

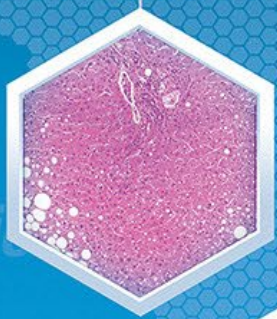
**XENOTECH**

A BioIVT Company

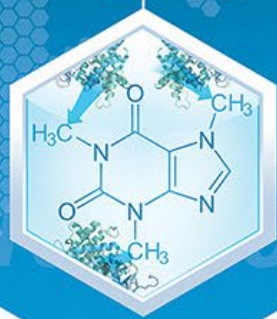
*25 Years*

# ADME | DRUG-DRUG INTERACTION | DMPK CONTRACT RESEARCH & TEST SYSTEM EXPERTISE

Cell & Tissue-Based Products



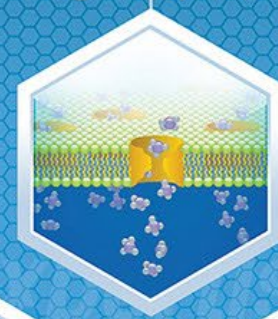
Reaction Phenotyping



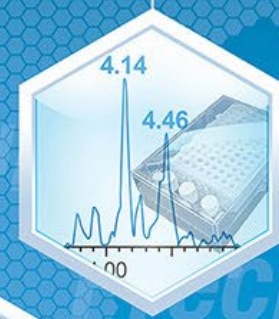
Enzyme Induction & Inhibition



Drug Transporters



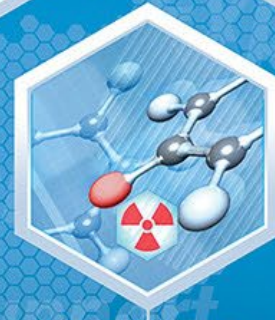
Metabolite ID & Production



Screening



Pharmacokinetics & QWBA



Radiolabeling



Bioanalytical

Welcome to the webinar...  
We will begin shortly

# Design Intricacies and Decision-Making Strategy for Drug-Drug Interaction Studies



**Andrew G. Taylor, Ph.D.**  
Manager of Technical Support for Services

# Design Intricacies and Decision-Making Strategy for DDI Studies

6 December 2022

Andrew G. Taylor, Ph.D.



# Outline



- **Why run these studies?**
- **Predominant types of *in vitro* ADME & Drug-Drug Interaction (DDI) studies & test systems**
- **Importance of proper study design & data interpretation**
- **When to conduct *in vitro* DDI ADME studies?**
- **Additional highlights and information**

# Regulatory Guidance

## FDA: Final January 2020


**In Vitro Drug  
Interaction Studies —  
Cytochrome P450  
Enzyme- and  
Transporter-Mediated  
Drug Interactions  
Guidance for Industry**

Additional copies are available from:  
Office of Communications, Division of Drug Information  
Center for Drug Evaluation and Research  
Food and Drug Administration  
10001 New Hampshire Ave., Hillandale Bldg., 4<sup>th</sup> Floor  
Silver Spring, MD 20993-0002  
Phone: 855-543-3784 or 301-796-3400; Fax: 301-431-6353  
Email: [druginfo@fda.hhs.gov](mailto:druginfo@fda.hhs.gov)  
<https://www.fda.gov/Drugs/GuidanceComplianceRegulatoryInformation/Guidances/default.htm>

**U.S. Department of Health and Human Services  
Food and Drug Administration  
Center for Drug Evaluation and Research (CDER)**

**January 2020  
Clinical Pharmacology**

## EMA: Final 2013

  
**EUROPEAN MEDICINES AGENCY**  
SCIENCE · MEDICINES · HEALTH

21 June 2012  
CPMP/EWP/560/95/Rev. 1 Corr. 2\*\*  
Committee for Human Medicinal Products (CHMP)

**Guideline on the investigation of drug interactions**

Discussion in the Efficacy Working Party (EWP)	June/October 1996 February 1997
Transmission to the CPMP	March 1997
Transmission to interested parties	March 1997
Deadline for comments	September 1997
Re-submission to the EWP	December 1997
Approval by the CPMP	December 1997
Date for coming into operation	June 1998
Draft Rev. 1 Agreed by the EWP	April 2010
Adoption Rev. 1 by CHMP for release for consultation	22 April 2010
End of consultation Rev. 1 (deadline for comments)	31 October 2010
Agreed by Pharmacokinetics Working Party	February 2012
Adopted by CHMP	21 June 2012
Date for coming into effect	1 January 2013

This guideline replaces guideline CPMP/EWP/560/95.

**Keywords** *Interaction, guideline, metabolism, inhibition, induction, transport, enzyme, transport protein, transporter, absorption, food, distribution, PBPK, herbal, SmPC*

\* The correction concerns section 3.3.4.3 (p. 28) and the corresponding decision tree no. 6 (p. 61) to read "if the observed Ki value is lower or equal to  $f \cdot f_i$ "; Appendix VII, Table 5 to read "See section 5.4.2." Decision tree 4.

30 Churchill Place • Canary Wharf • London E14 5EU • United Kingdom  
Telephone +44 (0)20 3660 6000 Facsimile +44 (0)20 3660 5555  
Send a question via our website [www.ema.europa.eu/contact](http://www.ema.europa.eu/contact)  
An Agency of the European Union  
© European Medicines Agency, 2015. Reproduction is authorised provided the source is acknowledged.

## PMDA: Final 2019

事務連絡  
平成 31 年 2 月 8 日

各都道府県衛生主管部（局）薬務主管課 御中

厚生労働省医薬・生活衛生局医薬品審査管理課

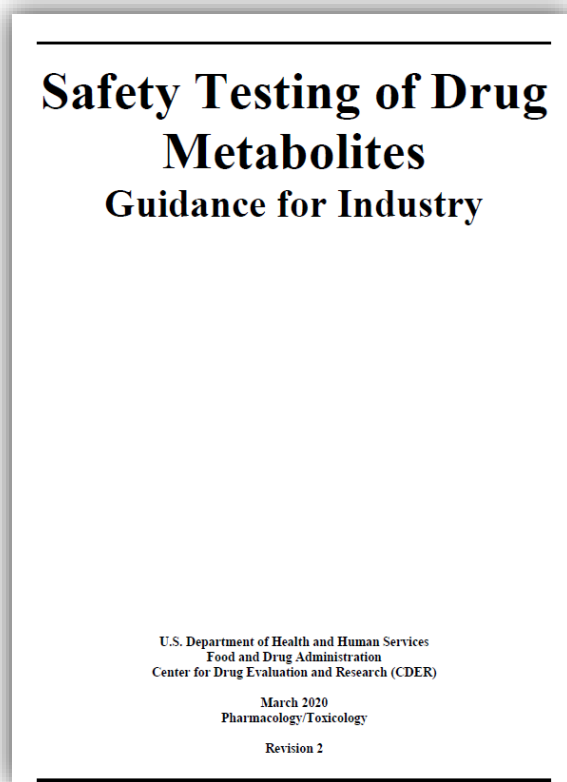
「医薬品開発と適正な情報提供のための薬物相互作用ガイドライン」等の英文版の送付について

標記について、別添1及び2のとおり取りまとめましたので、貴管下関係業者に対して周知方お願いします。

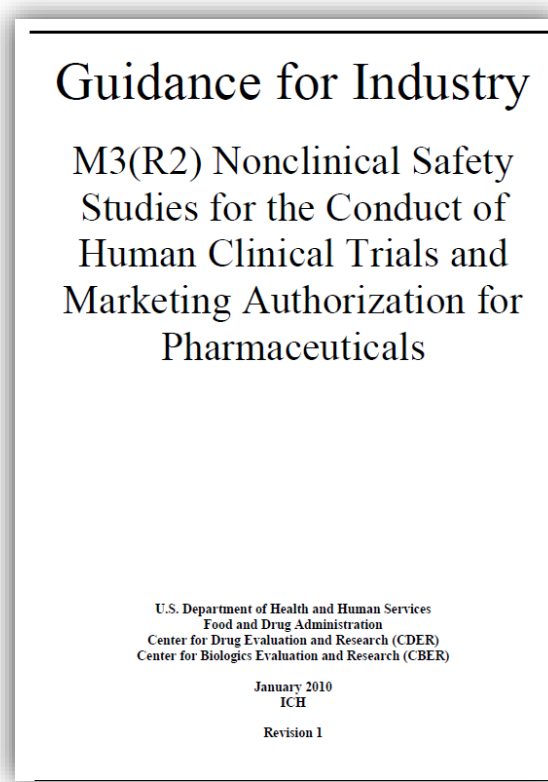
別添1 Guideline on drug interaction for drug development and appropriate provision of information  
別添2 Question and Answer for the "Guideline on drug interaction for drug development and appropriate provision of information"

# Additional Guidance

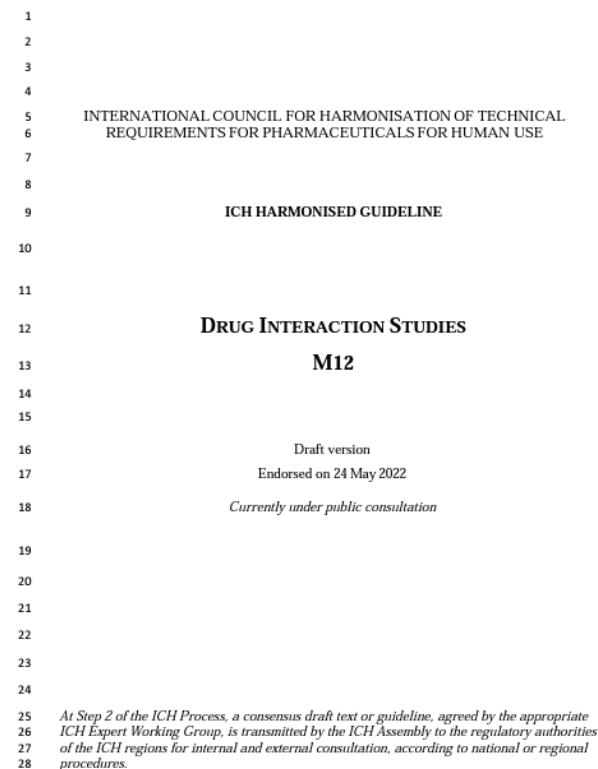
FDA “MIST”: Rev 2  
March 2020



FDA / ICH M3: Final 2010



ICH M12: Draft



# In vitro ADME & DDI study types

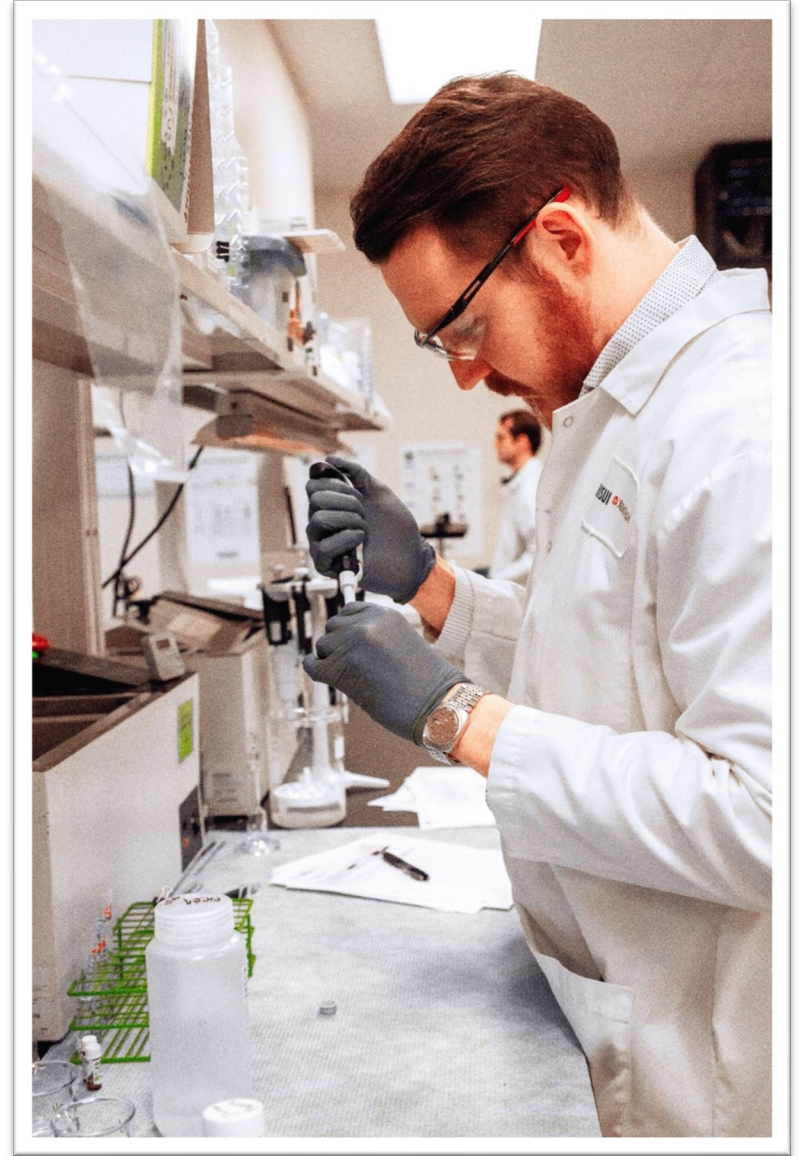
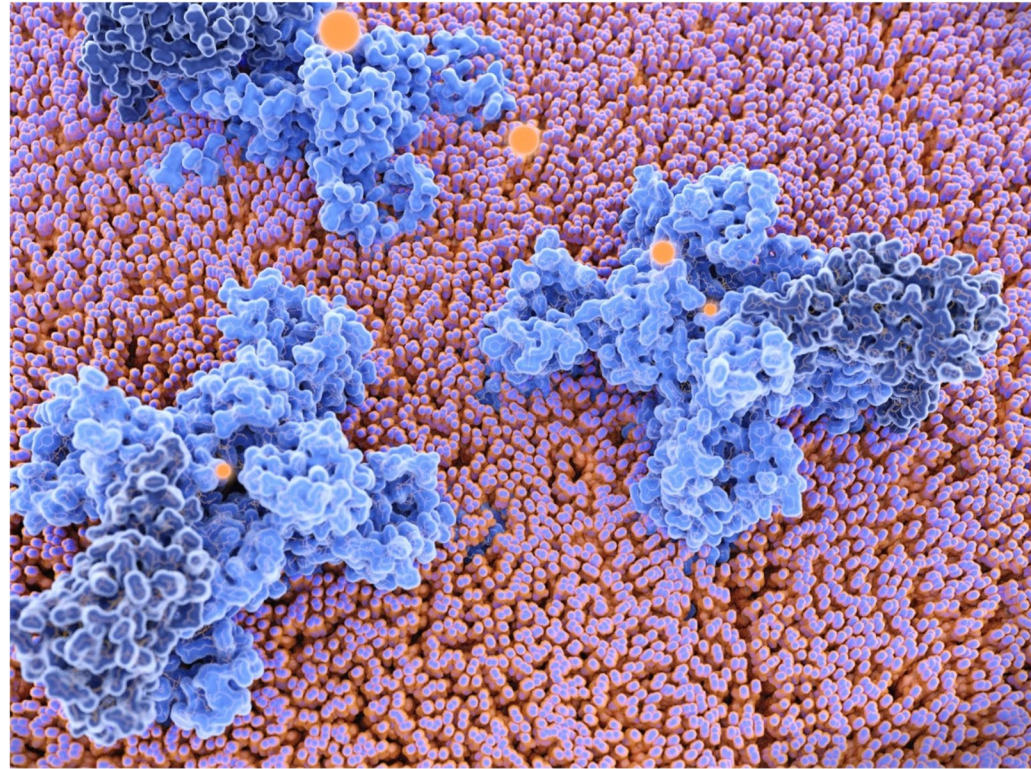
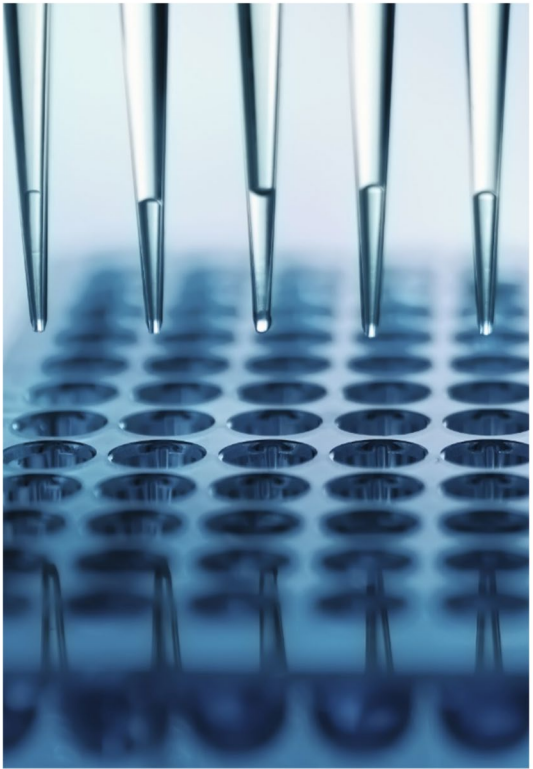
ADME component	Type of in vitro study
Drug Metabolism (M, E)	1. Inter-species comparative metabolism – $C_{\text{lint}}$ & $t_{1/2}$
	2. Metabolite ID – Qualitative analysis of metabolite profile
	3. Reaction phenotyping – Determine which CYPs are metabolizing
Drug Metabolizing Enzymes (M, E)	1. CYP Inhibition – Profile specific CYP inhibitions
	2. CYP induction – Induction potential for specific CYPs
Drug Transporters (A, D, E)	1. Transporter substrate – Determine Transporter substrate profile
	2. Transporter inhibition – Profile specific inhibition of major Transporters

# Test Systems for in vitro DDI Experiments

- **Subcellular fractions** – non-living, cell/tissue lysates that have been fractionated to enrich for certain enzyme activities.
  - **S9**
  - **Microsomes**
  - **Cytosol**
  - These subcellular fractions are derived from many relevant drug metabolizing organs/tissues. May require co-factors.
- **Cryopreserved primary hepatocytes** – living, isolated directly from living liver tissue and frozen for use at a later time. From multiple small animal models and human livers.

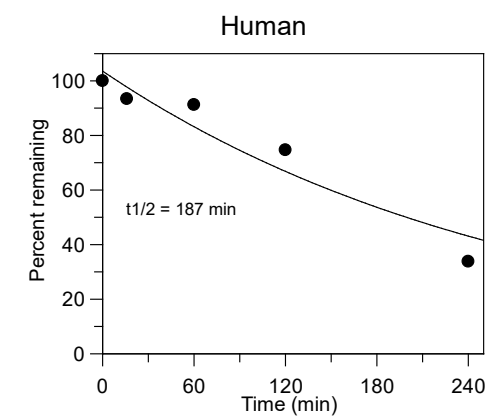
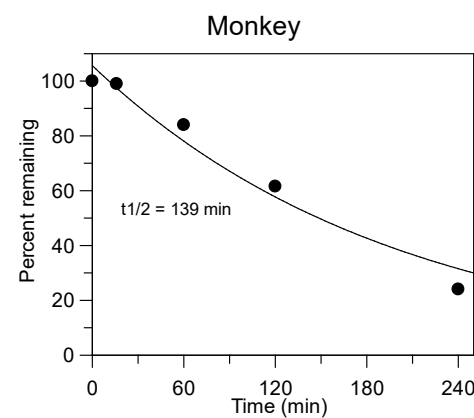
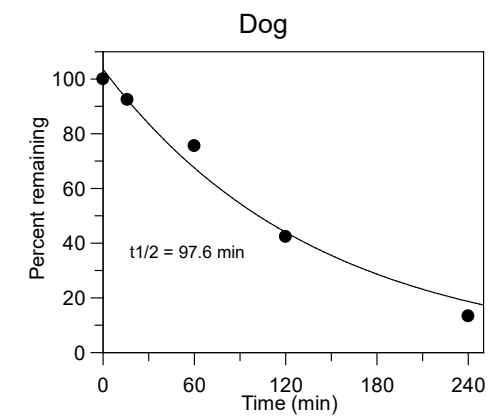
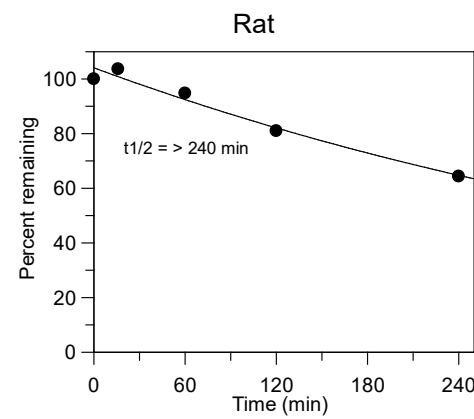
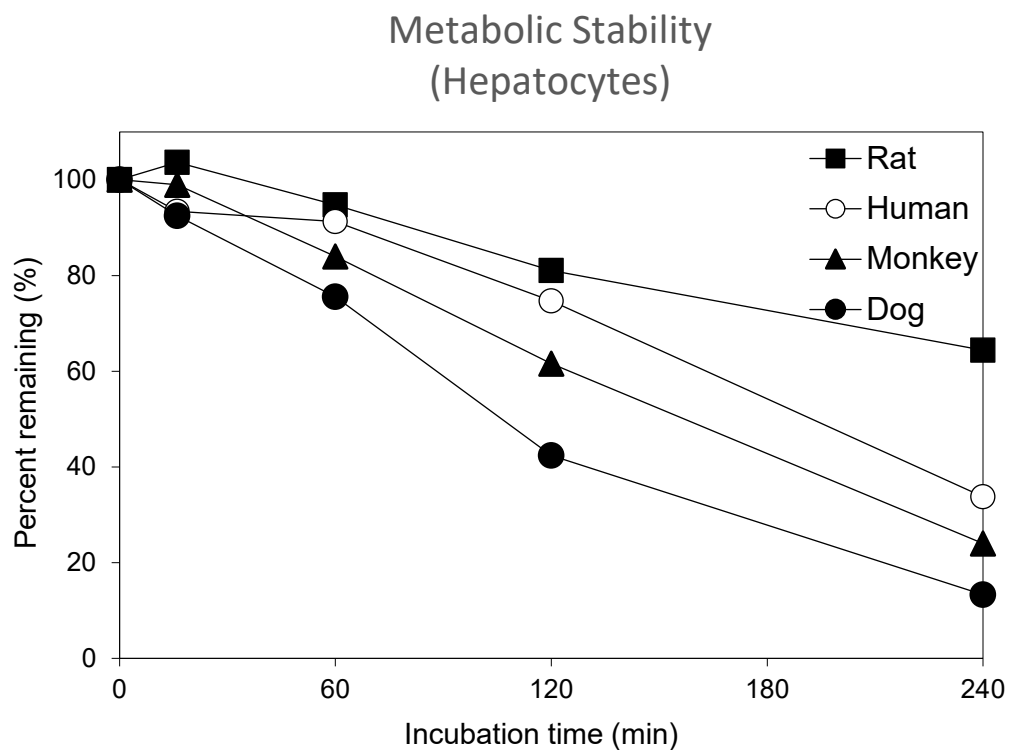


# Drug Metabolism Studies



# Drug Metabolism: Inter-Species Comparative Metabolism

- Design: Drug incubations with hepatocytes or subcellular fractions from various species
- **Typical species:** Human, Rat, Mouse, Dog, Rabbit, Monkey, Pig

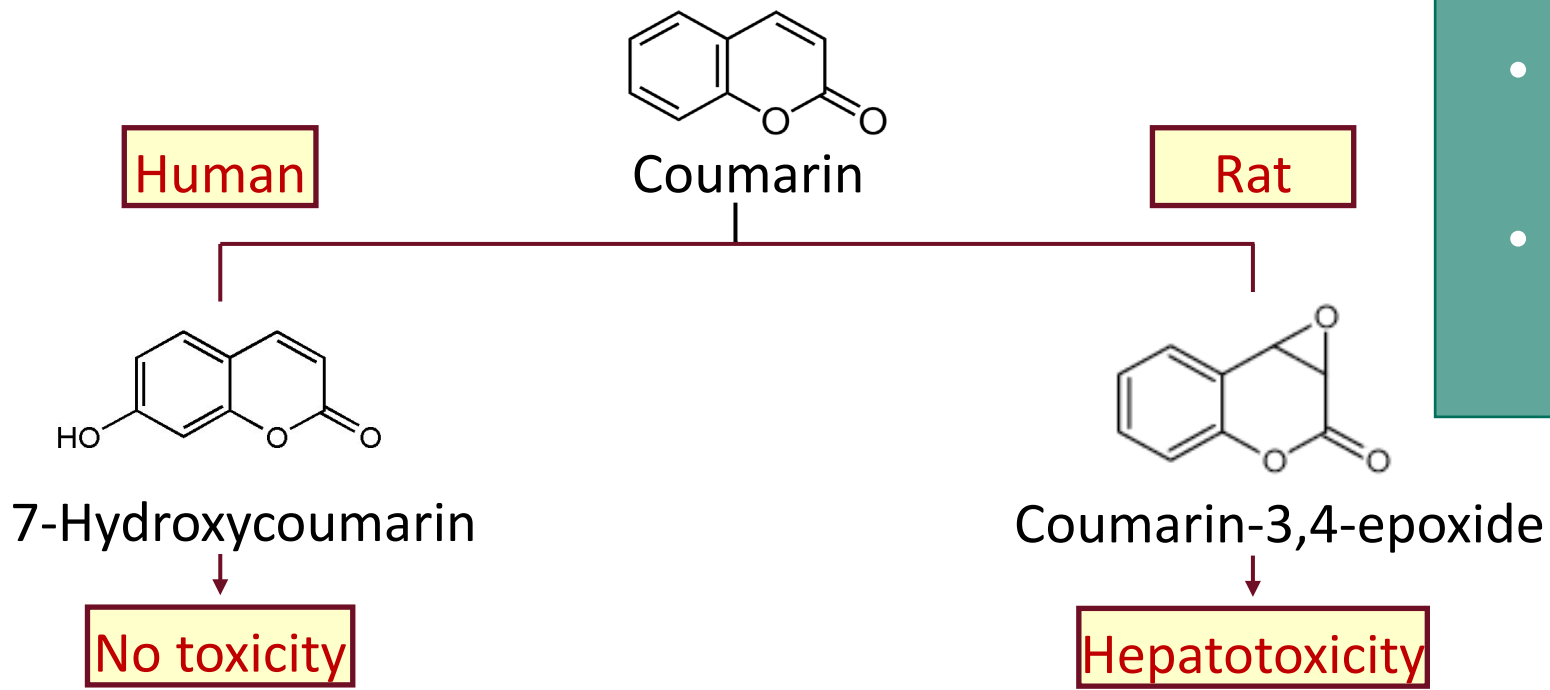




# Drug Metabolism: Inter-Species Comparative MetID

## Goals:

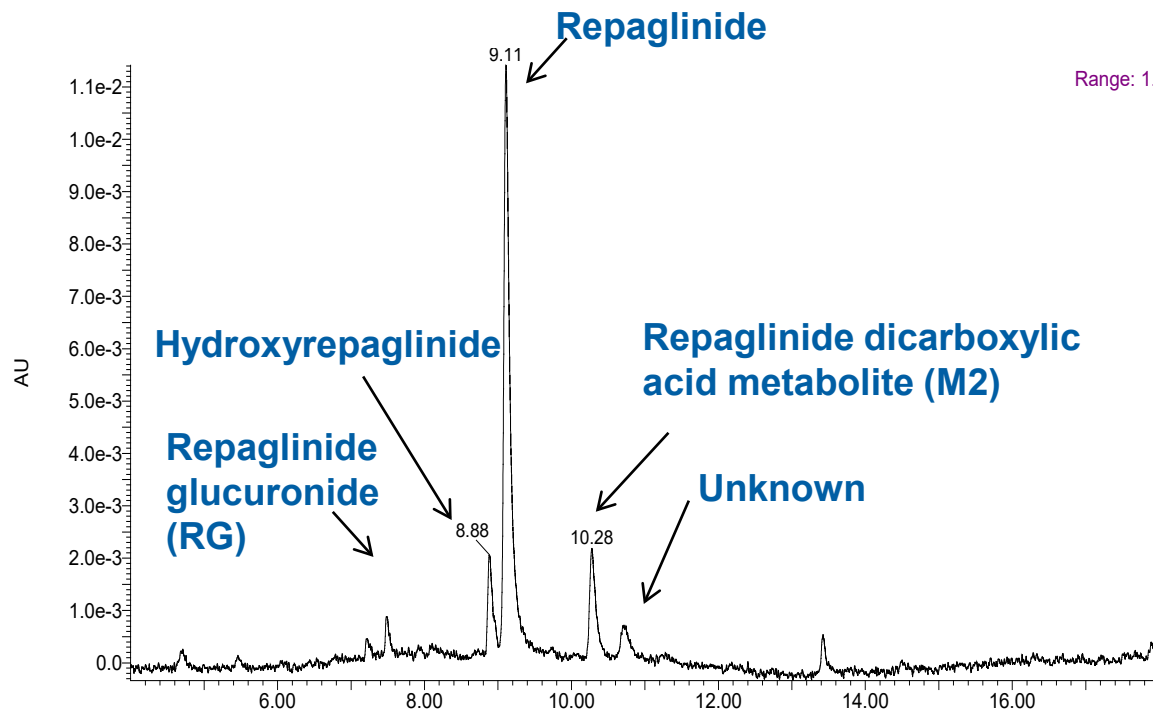
- Complete profile of metabolites
- Are there human specific or disproportionate metabolites?
- Which other species have a similar metabolic profile?



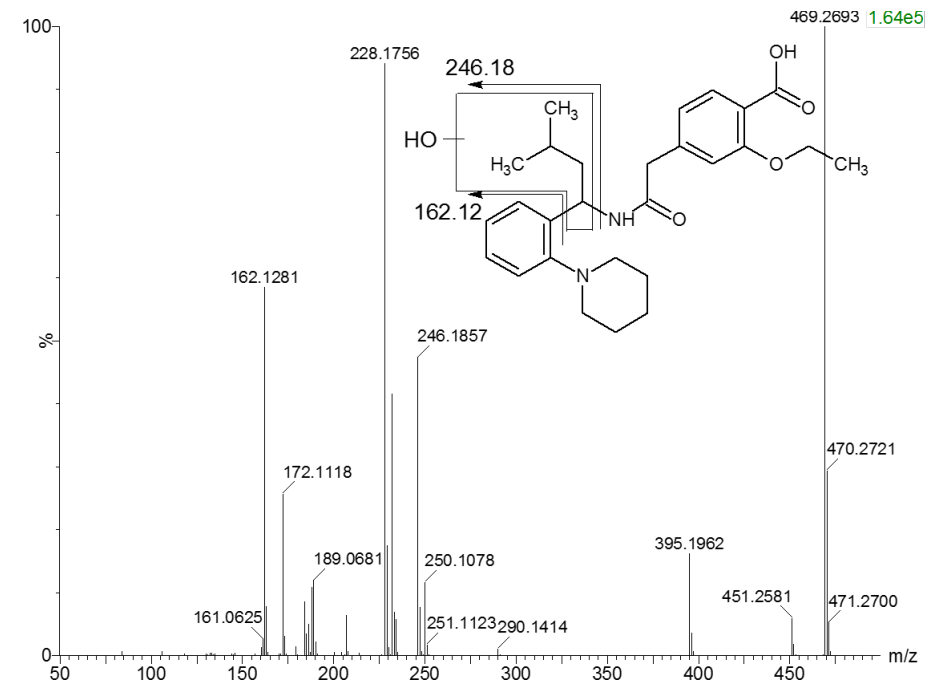
# Metabolite Characterization & ID

LC-MS/MS analysis – Qualitative identification of the metabolites

50  $\mu$ M Repaglinide; Human hepatocytes; 60 minutes; 37°C



Hydroxyrepaglinide LC-MS/MS





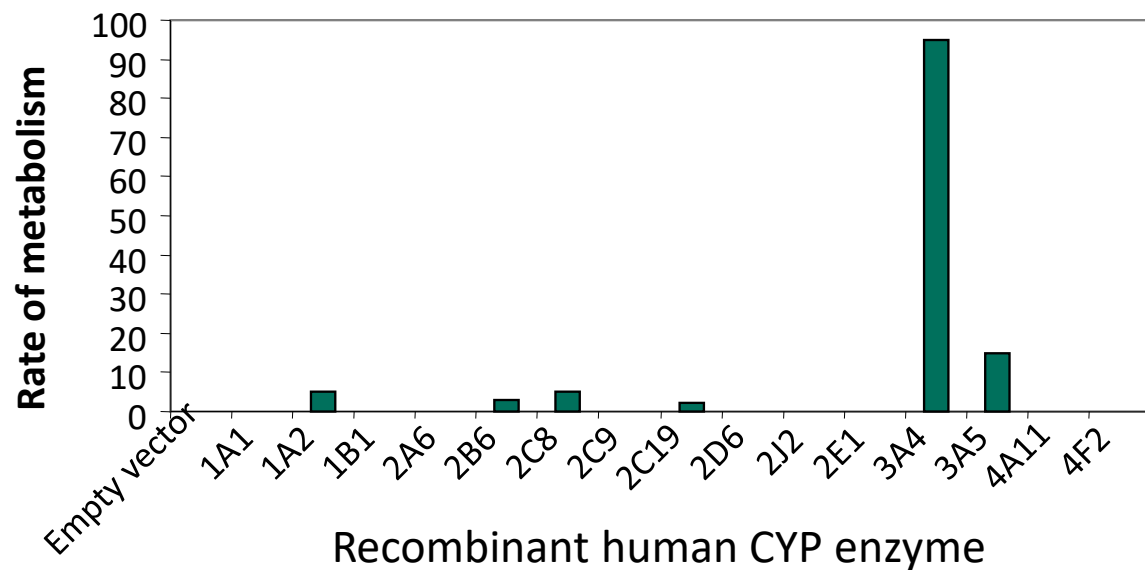
# Cross-Species Met ID

Component	Retention time (min)	Mass shift	Proposed biotransformation	Mouse	Rat	Dog	Pig	Human
C1	3.43	255.9889	Sulfation + glucuronidation	+	+	+	+	+
C2	3.63	354.0783	Di-glucuronidation + hydrogenation	+	+	+	+	+
C3	3.78	159.9135	Di-sulfation	+	+	+	+	+
C4	4.00	258.0045	Sulfation + glucuronidation + hydrogenation	+	+	+	+	+
C5	4.41	161.9298	Di-sulfation + hydrogenation	+	+	+	+	+
C6	4.44	194.0428	Glucuronidation + oxygenation + hydrogenation	ND	ND	ND	+	+

# Drug Metabolism: CYP Reaction Phenotyping (Victim potential)

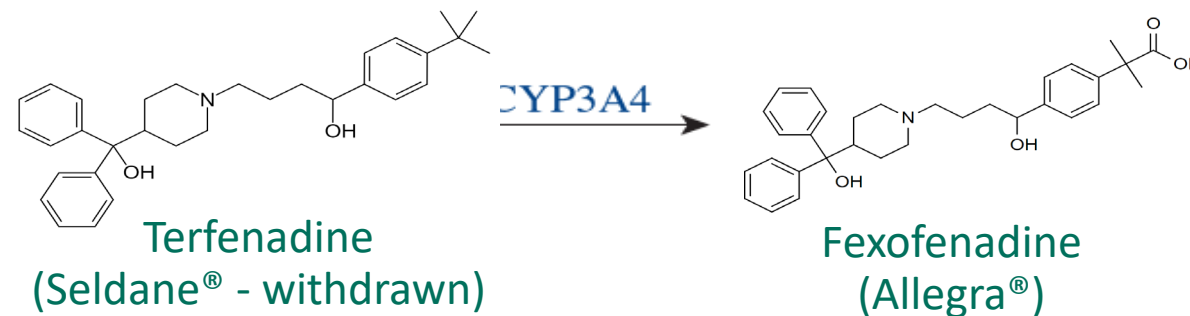
- Design: Incubate drug + recombinant human CYPs or human liver microsomes or hepatocytes ± selective inhibitors
- Goal: Determine which CYPs drive the metabolism of the drug
- Unique CYP metabolism is of concern

High DDI potential: few enzymes involved; Avoiding DDI, multiple pathways are preferred



Follow-up studies:

- Confirm with selective inhibitors
- Evaluate non-CYP pathways in HLM or hepatocytes



# Drug Metabolizing Enzymes (Perpetrator Potential)

Enzyme Induction



Enzyme Inhibition

# Drug Metabolizing Enzymes: CYP Inhibition

## Design:

Drug incubations with  
HLM + marker substrate  $\pm$   
pre-incubation

## Goal:

Predict clinically  
relevant inhibition of  
CYP enzymes

## **CYP**

CYP1A2

CYP2B6

CYP2C8

CYP2C9

CYP2C19

CYP2D6

CYP3A4

CYP3A4

## **Activity Assay**

Phenacetin *O*-dealkylation

Bupropion hydroxylation

Amodiaquine *N*-dealkylation

Diclofenac 4'-hydroxylation

*S*-Mephenytoin 4'-hydroxylation

Dextromethorphan *O*-dealkylation

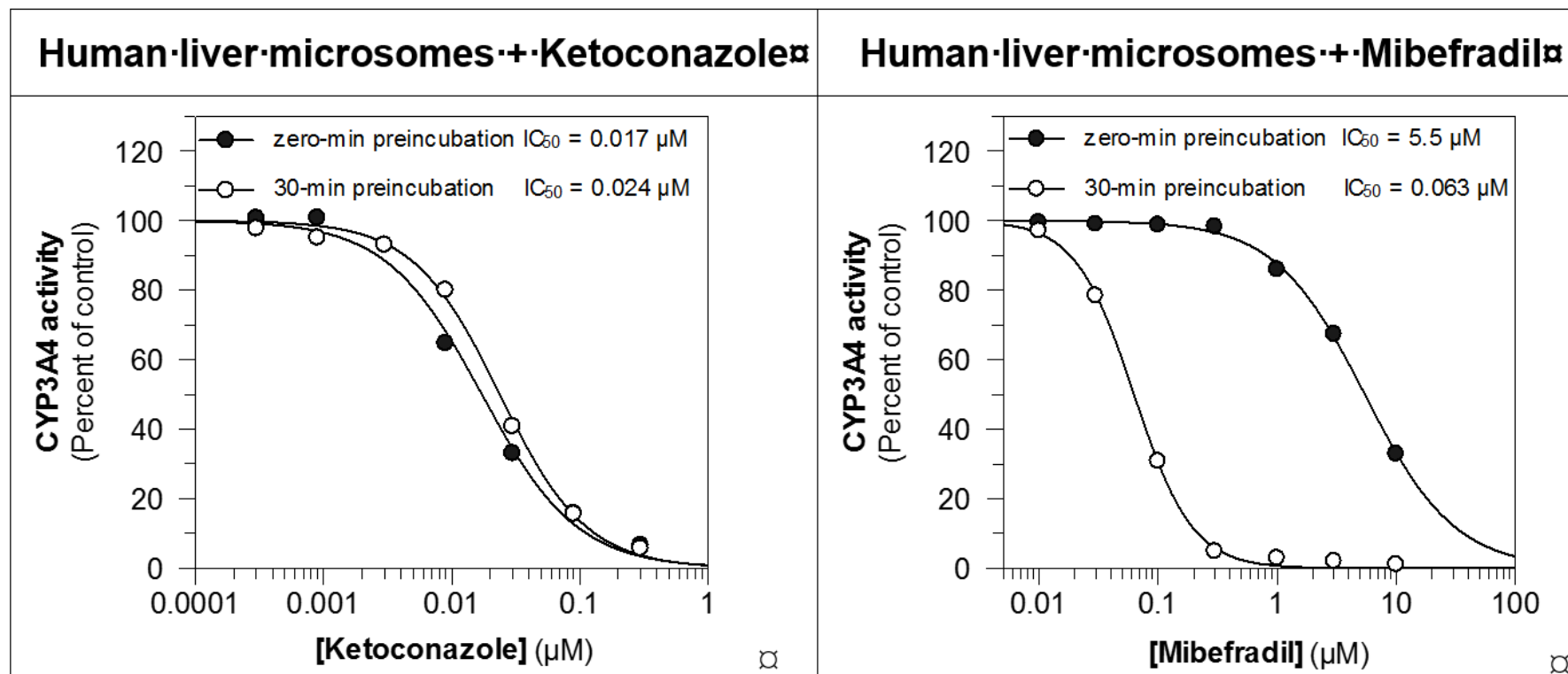
Testosterone 6 $\beta$ -hydroxylation

Midazolam 1'-hydroxylation



# Drug Metabolizing Enzymes: CYP Inhibition

- Design: Drug incubations with HLM + marker substrate  $\pm$  pre-incubation
- Goal: Assess inhibition of CYP enzymes



## Ketoconazole:

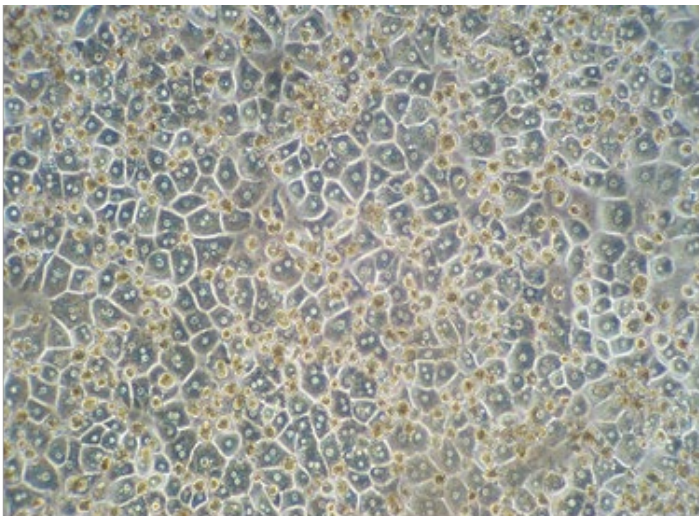
Potent inhibitor of 3A4  
precludes  
coadministration of  
other drugs

## Mibefradil

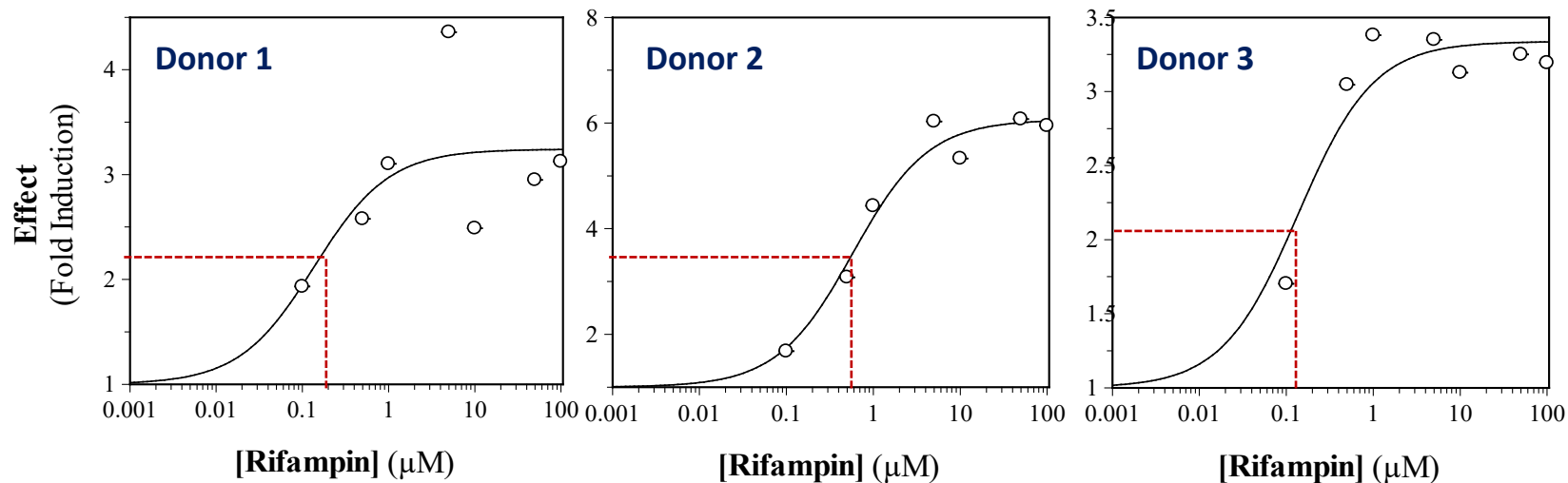
Removed from market  
in 1998 due to  
potential for fatal DDIs  
TDI has greater DDI  
risk

# Drug Metabolizing Enzymes: CYP Induction

- Design: Drug incubations in cultured human hepatocytes, measure mRNA (or activity) of various CYPs
- Goal: Assess induction of CYP enzymes
- For both inhibition and induction studies, basic or mechanistic models are further utilized to determine the clinical DDI potential of a drug and requirement of a clinical DDI study.



Cultured Human Hepatocytes



Parameter	Value	Std. Error
E <sub>max</sub>	3.2417	0.3163
EC <sub>50</sub>	0.1371	0.1427

Parameter	Value	Std. Error
E <sub>max</sub>	6.0649	0.1981
EC <sub>50</sub>	0.5795	0.1244

Parameter	Value	Std. Error
E <sub>max</sub>	3.3378	0.1200
EC <sub>50</sub>	0.1362	0.0528

# Drug Transporters



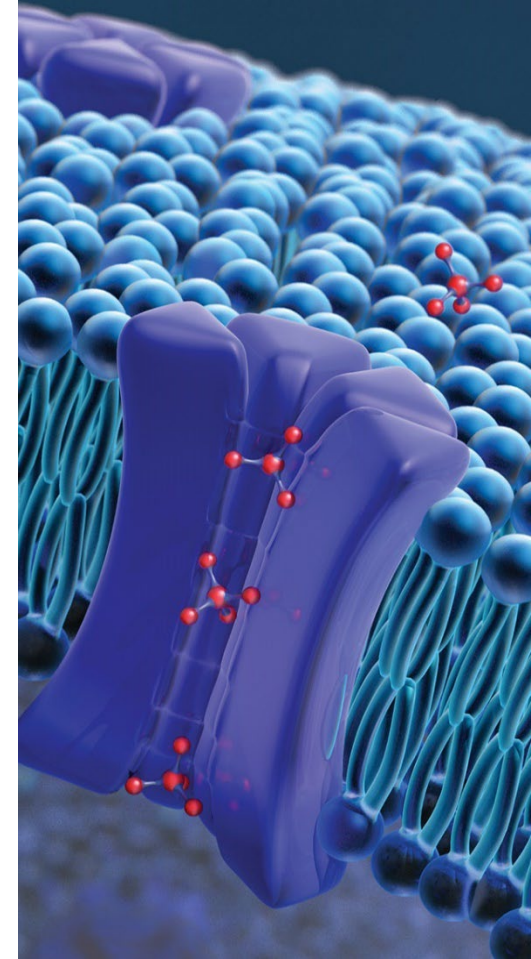
**Substrate Potential  
(Victim)**

**Transporter Inhibition  
(Perpetrator)**

# “Why conduct drug transporter DDI studies?”

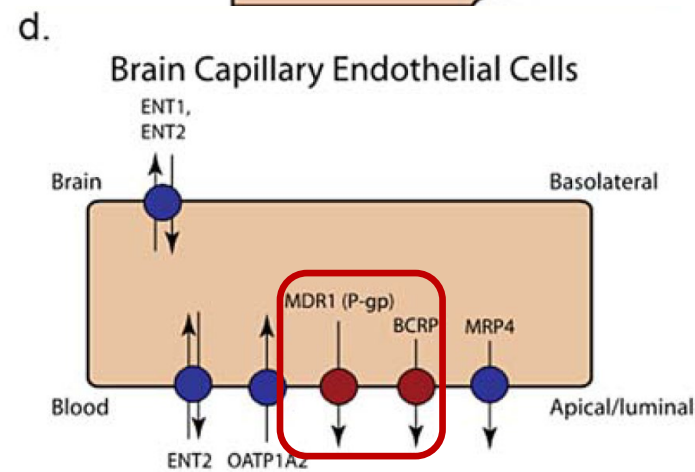
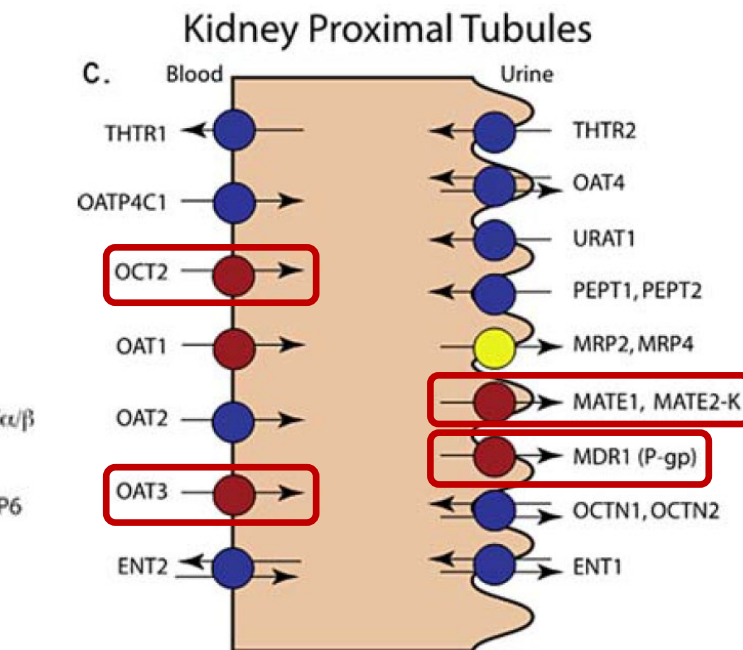
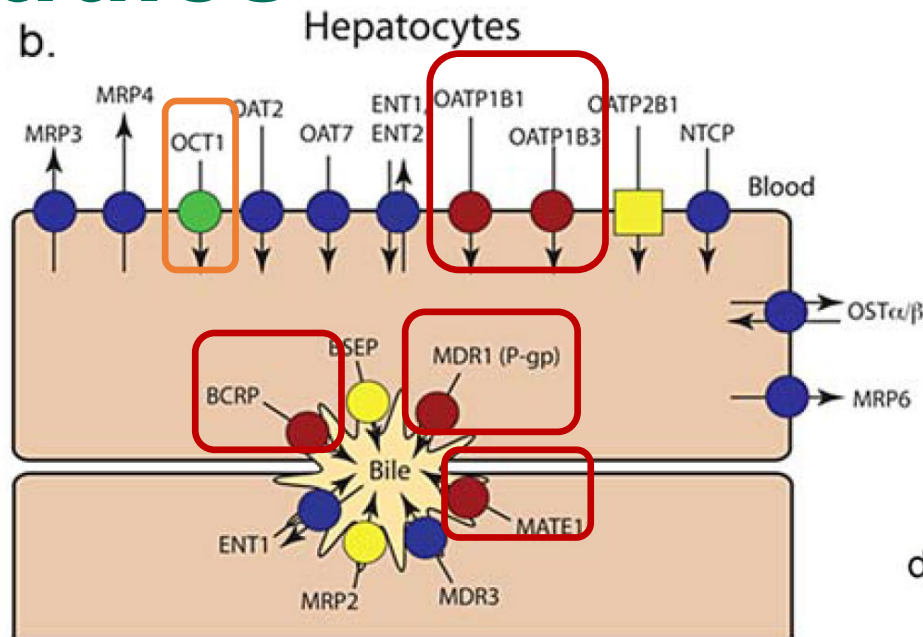
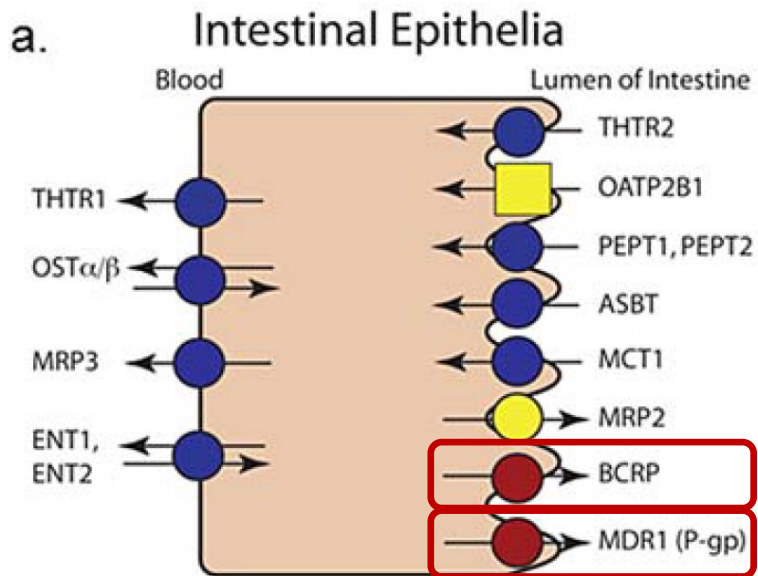
## ADME

- Drug absorption, distribution, tissue-specific drug targeting, and elimination
- Drug-drug interactions
  - Clearance of transporter substrates (Victims) can be impacted by transporter inhibitors or inducers (Perpetrators)
  - Toxicity or loss of efficacy
- Real world example - Statins
  - Hepatic uptake transporter (OATPs) substrates: taken up in the liver, reduce cholesterol
  - Cyclosporine inhibits OATPs: up to 10-fold increase in statin exposure
  - Toxic side effect: rhabdomyolysis (skeletal muscles break down, cells released into bloodstream, can lead to kidney failure and possibly death)





# Transporter studies



Inhibition for all in **red** (FDA & PMDA); **orange** (EMA)

Substrate potential:

P-gp and BCRP (all **orally administered** drugs)

Hepatic uptake: If hepatic metabolism or biliary secretion  $\geq 25\%$

Renal: If active renal secretion  $\geq 25\%$  of total clearance

Figures from Zamek-Gliszczyński et al. ITC3 (2018) *CPT* 104:890-899

# Importance of proper study design & data interpretation for in vitro DDI studies



# Importance of study design: CYP inhibition study

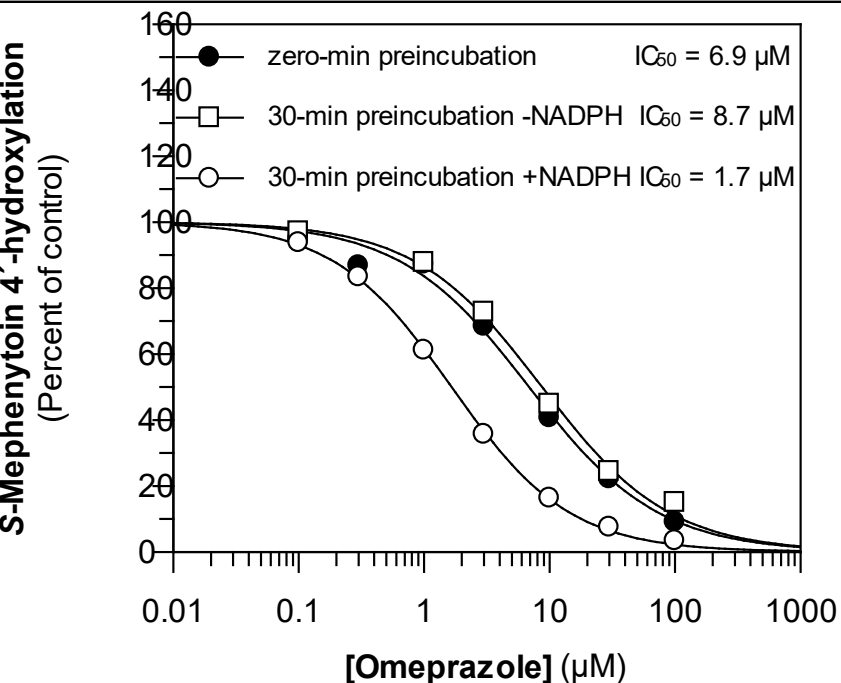
False negative results arise from poorly designed studies

- Example: Clinically relevant time-dependent inhibition of CYP2C19 by omeprazole missed with high [protein] and long marker substrate incubation

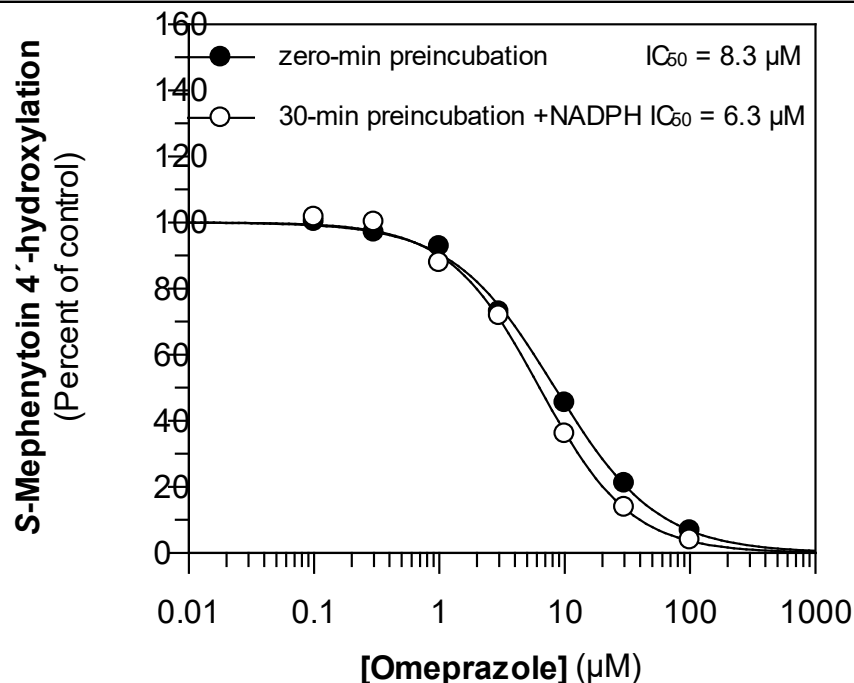
**Detected**

**Missed**

0.1 mg/mL protein, 5 min substrate incubation



1 mg/mL protein, 30 min substrate incubation

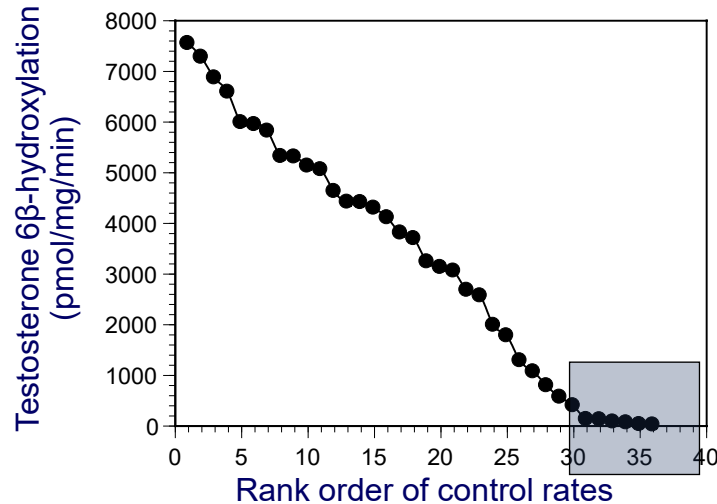


Time-dependent inhibition of CYP2C19 by omeprazole is readily detectable with HLM at 0.1 mg/mL with a 5-min substrate incubation period (**left**) but not at 1.0 mg/mL with a 30-min incubation period (**right**)

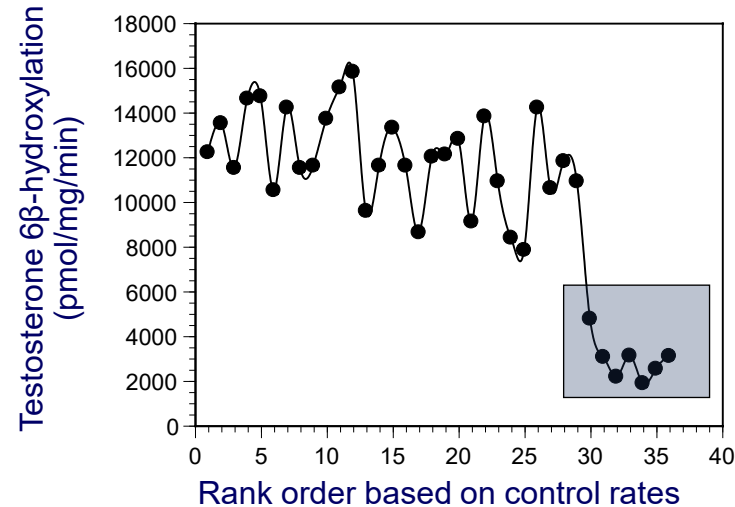
# Proper test system selection: CYP Induction study

CYP induction studies: positive controls with very large induction

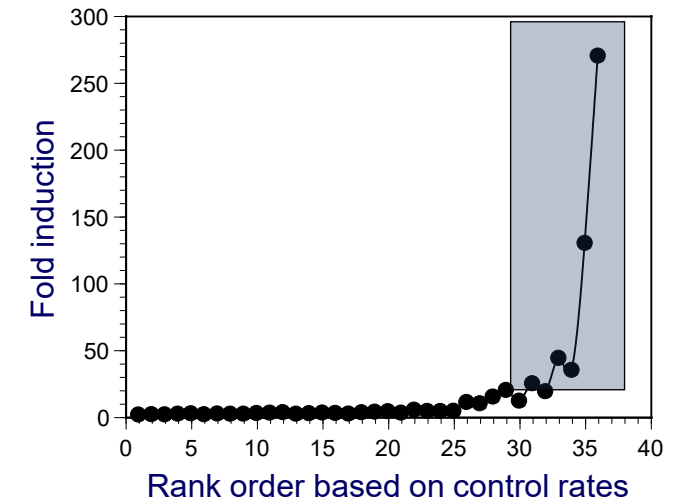
Control hepatocytes  
CYP3A4 activity



Rifampin-treated (20  $\mu$ M) hepatocytes  
CYP3A4 activity



Fold induction



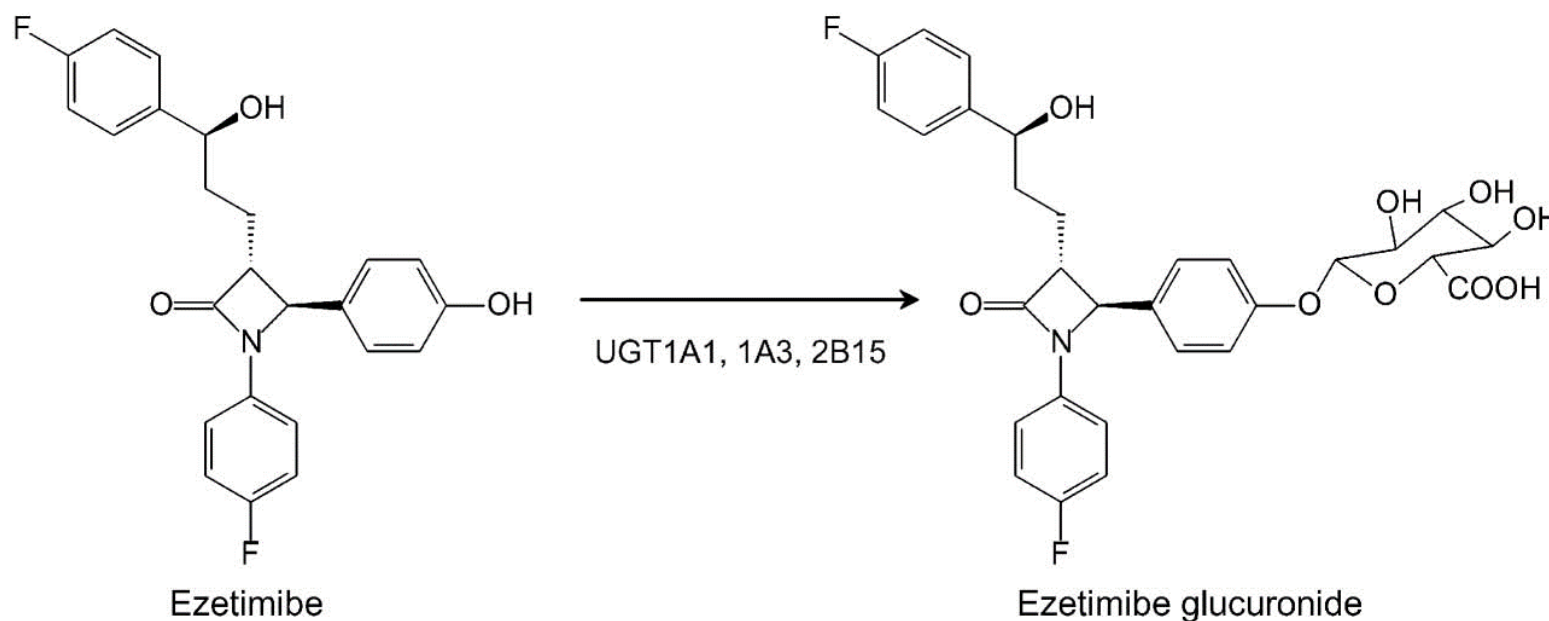
- When it comes to induction, more is not always better
- A high fold-induction (>20 fold) of CYP3A4 activity by rifampin is a sign of incomplete hepatocellular differentiation of the cultured human hepatocytes



# Proper test system selection : Reaction phenotyping

Metabolism studies: Choose the right test system based on the structure

- Ezetimibe is oxidized by CYP3A4 however results with HLM & NADPH alone can be misleading.



Oxidation does not occur clinically due to rapid phenolic glucuronidation. Recombinant human UGTs or human hepatocytes would be a better test system. CYPs are not the only enzyme system.

# Timing of ADME studies



# Drug Development Pipeline: Timing *in vitro* DDI studies

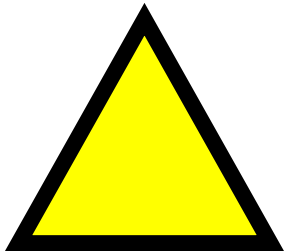
Priority depends on strategy for each drug & need for de-risking at each stage

Type of drug	Lead optimization	Pre-IND	Phase I to NDA
Typical small molecule	<ol style="list-style-type: none"> <li>1. Comparative metabolism</li> <li>2. Metabolite ID</li> <li>3. Screening for others</li> </ol>	<ol style="list-style-type: none"> <li>1. CYP inhibition</li> <li>2. CYP Induction</li> <li>3. Transporter inhibition</li> <li>4. Limited transporter substrate</li> </ol>	<ol style="list-style-type: none"> <li>1. Reaction phenotyping</li> <li>2. Additional transporter substrate (dependent on routes of elimination)</li> </ol>
Small molecule with orphan, breakthrough status, etc.	<ol style="list-style-type: none"> <li>1. Comparative metabolism</li> </ol>	May be able to defer	<ol style="list-style-type: none"> <li>1. Metabolite ID</li> <li>2. CYP inhibition</li> <li>3. Transporter inhibition</li> <li>4. Reaction phenotyping</li> <li>5. CYP induction</li> </ol>
Peptides, oligos, ADCs, other biologics	May be able to defer	May be able to defer	<ol style="list-style-type: none"> <li>1. Metabolite ID</li> <li>2. CYP inhibition</li> <li>3. Transporter inhibition</li> <li>4. Reaction phenotyping</li> <li>5. CYP induction</li> </ol>

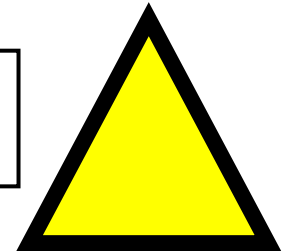
# “Why conduct these studies? Is this just box checking?”

**No.** The information in aggregate has real utility:

1. Provide deeper understanding of the molecule
  - Metabolism, enzymes involved in metabolism, etc.
  - The information generated from DDI studies goes on the drug label
  - From the pharma company’s perspective these studies can inform go/no-go decisions for a drug candidate
  - Predictive toxicology and dose selection for certain non-clinical *in vivo* studies
2. Prepare for clinical studies
  - Prediction of FIH dose and DDI risk
3. Satisfy regulatory expectations and comply with regulatory guidance

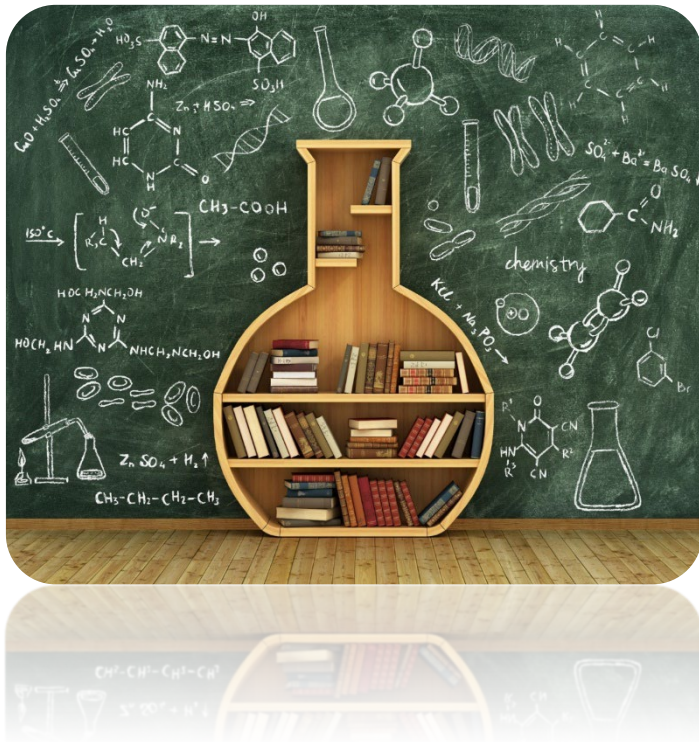


**These studies may appear deceptively simple – maximizing insight can be complex**



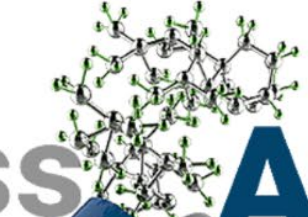


# For questions or further resources:



**XENOTECH**

[How We Help](#) [Contract Services](#) [Test Systems](#) [Scientific Resources](#) [About](#) [Contact Us](#)

access  **ADME**

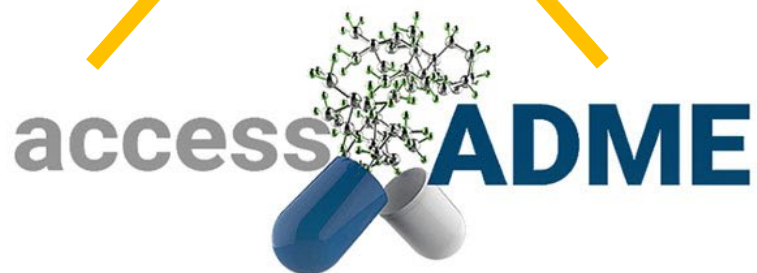


**How Can We Help?**

# Further Resources

<https://www.xenotech.com/access-adme-research-resources/>

- Videos
- Webinars
- Blogs
- And more!



Scientific Resources Related to  
ADME, DMPK in Drug  
Development



## Webinar Topic Request Form:

[www.xenotech.com/scientific-resources/upcoming-webinars](http://www.xenotech.com/scientific-resources/upcoming-webinars)

## XenoTech's Services & Products:

### **In Vitro ADME/PK & DDI**

- Drug Transport
- Drug Metabolism
- Enzyme Inhibition & Induction
- Protein Binding
- Metabolite Identification
- ADME Screening

### **In Vivo ADME/PK & Distribution**

- QWBA
- Microautoradiography
- Excretion / Mass Balance
- Tissue Distribution
- Blood / Plasma & Lymphatic Partition Rate

### **Bioanalytical**

### **Pharmacology**

- In Vitro Ligand Binding & Radioreceptor Assays
- Immunoassays

### **Chemical Synthesis**

- Radiolabeled Synthesis
- Metabolite Synthesis
- Peptide Synthesis

### **Consulting...**

### **Cellular Products**

- Hepatocytes (Cryo/Fresh, Genotyped...)
- Non-Parenchymal Cells (Kupffer Cells)

### **Subcellular Fractions**

- Liver Microsomes
- S9 Fractions
- Cytosol
- Homogenate
- Lysosomes & Tritosomes
- Mitochondria
- Extrahepatic Fractions

### **Custom Products**

- Various Species, Tissues & Preparations

### **Research Biobank**

- Normal & Diseased Tissue Samples

### **Recombinant Enzymes**

### **Substrates & Metabolites**

### **JCRB Cell Lines...**

