



## XenoTech's Products - Supporting Preclinical DMPK

Available products by *in vitro* study type

Applications	In Vitro Test Systems								
	Plateable Cryopreserved Hepatocytes	Pooled Cryopreserved Hepatocytes	Individual Cryopreserved Hepatocytes	Kupffer Cells	Pooled S9/Cytosol/Homogenate	Pooled Microsomes	Individual Microsomes/S9/Cytosol	Recombinant Enzymes (rCYPS)	Lysosomes/Tritosomes
Species Differences	+	+++	++	-	++	++	+	++	++
Metabolic Stability	+++	+++	++	-	++	++	+	+	+++
In Vitro Toxicity	++	-	-	+	+++	+	-	-	-
Enzyme Induction	+++	-	-	-	-	-	-	-	-
P450 Inhibition	+	++	+	-	+	+++	+	++	-
Reaction Phenotyping	+	+	+	-	++	+++	+	++	-
Integrated Metabolism	+++	+++	++	-	++	++	+	-	+
Uptake Transporter Assays	+++	+++	++	-	-	-	-	-	-
Efflux Transporter Assays	+++	-	-	-	-	-	-	-	-
Genetic Polymorphisms	+	+++	+	-	-	-	++	+++	-
3D Hepatic Models	++	-	++	++	-	-	-	-	-

+++ = Preferred/best  
 ++ = Good  
 + = Acceptable  
 - = Not applicable

Order these products online at [www.xenotech.com/products](http://www.xenotech.com/products) or contact a customer service representative at **913.438.7450**



## XenoTech's Products - Supporting Preclinical DMPK

### Available Products by *In Vitro* Study Type

<i>In Vitro</i> Study Type	Available Sekisui XenoTech Products
Enzyme Induction	Cryoplateable human/animal hepatocytes, <i>CryostaX</i> <sup>TM</sup> Plateable Pools, fresh human/animal hepatocytes, Ready-Plate <sup>TM</sup> hepatocytes, treated animal liver microsomes
Hepatic Uptake	<i>HepatoSure</i> <sup>TM</sup> pooled human hepatocytes (n=100) or <i>CryostaX</i> <sup>TM</sup> (n=10/20), <i>CryostaX</i> <sup>TM</sup> Plateable Pools (n=5/10), cryoplateable human hepatocytes, Transporter certified human hepatocytes
Hepatic Efflux	Transporter certified human hepatocytes, <i>CryostaX</i> <sup>TM</sup> Plateable Pools, cryoplateable human hepatocytes
Enzyme Inhibition	Human liver microsomes: <i>XTreme 200</i> -donor pool Human S9, cytosol: <i>XTreme 200</i> -donor pool Hepatocytes: <i>HepatoSure</i> <sup>TM</sup> , <i>CryostaX</i> <sup>TM</sup> , cryoplateable human/animal hepatocytes, Ready-Plate <sup>TM</sup> hepatocytes Recombinant enzymes: Bactosomes (rCYPs) CYP inhibitors/substrates
Metabolic Stability/Clearance	Human liver microsomes, S9: <i>XTreme 200</i> or 50-donor pool Human liver homogenate Hepatocytes: <i>HepatoSure</i> <sup>TM</sup> , <i>CryostaX</i> <sup>TM</sup> , <i>CryostaX</i> <sup>TM</sup> Plateable Pools, cryoplateable human/animal hepatocytes, Ready-Plate <sup>TM</sup> hepatocytes Animal microsomes, S9, extrahepatic subcellular fractions Lysosomes/tritosomes
Metabolite Characterization	Human liver microsomes, S9: <i>XTreme 200</i> or 50-donor pool, human liver homogenate Hepatocytes: <i>HepatoSure</i> <sup>TM</sup> , <i>CryostaX</i> <sup>TM</sup> , animal cryopreserved/fresh hepatocytes
Species Comparison	Human, monkey, dog, minipig, rabbit, guinea pig, hamster, rat, mouse, etc. microsomes, S9, hepatocytes
Reaction Phenotyping	Human liver microsomes: <i>XTreme 200</i> or 50-donor pool, reaction phenotyping kit Hepatocytes: pooled human hepatocytes Bactosomes (rCYPs)
Reactive Metabolites	Human liver microsomes, S9: <i>XTreme 200</i> or 50-donor pool, human liver homogenate Hepatocytes: <i>HepatoSure</i> <sup>TM</sup> , <i>CryostaX</i> <sup>TM</sup> , human/animal cryopreserved/fresh hepatocytes
Lysosomal Trapping	<i>HepatoSure</i> <sup>TM</sup> , cryoplateable human hepatocytes, <i>CryostaX</i> <sup>TM</sup> Plateable Pools, Fa2N-4 immortalized hepatocytes
Kupffer Cell Mediated Toxicity/Macrophage Stimulation	Kupffer cells/Kupffer cell thawing and culture media
Polymorphic Enzymes	Geneknown <sup>TM</sup> human hepatocytes – CYP2C8, CYP2C9, CYP2C19, CYP2D6, CYP3A5, UGT1A1, OATP1B1 Genotyped human liver microsomes – CYP2C9, CYP2C19, CYP2D6, CYP3A5, UGT1A1, UGT1A9
Gender Differences/ Demographic Differences	Male/female specific lots of liver microsomes, across all species Individual donor human liver microsomes and hepatocytes Ethnic and age specific pools available
Enzyme Cofactors	<i>RapidStart</i> <sup>TM</sup> NADPH regenerating system
Hepatocyte Media	Cryopreservation thawing kits, resuspension medium, culture medium and incubation medium
Fresh hepatocytes	Fresh plated human/animal hepatocytes 6, 12, 24, 48 and 96-plate well formats Fresh human/animal suspensions
Custom Products	Human/animal cellular and subcellular products, various tissues/organs. By request only.

Order these products online at [www.xenotech.com/products](http://www.xenotech.com/products) or contact a customer service representative at **913.438.7450**