

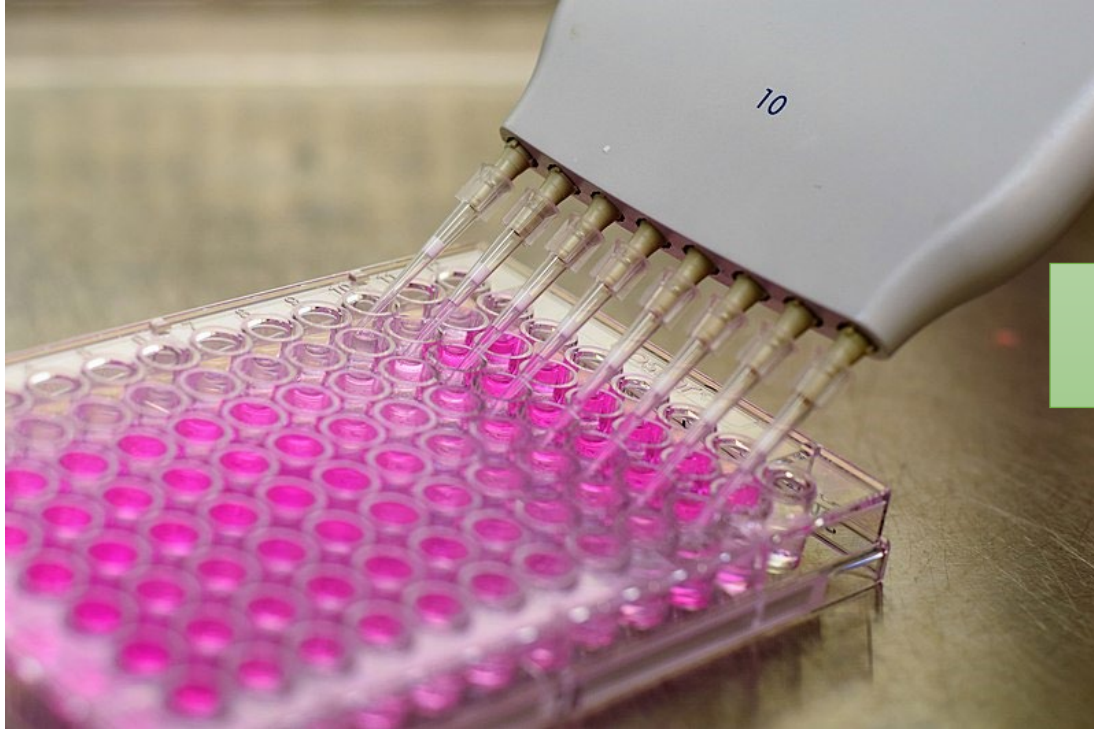
Model-Based Approaches to DDI Risk Prediction: Navigating the Transition from In Vitro Data to In Silico Modeling



Pallavi Limaye, Ph.D., DABT
Director, Scientific Consulting
XenoTech



In Vitro to In Vivo Extrapolation IVIVE



Not a stand alone study
A follow up analysis to routine in vitro ADME-DDI studies



FDA Guidance Verbiage - Background

“The evaluation of DDI potential often starts with **in vitro** experiments to identify potential factors influencing drug disposition to elucidate potential DDI mechanisms and to yield kinetic parameters for use in further studies. Results of in vitro experiments, **along with clinical pharmacokinetic (PK) data, provide mechanistic information** that can inform the need for and proper design of potential future clinical studies.”

— FDA 2020, “Drug–Drug Interaction Assessment for Therapeutic Proteins Guidance for Industry”

ADME & DDI study types for in vitro to in vivo prediction

Models focus on DDI via perpetrator potential of a drug

ADME component	Type of in vitro study to determine perpetrator potential
CYP450 Enzymes	<ol style="list-style-type: none">1. CYP Inhibition – Identification of specific CYP enzymes inhibited by the drug2. CYP induction – Induction potential for specific CYPs
Drug Transporters	<ol style="list-style-type: none">1. Transporter inhibition – Profile specific inhibition of major Transporters

Models for DDI Prediction

- Basic static models
- Mechanistic static models
- Dynamic mechanistic models (PBPK)



What we can provide:

- Basic static models
- Mechanistic static models
- Dynamic mechanistic models (PBPK)

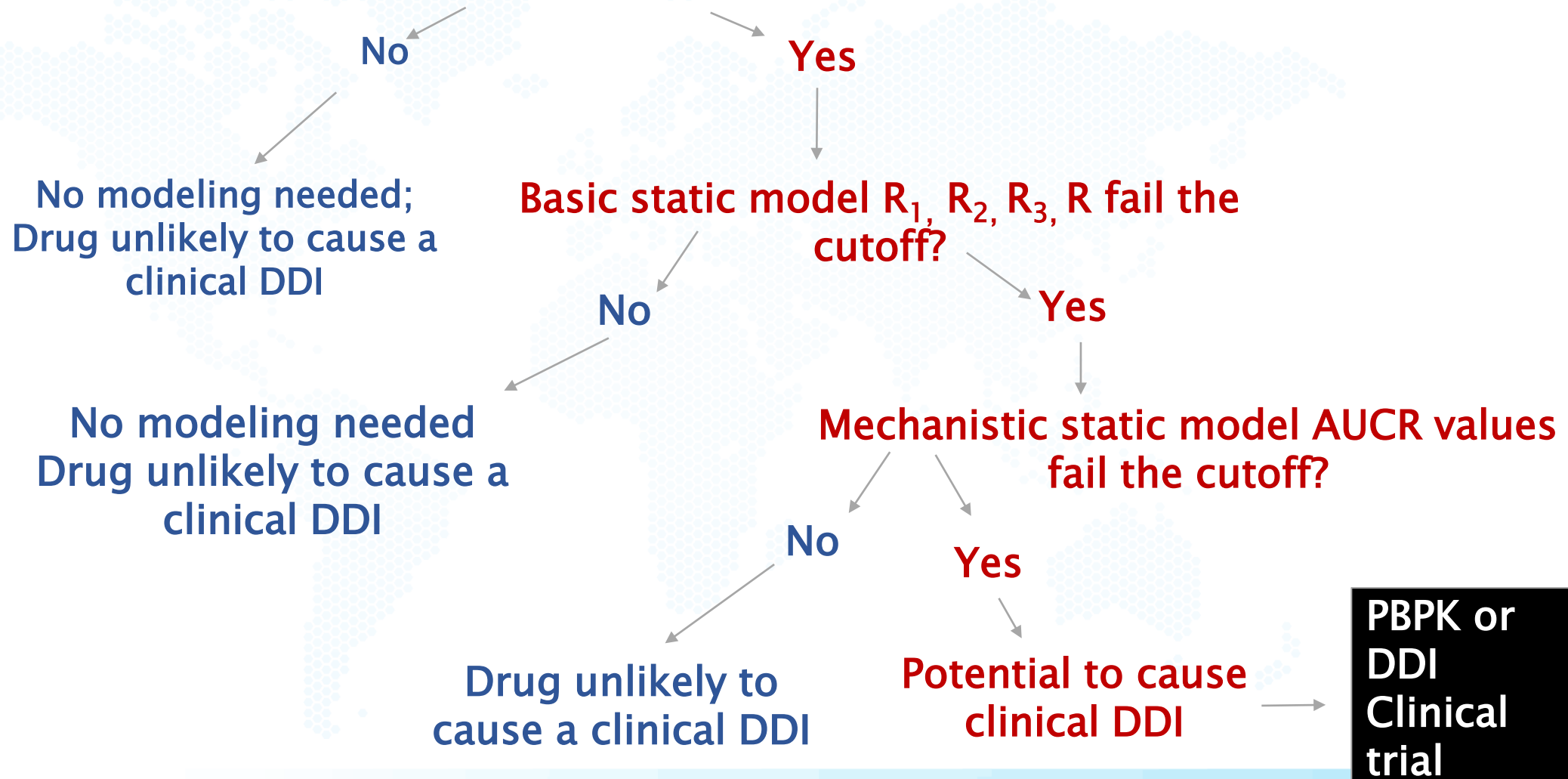


Predictive Models Based on Study Types

Study	Basic Static		Static Mechanistic
CYP inhibition	Hepatic:		AUCR
	Direct	TDI	
	R_1	R_2	
CYP induction	R_3 , Relative Induction Score (RIS)		
Transporter inhibition	R		-

In vitro CYP Inhibition/Induction or Transporter Inhibition?

Decision Tree



Basic Static Models: CYP Inhibition and Induction

Type of Data	Equation	Cut Off Values
Hepatic CYP reversible inhibition	$R_1 = 1 + (I_{\max,u} / K_{i,u})$	Potential to inhibit if $R_1 \geq 1.02$
Intestinal CYP inhibition	$R_{1,\text{gut}} = 1 + (I_{\text{gut}} / K_i)$	Potential to inhibit if $R_{1,\text{gut}} \geq 1.1$
Hepatic CYP irreversible/TDI	$R_2 = (K_{\text{obs}} + k_{\text{deg}}) / k_{\text{deg}}$	Potential to inhibit if $R_2 \geq 1.25$
CYP induction	$R_3 = 1 / [1 + d \times ((E_{\max} \times 10 \times I_{\max,u}) / (EC_{50} + 10 \times I_{\max,u}))]$ RIS = 1. $E_{\max} \times I_{\max,u} / EC_{50} + I_{\max,u}$ 2. $I_{\max,u} / EC_{50}$ values	Potential to induce if $R_3 \leq 0.8$ Potential to induce based on AUC decrease of victim drug depending on RIS

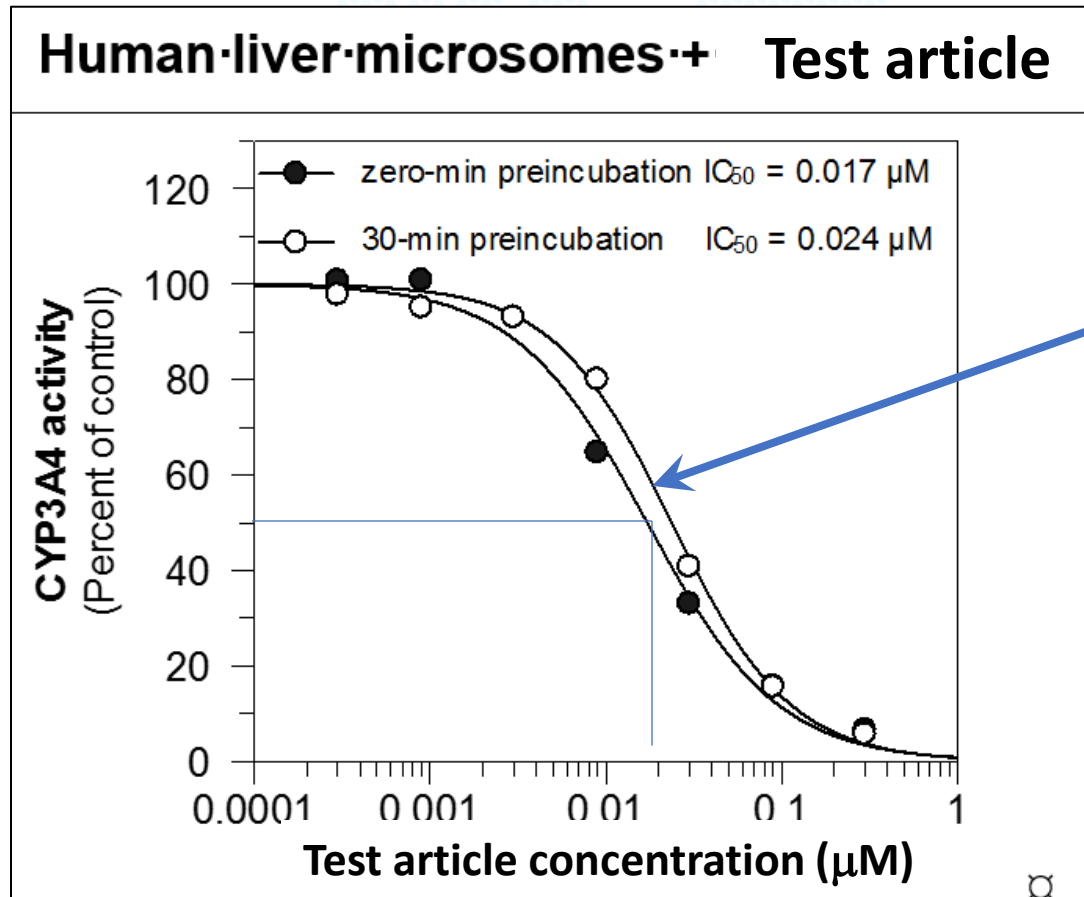
Basic Models: Transporter Inhibition

Transporter	Ratio	Potential to inhibit if
P-gp	I_{gut} / IC_{50}	Ratio ≥ 10
BCRP	I_{gut} / IC_{50}	Ratio ≥ 10
OATP1B1	$1 + (I_{\text{in,max,u}}) / IC_{50}$	Ratio ≥ 1.1
OATP1B3	$1 + (I_{\text{in,max,u}}) / IC_{50}$	Ratio ≥ 1.1
OAT1	$I_{\text{max,u}} / IC_{50}$	Ratio ≥ 0.1
OAT3	$I_{\text{max,u}} / IC_{50}$	Ratio ≥ 0.1
OCT2	$I_{\text{max,u}} / IC_{50}$	Ratio ≥ 0.1
MATE1	$I_{\text{max,u}} / IC_{50}$	Ratio ≥ 0.1
MATE2-K	$I_{\text{max,u}} / IC_{50}$	Ratio ≥ 0.1

Predictive Models Based on Study Types

Study	Basic Static			Static Mechanistic
CYP inhibition	Hepatic:		Intestinal	AUCR
	Direct	TDI	Direct	
	R_1	R_2	$R_{1 \text{ gut}}$	
CYP induction	R_3 , Relative Induction Score (RIS)			
Transporter inhibition	R			-

CYP Direct Inhibition – In Vitro Data



Superimposable curves
indication direct or reversible
inhibition

Kinetic Constants

$IC_{50} (\mu M)$	0.017
$K_i (\mu M) = IC_{50}/2$	0.0085

Study shows that the drug is a
CYP3A4 direct inhibitor

Follow up Prediction Model: R_1 Value Determination

Hepatic CYP Direct inhibition

$$R_1 = 1 + (I_{\max,u} / K_{i,u})$$

Potential to inhibit if $R_1 \geq 1.02$

Parameters required

- $I_{\max,u}$ – Provided by sponsor ($I_{\max,u}$ is unbound C_{\max})
- $K_{i,u}$ – In vitro inhibition study data



R_1 Calculation for Hepatic Reversible Inhibition

Where $I_{\max,u} = 0.025 \mu\text{M}$

$K_{i,u} = 0.0084 \mu\text{M}$

$$R_1 = 1 + (I_{\max,u} / K_{i,u})$$

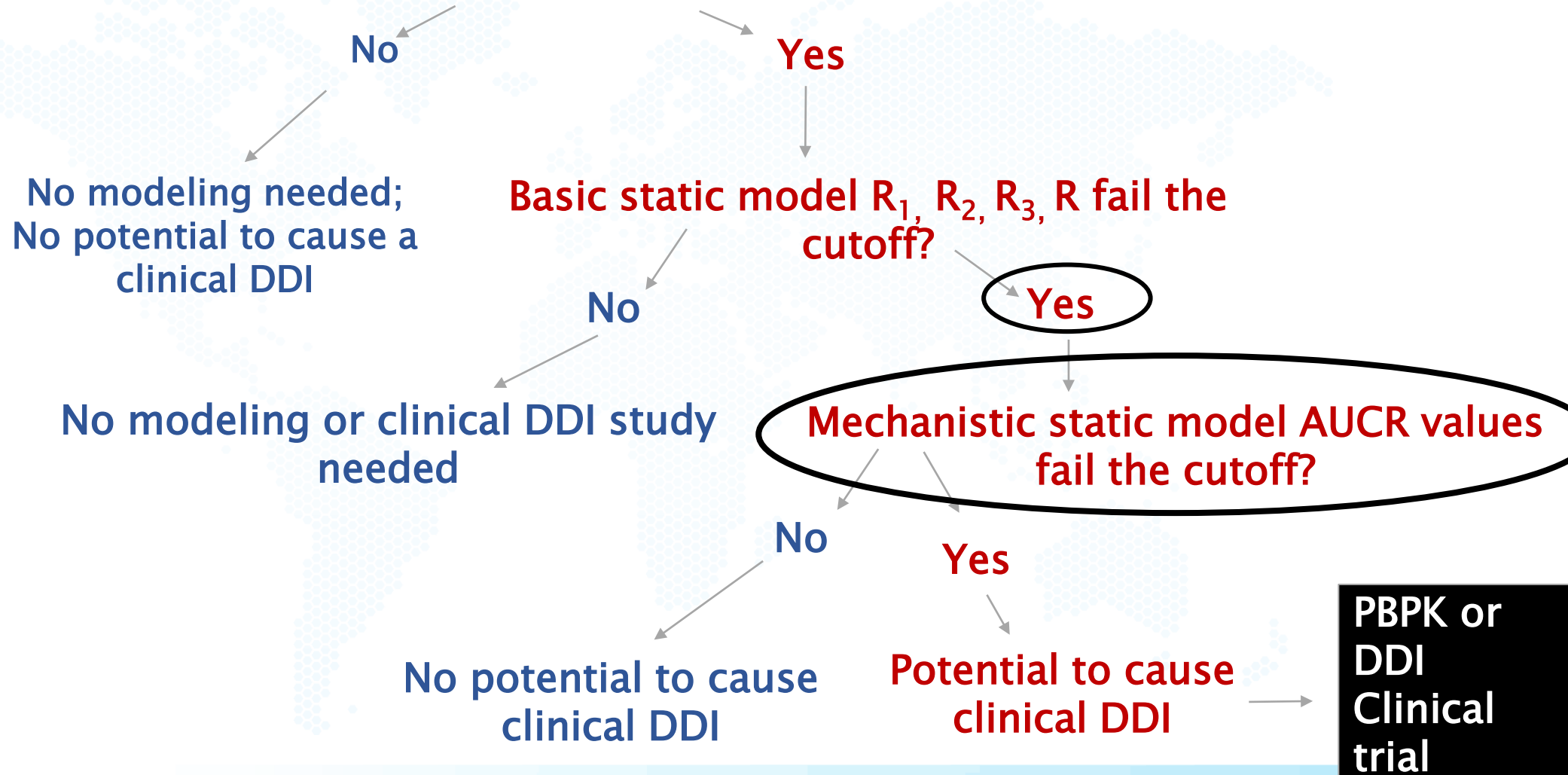
$$R_1 = 1 + (0.025 / 0.0084)$$

$$R_1 = 3.98$$

Potential to inhibit clinically if $R_1 \geq 1.02$

In vitro CYP Inhibition/Induction or Transporter Inhibition?

Decision Tree



Parameters Needed for Static Mechanistic Model

- Dose (μM)
- Maximal unbound total systemic ($I_{\text{max},u}$)
- Fraction of metabolism of a victim drug (f_m)
- Hepatic blood flow (Q_h)
- Blood flow through enterocytes (Q_{en})
- Fraction of absorption (F_a)
- Intestinal availability (F_g)
- Hepatic inlet ($I_{\text{in,max}}$) and enterocytic (I_g) drug concentrations
- Extent of binding to plasma proteins ($f_{u,p}$)
- Blood-to-plasma concentration ratio (R_b)

Static Mechanistic Model

When the same CYP enzyme is **inhibited and induced** by the drug, a **net effect** is calculated.

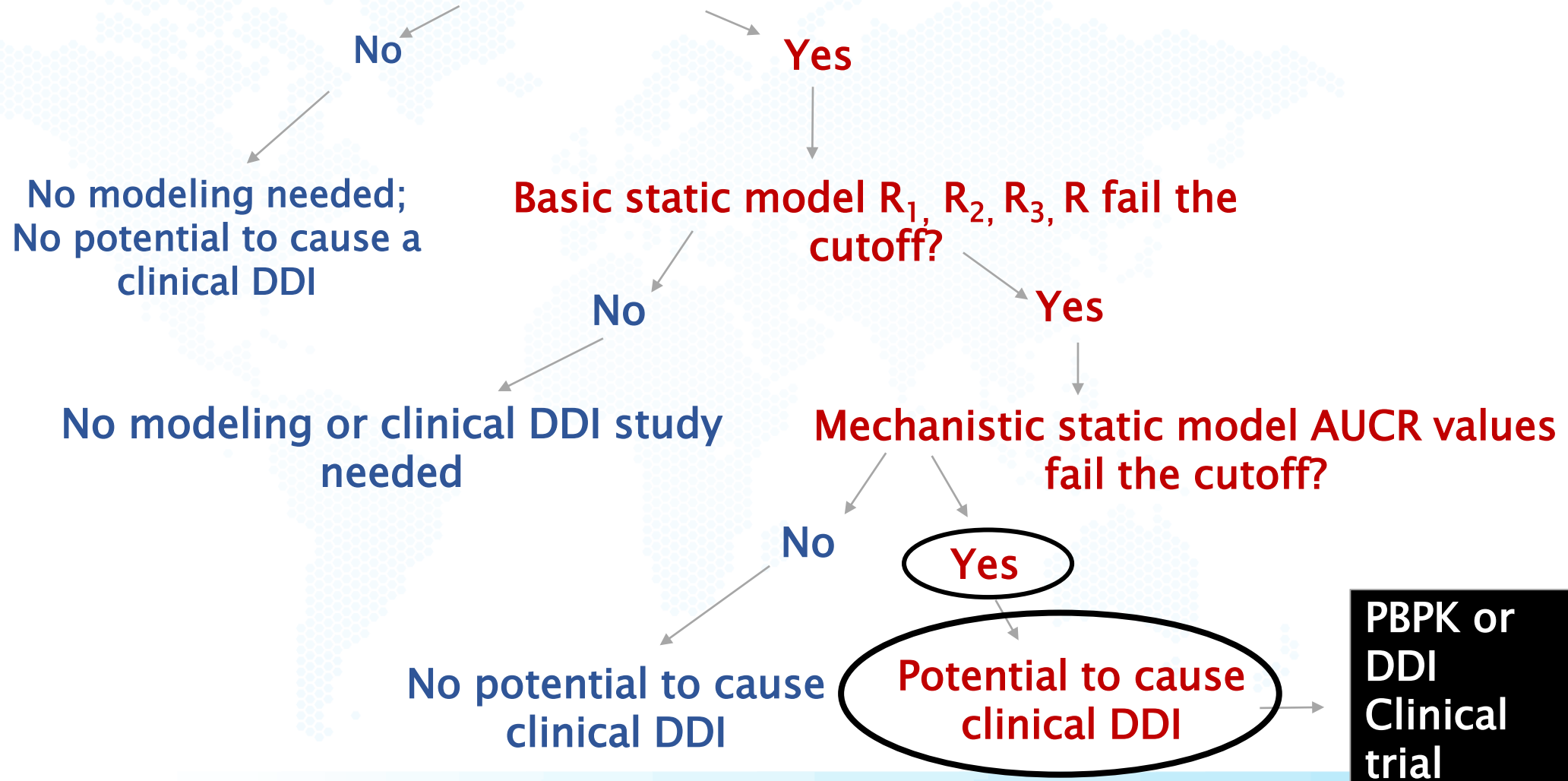
$$AUCR = [1 / ((A_g \times B_g \times C_g) \times (1 - F_g) + F_g)] \times [1 / ((A_h \times B_h \times C_h) \times f_m + (1 - f_m))]$$



Potential to cause induction if $AUCR \leq 0.8$
Potential to cause inhibition if $AUCR \geq 1.25$

In vitro CYP Inhibition/Induction or Transporter Inhibition?

Decision Tree



Previous Consulting Project Example- Overall Modeling Strategy

Enzyme	Substrate (CYP inhibition)	R1 Value (Hepatic direct inhibition)	In vitro hepatocyte culture ID (CYP induction)	R3 Value (CYP induction)	MSM: AUCR Value
CYP2B6	Efavirenz	2.51	HC10-1	0.726	1.002
			HC5-30	0.756	1.002
			HC7-4	0.637	1.002
		≥1.02			≤0.8
				≤0.8	≥1.25

Conclusions:

- Based on the FDA guidance, CYP2B6 was **not predicted to be affected** by the drug to a clinically significant extent.
- C_{max} plasma concentrations would need to be >115-times greater** than the observed C_{max} in patients before AUCR values will fall between the cutoff.

Summary

- Model-based approach to be offered as the follow up to the routine perpetrator potential studies i.e. CYP inhibition, CYP induction, and transporter inhibition
- Great value to the sponsors in assessing the clinical potential that may eliminate the need of conducting clinical studies
- Step-wise approach brings robustness to the prediction

Thank you for watching!

Questions? Get in touch through the **Contact Us** tab on our website

Please contact your regional account manager if you are interested in placing a contracted study or have interest in high-quality test systems for your assays

