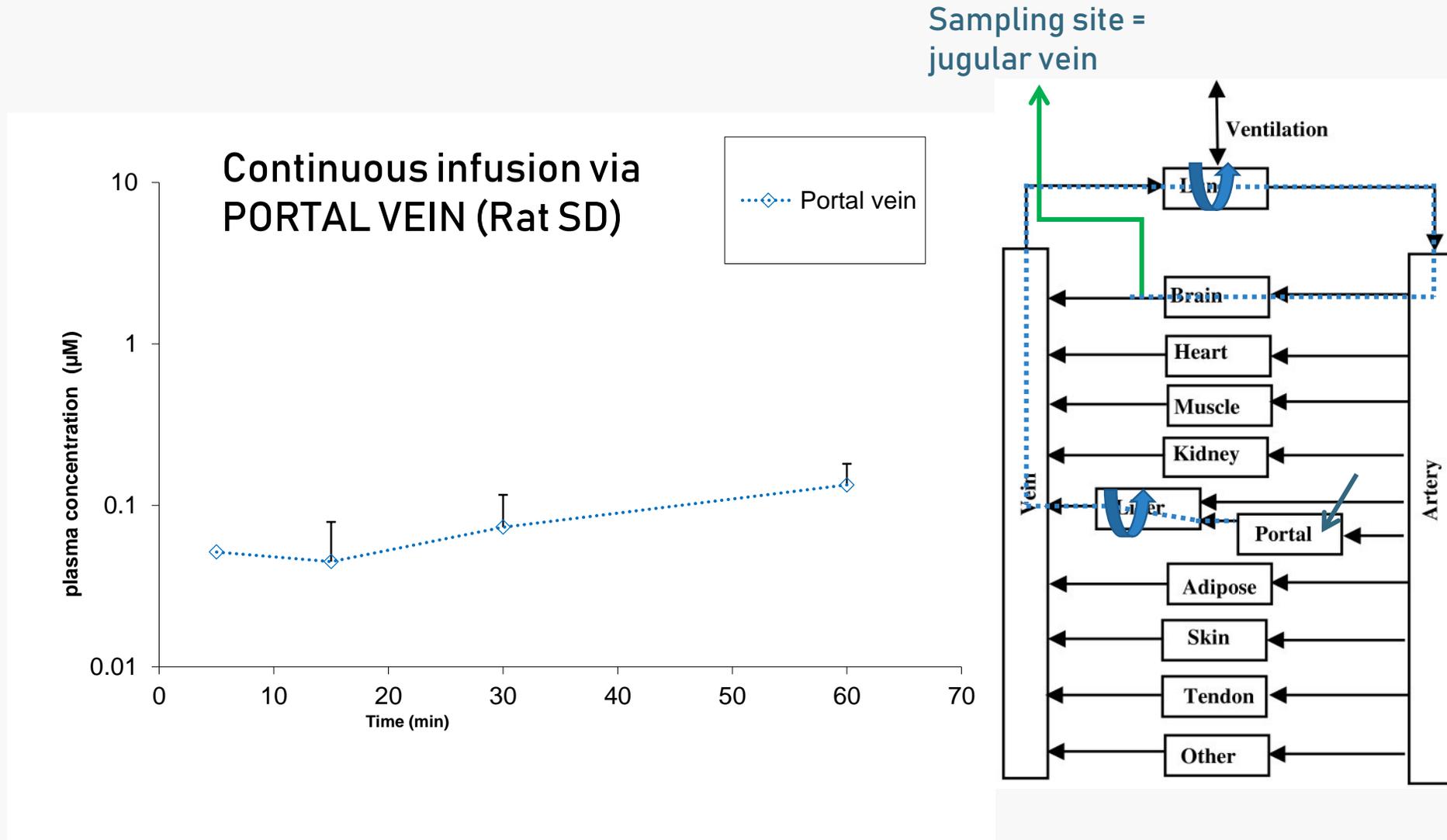
rat CYP phenotyping for cpd 10

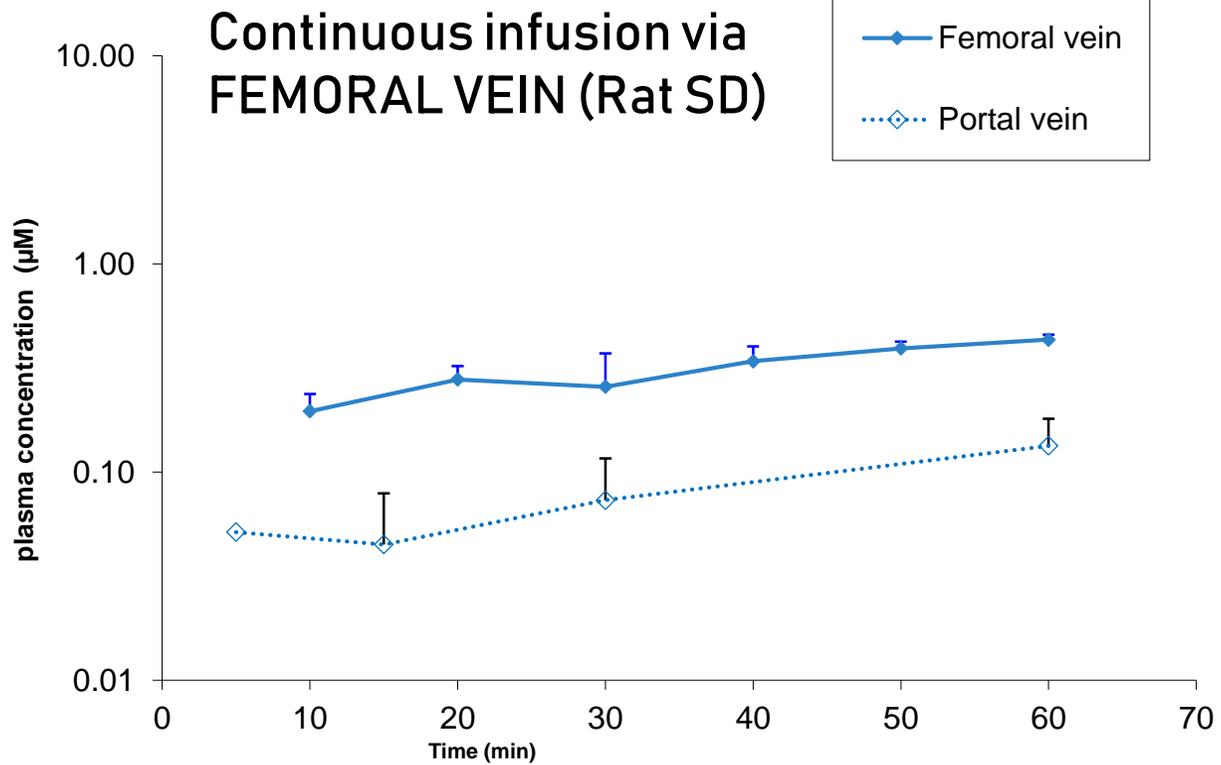


compd	RLM $Cl_{int}$ (( $\mu$ L/min)/ mgP)	HLM $Cl_{int}$ (( $\mu$ L/min)/ mgP)	CYP1 a1 $Cl_{int}$ (( $\mu$ L/min)/ mgP)	rat Cl ((mL/min)/ kg)
8	123	138	ND	30
37	177	1	8.7	450
39	29	3	3.8	107
41	34	2	<0.1	8
43	28	3	6	220
47	18	1	0.8	131
54	32	5	5.4	69
48	9	1	2.2	58
50	11	<1	0.4	30
13	52	4	0.7	87
14	36	3	0.7	30
15	22	4	2.5	ND
16	11	1	0.3	47
17	15	5	<0.1	ND
56	16	4	0.3	28
57	7	3	<0.1	24

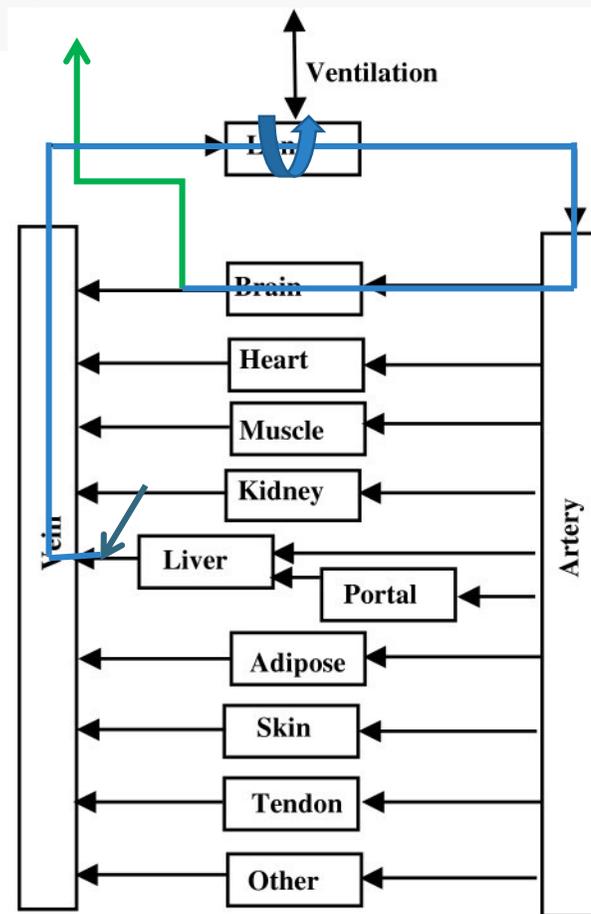


# Mechanistic in vivo PK to confirm hepatic and liver organ extraction

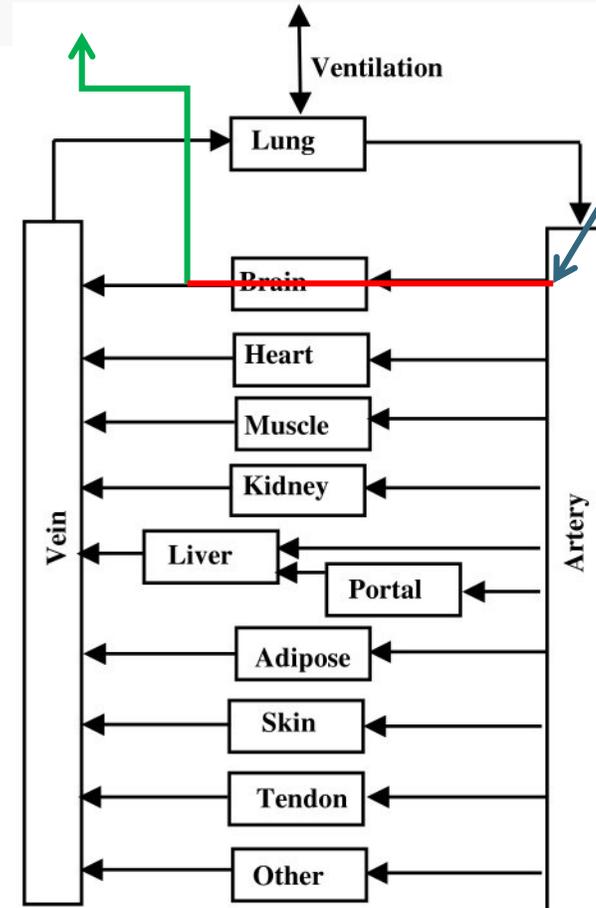
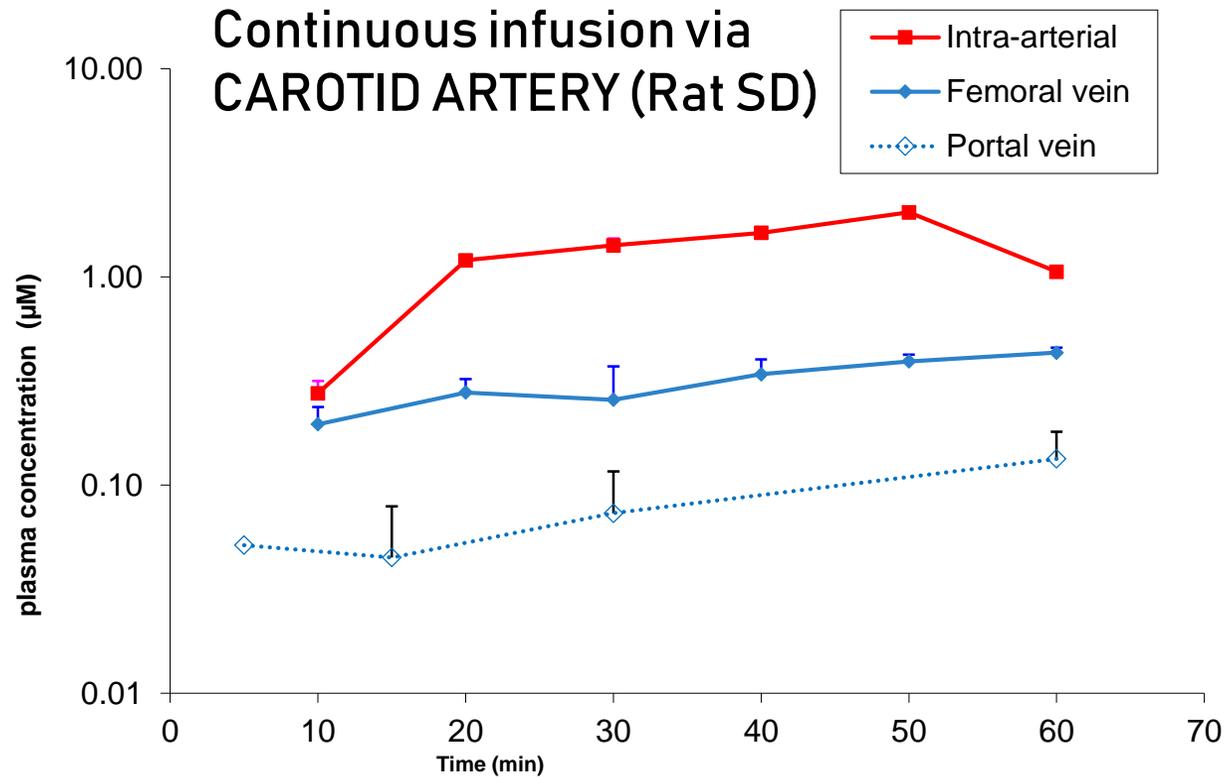




Sampling site =  
jugular vein

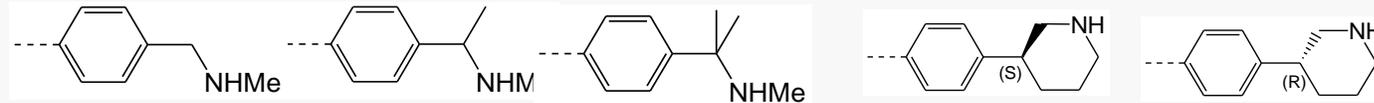
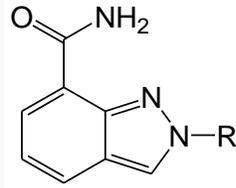


Sampling site =  
jugular vein



$C_{SS}^{femoral}$	$\mu\text{M}$	$0.52 \pm 0.06$
$Cl_p^*$	$\text{ml}/\text{min}/\text{kg}$	$140 \pm 16$
$C_{SS}^{portal}$	$\mu\text{M}$	$0.21 \pm 0.06$
$E\%_{hepatic}$	%	$59 \pm 19$
$C_{SS}^{arterial}$	$\mu\text{M}$	$2.0 \pm 0.5$
$E\%_{lung}$	%	$78 \pm 23$

# SPR *in vitro* & *in vivo*



CL<sub>int</sub>  
rCYP1a1  
(uL/min/nmol)

6000

2200

400

300

<100

CL<sub>b</sub>  
(mL/min/kg)

220

58

30

46

24

PARylation EC<sub>90</sub> (nM)

30

280

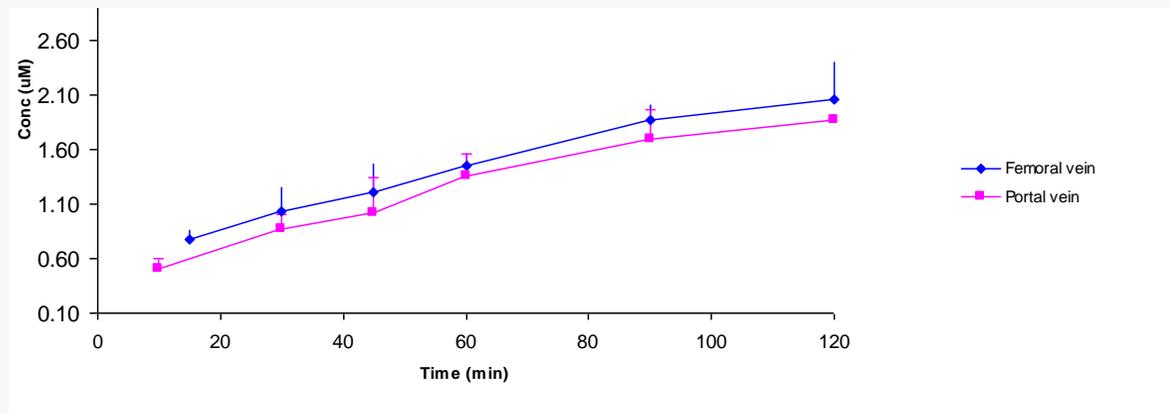


MK-4827 (Niraparib)

[J Med Chem.](#) 2009 Nov 26;52(22):7170-85.

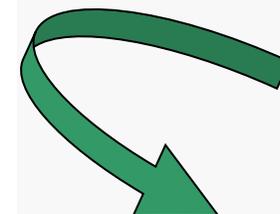
# MK-4827 IVIVC: Rat liver data

TOTAL blood CL = 46 mL/min/Kgbw – F% = 92%



Hep Intrinsic Clearance  
(µl/min/million cells)

Rat 2.2  
Hum 1.6



Predicted Rat hepatic CL<sup>1</sup>

ER<sub>H</sub> = 11%  
CL<sub>H</sub> = 7.1 (ml/min)/kg



Predicted hum hepatic CL<sup>1</sup>

ER<sub>H</sub> = 14%  
CL<sub>H</sub> = 2.8 (ml/min)/kg

**Infusion via Femoral vein**

Time (min)	Concentration (µM)			Average	Std. Dev.
	Rat 1	Rat 2	Rat 3		
15	0.80	0.85	0.67	0.78	0.09
30	1.16	1.17	0.79	1.04	0.22
45	1.36	1.36	0.90	1.21	0.26
60	1.53	1.52	1.30	1.45	0.13
90	1.89	2.00	1.69	1.86	0.15
120	2.29	2.22	1.64	2.05	0.36

**Infusion via Portal vein**

Time (h)	Concentration (µM)			Average	Std. Dev.
	Rat 4	Rat 5	Rat 6		
15	0.51	0.41	0.59	0.50	0.09
30	0.84	0.78	0.97	0.86	0.09
45	1.02	0.87	1.15	1.01	0.14
60	1.61	0.99	1.48	1.36	0.33
90	1.82	1.47	1.79	1.69	0.19
120	1.85	1.60	2.15	1.87	0.28

AUC (µM \* h)      174.8    177.9    140.5    **164.4**    20.8

AUC (µM \* h)      152.9    121.7    159.2    **144.6**    20.1

C<sub>ss</sub> (µM)  
Average 60 and 120 min      **1.8**

C<sub>ss</sub> (µM)      **1.6**

ER<sub>H</sub> = 10%

Measured hepatic CL (*in vivo*)  
CL<sub>H</sub> = ER<sub>H</sub> \* Q<sub>H</sub> = 6.8 ((ml/min)/kg)<sup>1</sup>

<sup>1</sup> Rat Q<sub>H</sub> = 68 ((ml/min)/kg) - Hum Q<sub>H</sub> = 20 ((ml/min)/kg).

Ring B.J.et al. *Pharmaceutical Sciences*, 2011 DOI 10.1002/jps

# Mechanism of Clearance in rats

TOTAL blood CL = 46 mL/min/Kgbw – F% = 92%

- 35% urinary excretion

=>  $CL_R = 0.35 * CL_B = 0.35 * 46 = 16.1$  ml/min/kg ; >  $GFR_{rat}$  (5-6 ml/min/kg)-Rat Kidney BF (40-52 ml/min/kg)

- 2 % biliary excretion
- 10 % hepatic extraction (determined by portal/femoral infusion)
- ~ 50 % extrahepatic metabolism

# Transporters Data

Compound	$P_{app}^{B-A}/P_{app}^{A-B}$					
	LLC-PK1 Papp (cm*10E-6/s)	Human MDR1	Mouse Mdr1a	Rat Mdr1a	Beagle MDR1	Monkey MDR1
MK-4827 (5 $\mu$ M)	3	12	22	17	10	ND
Verapamil (1 $\mu$ M)	1	6	12	7	4	ND

**MK-4827 is a Pgp substrate**

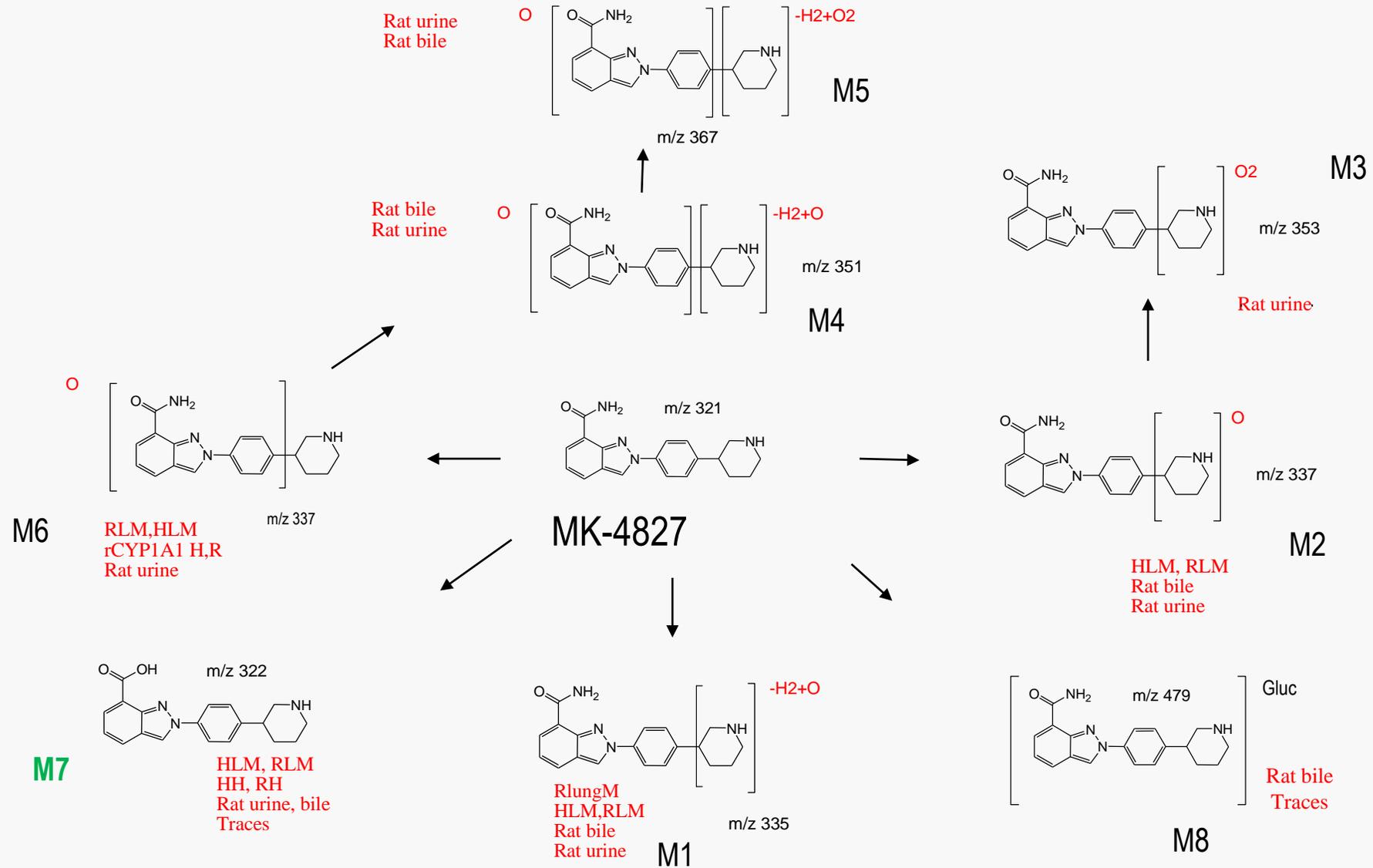
**Solubility >20 mg/mL**

# Turnover in lung and kidney preparations

MK-4827 in NADPH-supplemented preparations:

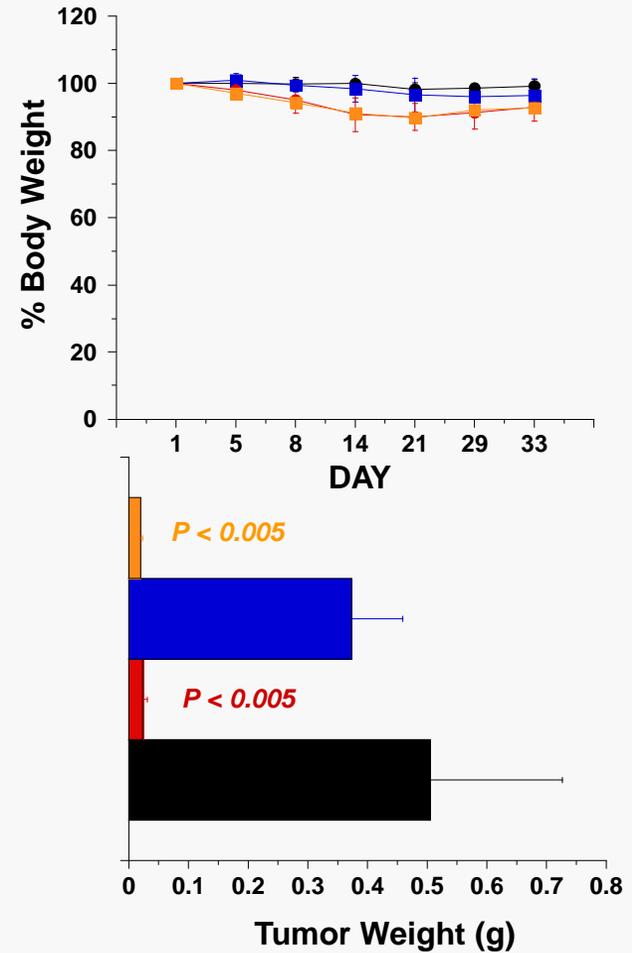
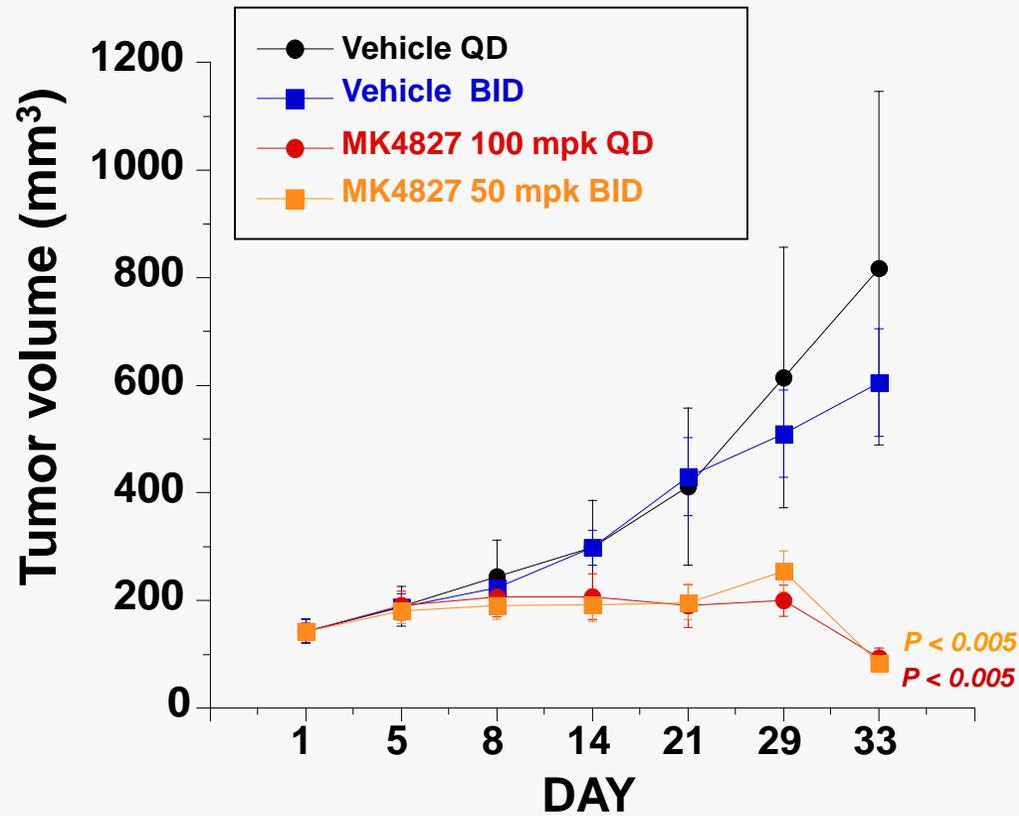
- rat lung microsomes (40 % turnover in 90 min)
  - rat kidney S9 fraction (50% turnover in 90 min)
  - human lung microsomes, from both non-smokers and smokers (stable)
  - human kidney S9 fraction (stable).
- 
- These results suggest low contribution of extra-hepatic metabolism in humans.
  - Predicted low hepatic CL in human (IVIVE using HH)
  - Potential renal CL in human due to Low PPB – Mdr1 active efflux
  - Expected good F%

# Proposed Metabolite Scheme (Rat and Human)



# Activity in MDA-MB-436 (BRCA1-) xenografts

Administered orally



Approved as PCC April 2007 (MK-4827)

## MK-4827 (Niraparib) preclinical profiling - PK

IV (male, n=3, saline)

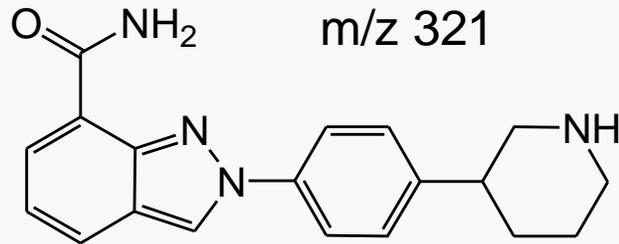
Species	Dose (mg/kg)	Cl <sub>b</sub> (mL/(min*kg))	Vd (L/kg)	T <sub>1/2</sub> (hr)	AUC (μM*hr) <sub>0-x</sub>
Rat	3	46±3	22.8±8.7	5.1±2.0	2.9±0.1
Dog	1	89±6	29.3±5.2	5.3±0.3	0.6±0.0
Monkey	1	17.5±6	17.1±2.5	6.8±1.0	1.5±0.2
		±	±	±	±

P.O. (male, n=3, water)

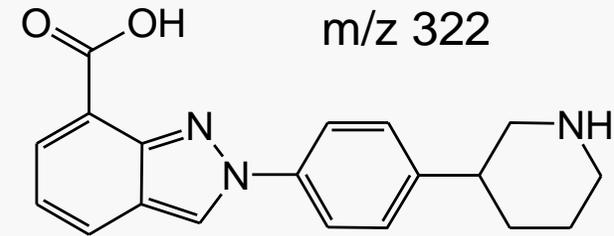
Species	Dose (mg/kg)	C <sub>max</sub> (μM)	T <sub>max</sub> (hr)	AUC (μM*hr) <sub>0-x</sub>	%F
Rat	4	0.3±0.1	4.0±1.0	3.7±1.8	92±29
Dog	2	0.2±0.0	0.9±0.3	0.8±0.1	80±14
Monkey	2	0.3±0.1	1.3±0.6	1.9±0.5	70±17
		±	±	±	±

**Dose proportional PK in rat – MK-4827**

# Proposed Metabolite Scheme (Dog)



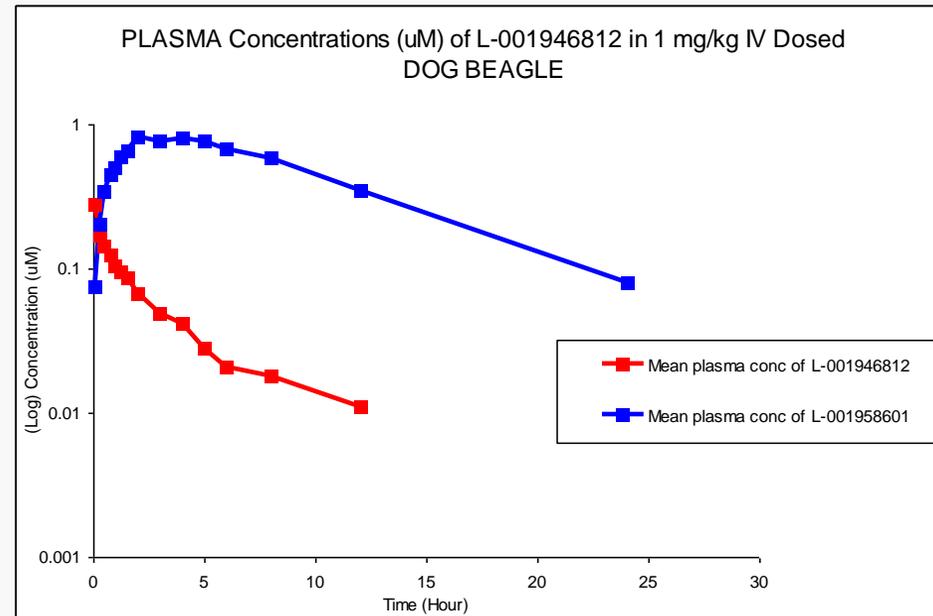
MK-4827



L-00958601

M7

hepatocytes  
microsomes  
urine  
plasma



# Rat/Dog Excretion Study 3 mg/kg iv (n=3)

	Percent of Dose Recovered (0-48h)		
	urine	bile (0-24h)	total
Total	48±1	32 ±4	80±4
Parent	35±2	2±2	37±3
Water	2±1	2±1	

Data obtained using MK4827 labeled with tritium in position 3 of the indazole ring.  
Low tritium loss was detected both in vivo and in vitro

In dogs only 2 % of parent excreted in urine



# Hum CL – t1/2 prediction for MK4827

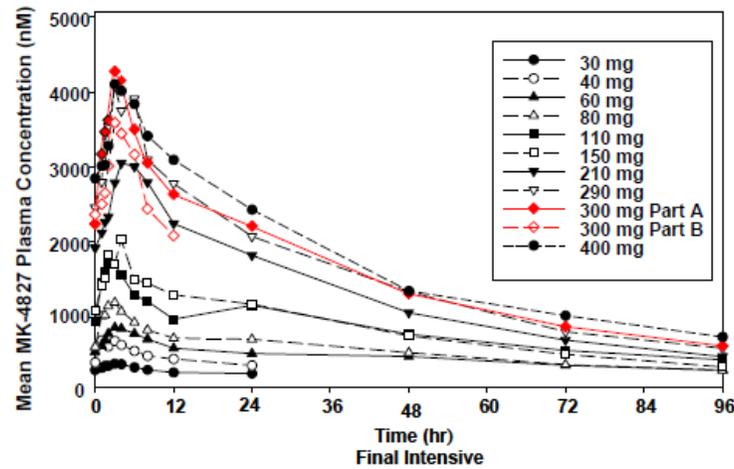
- Total hum CL =  $CL_{\text{metab liver}} + CL_R + CL_{\text{metab extrahep}}$ 
  - $CL_{\text{metab liver}}$ : IVIV scaling from HH = 2.8 ml/min/kg (range 1.8-3.8)
  - $CL_R = CL_{R\text{-rat}} * (f_{u\text{hum}} / f_{u\text{rat}}) * KBF_{\text{hum}} / KBF_{\text{rat}} * ER_{\text{hum}}/ER_{\text{rat}} = 4.5 \text{ ml/min/kg}$   
(range 3.5-7)
  - $CL_{\text{metab extrahep}}$ : expected to be negligible in human
  - $Vd_{ss} = Vd_{ss \text{ hum}} = Vd (\text{animal}) * f_u(\text{hum})/f_u(\text{animal}) = 26 \text{ (L/kg)}$  range (17-37)

		Vd (L/kg)		
		17	26	37
total		min	avrg	max
CL	5.5	28	43	61
(mL/min/kg)	8.3	21	<b>36</b>	46
	11	16	25	36

Predicted half-life (hr)

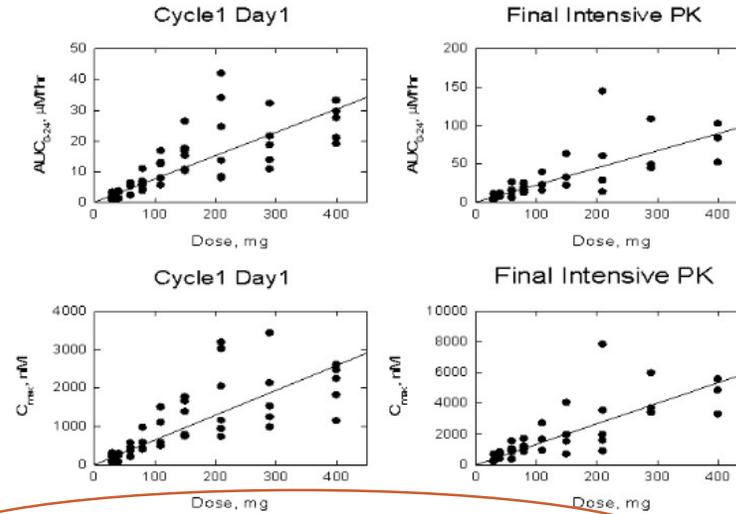
# Human PK data (Phase II)

Fig 1: Pharmacokinetic (PK) profile



- Bio-Exponential PK; PK parameters were dose proportional
- The terminal elimination  $t_{1/2}$  was estimated at 30-40 hours
- Two to four-fold accumulation was observed with once daily dosing

## Pharmacokinetic (PK) Profile



- Terminal half-life of MK-4827 estimated to be 37-42 hours
- Steady state trough levels approached following one 21-day cycle
- Two-four-fold accumulation in AUC0-24, Cmax and C24 over the 1st cycle
- Dose proportional pharmacokinetics

<http://meetinglibrary.asco.org/content/62393?format=posterImg>

[Niraparib](#) was approved in 2017

# Conclusion

Case studies in drug discovery have been reported showing relevant lungs and generalized extrahepatic metabolism in rats. The large extrahepatic contribution appeared to be rat specific with no analogy to human metabolism. Rat CYP1a1 was identified as the major enzyme responsible

The extraction fraction (CL/Q) in the lung is in general small, as a result of its modest metabolic activity but due to the fact that the lung blood flow is equivalent to the total cardiac output, in some cases, drug metabolism in the lungs can be substantial.

In general, xenobiotic metabolizing enzymes (XMEs) are expressed in lower levels in the extrahepatic tissues than in the liver, making the former less relevant for the clearance of xenobiotics. **Local metabolism, however, may lead to tissue-specific adverse responses, e.g. organ toxicities, allergies or cancer.\***

Fonsi M - Drug Metab Lett. 2014;8(1):51-66.

\* U. Gundert-Remy et al. Drug Metab Rev, Early Online: 1-34

**Thank you very much!**