



XenoTech is a premier provider of custom products to over 200 companies worldwide

XenoTech's Custom Products Division has over 15 years of experience preparing hepatocytes and subcellular fractions from a variety of different species. Every year, our dedicated custom products team prepares more than 50 cellular and subcellular products, tailor-made according to our customers' requests. If your ADME research demands quality data from a unique test system, let XenoTech's experienced team prepare your custom product.

Typical Custom Preparations:

- Fresh animal hepatocytes
- Individual human subcellular fractions (poor & extensive metabolizers, age & race-specific, etc)
- Customized human liver microsome pools (activity, etc)
- Animal hepatic & extrahepatic subcellular fractions
- Human extrahepatic subcellular fractions
- Pre-lysate tissue samples
- Fraction preparation from customer provided tissue

Why Choose XenoTech?

- More than 15 years experience
- Dedicated custom products department
- Over 40 characterization assays available
- No minimum set-up fee
- Validated LC-MS/MS characterization methods
- Custom-designed preparation methods

Surplus Custom Product Inventory

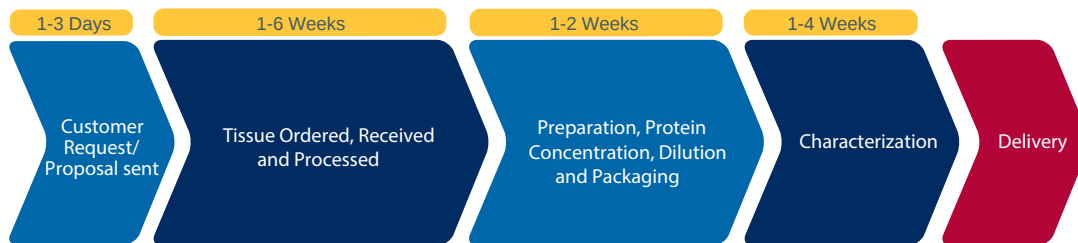
XenoTech routinely prepares custom batches of hepatic, intestinal, pulmonary, and renal subcellular fractions from various species and strains in accordance with client specifications. We have a surplus inventory of products remaining from these custom preparations. You can find a complete list of surplus inventory on our website at www.xenotechllc.com/Products/Custom-Products.

Assurance of Quality

XenoTech's custom products are manufactured and characterized following the principles of GLP regulations and management-approved standard operating procedures. The product proposal, design, preparation and characterization are managed by a dedicated Senior Production Scientist to ensure the highest of quality for each custom product

Typical Custom Product Timeline

Typical custom preparations vary in time to completion. This variation is caused by a number of factors including how long it takes to receive the specific tissue requested and the number of characterization assays requested. A typical custom product preparation usually takes around 4-12 weeks. Below is an example of a typical timeline.



Timeline represents a typical custom product preparation. Timeline varies based on tissue availability and characterization requested

To get started with your custom product quote, contact us at info@xenotechllc.com or call **+913.438.7450 / 877.588.7530** toll-free in North America.

Custom Products At-A-Glance

Species

African Green monkey	Fly	Marmoset monkey
B6C3F1 mouse	Fox Chase SCID mouse	Mongolian gerbil
B6SJLF1 mouse	FVB mouse	New Zealand rabbit
Balb/c mouse	GK rat	NU/NU nude mouse
Beagle dog	Goat	Ob/ob Lean mouse
Brown Norway rat	Golden Syrian hamster	Ob/ob Obese mouse
C57BL/6 mouse	Gottingen minipig	Owl monkey
C57BL/6J mouse	Hanford Minipig	Pig (standard)
Cat	Hartley guinea pig	Rhesus monkey
CD-1 mouse	Horse	SCID mouse
Chicken	Human (many demographics)	Sheep
Cow	Japanese White rabbit	Swiss Webster mouse
Cynomolgus monkey	KK.Cg-A/J mouse	Trout
Db/db Lean mouse	KK/H1J mouse	Wistar Han rat
Db/db Obese mouse	Lewis rat	Yucatan minipig
Fischer 344 rat	Long-Evans rat	ZDF Lean & Obese rat

Tissues

Adrenal gland	Eye	Liver	Skin
Blood Plasma	Heart	Lung	Spleen
Brain	Ileum	Muscle	Stomach
Colon	Jejunum	Ovary	Testicle
Duodenum	Kidney	Pancreas	Etc.

Fractions

Cytosol	Lysosomes	Mitochondria
Hepatocytes	Mast Cells	Pre-lysate
Homogenate	Microsomes	S9 fraction

Characterization Assays

<p>Content Total P450 content by spectral analysis Cytochrome b₅ content by spectral analysis NADPH-cytochrome c reductase activity</p> <p>Cytochrome P450 activity 7-Benzoxoresorufin O-dealkylation 7-Ethoxoresorufin O-dealkylation 7-Methoxoresorufin O-dealkylation 7-Pentoxoresorufin O-dealkylation 7-Ethoxycoumarin O-dealkylation Phenacetin O-dealkylation Coumarin 7-hydroxylation Bupropion hydroxylation Amodiaquine N-dealkylation Diclofenac 4'-hydroxylation S-Mephenytoin 4'-hydroxylation Dextromethorphan O-demethylation Chlorzoxazone 6-hydroxylation Ebastine hydroxylation Testosterone 2α-hydroxylation Testosterone 2β-hydroxylation Testosterone 6β-hydroxylation Testosterone 7α-hydroxylation Testosterone 15α-hydroxylation Testosterone 16α-hydroxylation</p>	<p>Cytochrome P450 activity continued... Testosterone 16β-hydroxylation Midazolam 1'-hydroxylation Lauric acid 12-hydroxylation 4-Nitrophenol hydroxylation Tolbutamide hydroxylation</p> <p>UDP-Glucuronosyltransferase activity 4-Methylumbelliferone glucuronidation 17β-Estradiol 3-glucuronidation Trifluoperazine glucuronidation 1-Naphthol glucuronidation Propofol glucuronidation Morphine 3-glucuronidation Chenodeoxycholic acid 24-glucuronidation Testosterone 17-glucuronidation 7-Hydroxycoumarin glucuronidation 7-Hydroxycoumarin sulfonation</p> <p>Monoamine oxidase activity 5-Hydroxytryptamine oxidation 4 (Dimethylamino)benzylamine oxidation</p> <p>Carboxylesterase activity Clopidogrel hydrolysis Methylprednisolone 21-hemisuccinate hydrolysis</p>	<p>Glutathione S-transferase activity 1-Chloro- 2,4-dinitrobenzene-glutathione conjugation</p> <p>N-Acetyltransferase activity p-Aminobenzoic acid acetylation Sulfamethazine acetylation Dapsone acetylation</p> <p>Aldehyde Oxidase Phthalazine oxidation</p> <p>Lysosomal Enzymes Acid Phosphatase RNase Cathepsin B</p> <p>Marker Substrate Enzymatic Fold Induction CYP1A2 CYP2B6 CYP3A4</p> <p>Uptake Transporter OATP1B1 OATP1B3 OCT1 NTCP</p>
---	--	--

The list above represents products previously prepared by XenoTech. If you don't see your specific need in the list above, or to get started with your custom product quote, contact us at info@xenotechllc.com or call **913.438.7450**