



XENOTECH
A BioIVT Company

HepatoSure®

HepatoSure® pooled cryopreserved human hepatocytes offers the largest donor pool on the market.

HEPATOSURE®

HepatoSure® is XenoTech's new 100-donor pool of cryopreserved human hepatocytes. The largest pool on the market, HepatoSure® provides the best value when it comes to consistency of data reproducibility and the ability to produce assured minimum yields (AMY). HepatoSure® pushes the envelope to deliver higher quality, more consistency and a better overall value than any other pooled hepatocyte product currently available.

Pooled human hepatocytes are a preferred test system in many drug discovery and development applications which require intact cellular systems for *in vitro* testing. Intact hepatocytes contain the major hepatic drug-metabolizing enzymes and co-factors required to evaluate the metabolism and potential drug-drug interactions of drug candidates effectively.

Packaging

- 1.5 mL (Catalog #HCP100.H15)
- 5 million cells AMY (Assured Minimum Yield)
- HepatoSure™ OptiThaw Kit included (Catalog #K8500)

Characterization Provided

Enzyme	Marker Substrate Reaction
CYP1A2	Phenacetin <i>O</i> -dealkylation
CYP2A6	Coumarin 7-hydroxylation
CYP2B6	Bupropion hydroxylation
CYP2C8	Amodiaquine <i>N</i> -dealkylation
CYP2C9	Diclofenac 4'-hydroxylation
CYP2C19	<i>S</i> -Mephenytoin 4'-hydroxylation
CYP2D6	Dextromethorphan <i>O</i> -demethylation
CYP2E1	Chlorzoxazone 6-hydroxylation
CYP3A4/5	Testosterone 6β-hydroxylation
CYP3A4/5	Midazolam 1'-hydroxylation
UGT	7-Hydroxycoumarin glucuronidation
SULT	7-Hydroxycoumarin sulfonation
Transporter	Marker Substrate
OCT1	1-Methyl-4-phenylpyridinium iodide
OATP1B1	Estrone-3-sulfate
OATP1B3	Cholecystokinin-8
NTCP	Taurocholic Acid
Intrinsic Clearance Data	
Aldehyde Oxidase Activity	

- % Viability

HepatoSure® OptiThaw Kit

XenoTech's HepatoSure® pooled human cryopreserved hepatocytes are shipped with a HepatoSure® OptiThaw Kit. This kit contains a nutrient rich media to ensure consistent yield and high viability post-thaw. Each kit is sufficient to thaw and isolate up to three 1.5mL vials of HepatoSure®. When using the OptiThaw kit in conjunction with XenoTech's HepatoSure® pooled cryopreserved human hepatocyte product, you are guaranteed a post-thaw viability of 70% or higher*.

HepatoSure® Comparison

HepatoSure® was tested against a comparable product currently on the market and demonstrated similar viability, enzymatic activity and overall performance. Costing less than the competition, HepatoSure® includes 100 donors, provides the necessary media for thawing and isolation (HepatoSure® Optimization Kit) and delivers high quality, consistent hepatocytes.

Features and Benefits of HepatoSure®

- High quality, reproducible hepatocytes
- 100 donors (reduces interindividual variability)
- High viability and enzyme activity levels
- Maintains viability >50% up to 4 hours
- Extensively characterized using LC/MS/MS methods
- Best overall value on the market
- Study Manager, Study Director and QC review of data

Contact one of our Product Sales Specialists at [913.438.7450](tel:913.438.7450) or email them if you have questions or would like additional information on HepatoSure®.

allproductsales@xenotechllc.com

Order HepatoSure® online at www.xenotech.com.

*This product requires a density gradient after thawing the hepatocytes in order to achieve greater than 70% viability of the recovered hepatocytes. The yield and viability given above are based on experiments performed at XenoTech using our thawing protocol that includes a density gradient fractionation and the HepatoSure® Optimization Kit. Details of XenoTech's hepatocyte thawing protocol can be found at www.xenotech.com. It is recommended to use XenoTech's thawing protocol with this product to ensure high viability of the recovered cells.