



# Hepatic ADME-Tox Products & Research Services

# Portfolio of Products & Services to Advance Research

**ACCELERATE R&D | GAIN REGULATORY APPROVAL | ACHIEVE COMMERCIAL OBJECTIVES**

BioIVT is the leading provider of high-quality hepatic products. For over 25 years, we have set the standard for hepatocytes and offer leading brands including TRANSPORTER CERTIFIED™ and LIVERPOOL® human hepatocytes. Our product offerings also include INVITROCYP™ human liver microsomes, animal hepatocytes and microsomes and INVITROGRO™ hepatocyte media. These research tools are used in our *in vitro* hepatic modeling solutions by analyzing the ADME-Tox properties of New Chemical Entities (NCEs). Our PHASEZERO® Research Services team works collaboratively with clients to provide them with technical expertise and guidance to advance their drug development program.



**Safety**

**Clearance**

**Distribution**

**Interactions**

**Absorption**

**Excretion**

**Toxicity**

**Metabolism**



## Liver Products & Media

Hepatocytes, homogenate, subcellular fractions and tissue available along with related media products

## Custom Products

Providing cell culture services and custom products to support researcher's cell and tissue culture needs

## Application Kits

Duplicate the experimental conditions we have optimized

## Collaborative Research Services

Expertise to help our clients achieve their research objectives

# Answers for All Your Liver-Related Research Questions

We take a collaborative approach to designing and implementing *in vitro* hepatic research programs. Your project will be assigned to an experienced team that will design, implement, and interpret the study results. We provide consulting support to ensure that study results can be used for regulatory submissions and meet the needs of your internal and external stakeholders. We view our clients as partners and have structured our processes to be flexible and adaptable to meet changing project requirements, and at the same time ensure that we provide high-quality deliverables on-time and within budget.

## Innovative Technology



The C-DILI Assay is a novel *in vitro* method to evaluate a compound's potential for cholestatic drug-induced liver injury.

## B-CLEAR Technology

B-CLEAR® Technology is a patented methodology that allows quantitation of biliary clearance in an *in vitro* model which correlates to *in vivo* biliary clearance.



## Experts in Hepatic Research

Led by Kenneth Brouwer, Ph.D., RPh., Vice President, ADME-Tox, our team of scientists provides expertise in hepatic transporters, metabolism, gene regulation, and disease state models. We have developed unique and innovative technologies that improve IVIVE correlations and predict clinical results. Through our academic networks we access subject matter experts who can provide consulting support for research programs.

# ADME-Tox Research Products

BioIVT is the premier provider of hepatic products to help our clients advance their ADME-Tox research programs.

- TRANSPORTER CERTIFIED™ Human Hepatocytes
- LIVERPOOL® Cryoplateable & Cryosuspension Human Hepatocytes
- Cryoplateable Human & Animal Hepatocytes
- Fresh Human and Animal Hepatocytes
- INVITROCYP™ Human Liver Microsomes
- Animal Liver Microsomes & S9
- INVITROGRO™ Hepatocyte Media

## BioIVT has a reputation for excellence:

- ✓ Highest quality standards
- ✓ Ethical procurement of tissues
- ✓ Automated manufacturing processes
- ✓ Largest lots in the industry
- ✓ Extensive inventory options and choices



<b>TRANSPORTER CERTIFIED</b>	<i>The industry standard for hepatocytes having physiologically-relevant transporter function, metabolic competence &amp; appropriate pathway functionality.</i>	<ul style="list-style-type: none"> <li>• Evaluated for uptake &amp; efflux</li> <li>• Extensively tested</li> <li>• High viability &amp; confluence</li> </ul>
<b>LIVERPOOL</b>	<i>Patented methodology that produces multi-donor lots of high-quality, cryopreserved hepatocytes. LIVERPOOL cryoplateable &amp; cryosuspension human hepatocytes are available.</i>	<ul style="list-style-type: none"> <li>• Various donor configurations</li> <li>• Large lot sizes</li> <li>• Customized lots</li> </ul>
<b>Cryoplateable Hepatocytes</b>	<i>Standard isolation methods that consistently produce high-quality cryoplateable hepatocytes which receive enzyme characterization &amp; plateability assessment.</i>	<ul style="list-style-type: none"> <li>• Human &amp; animal</li> <li>• High viability and confluence</li> <li>• Extensive inventory</li> </ul>
<b>INVITROGRO Media</b>	<i>Media that is optimized for thawing, plating &amp; culturing of hepatocytes. Various sizes available.</i>	<ul style="list-style-type: none"> <li>• Animal and human formulations</li> <li>• TORPEDO™ Antibiotic Mix</li> <li>• INVITROGRO™ KHB</li> </ul>
<b>INVITROCYP Microsomes</b>	<i>Human liver microsomes that are designed to meet your unique requirements for inhibition &amp; metabolism studies.</i>	<ul style="list-style-type: none"> <li>• H-CLASS - high activity</li> <li>• M-CLASS - moderate activity</li> <li>• *C-CLASS - customer specified</li> </ul>



## Custom-Cell ADME-Tox Biologicals

We manufacture a variety of primary cells and subcellular products from client-supplied or BioIVT acquired tissues. These Custom-Cell products are prepared to your specifications which can include donor demographics, enzyme characterization, vial configuration, lot size, buffer or media formulations.

BioIVT can purchase and transfer any strain of rats, mice and other commercially available species to our facility to produce various ADME-Tox cellular and subcellular products.

## Common Species Requested

### Rat

- Sprague Dawley
- Han Wistar
- Fischer 344
- Cotton
- Lewis

### Mouse

- ICR / CD-1
- Transgenic Mice

### Guinea Pig

- Dunkin Hartley

### Canine

- Beagle

### Mini-Pig

- Gottingen
- Yucatan

### Primates

- Cynomolgus Monkey
- Rhesus Monkey
- Marmoset Monkey

### Rabbit

- New Zealand White
- Dutch Belted
- Himalayan

### Livestock

- Bovine
- Chicken
- Goat
- Horse
- Pig
- Sheep

### Other

- Bird
- Feline
- Hamster
- Ferret
- Fish
- Woodchuck
- Customer Supplied

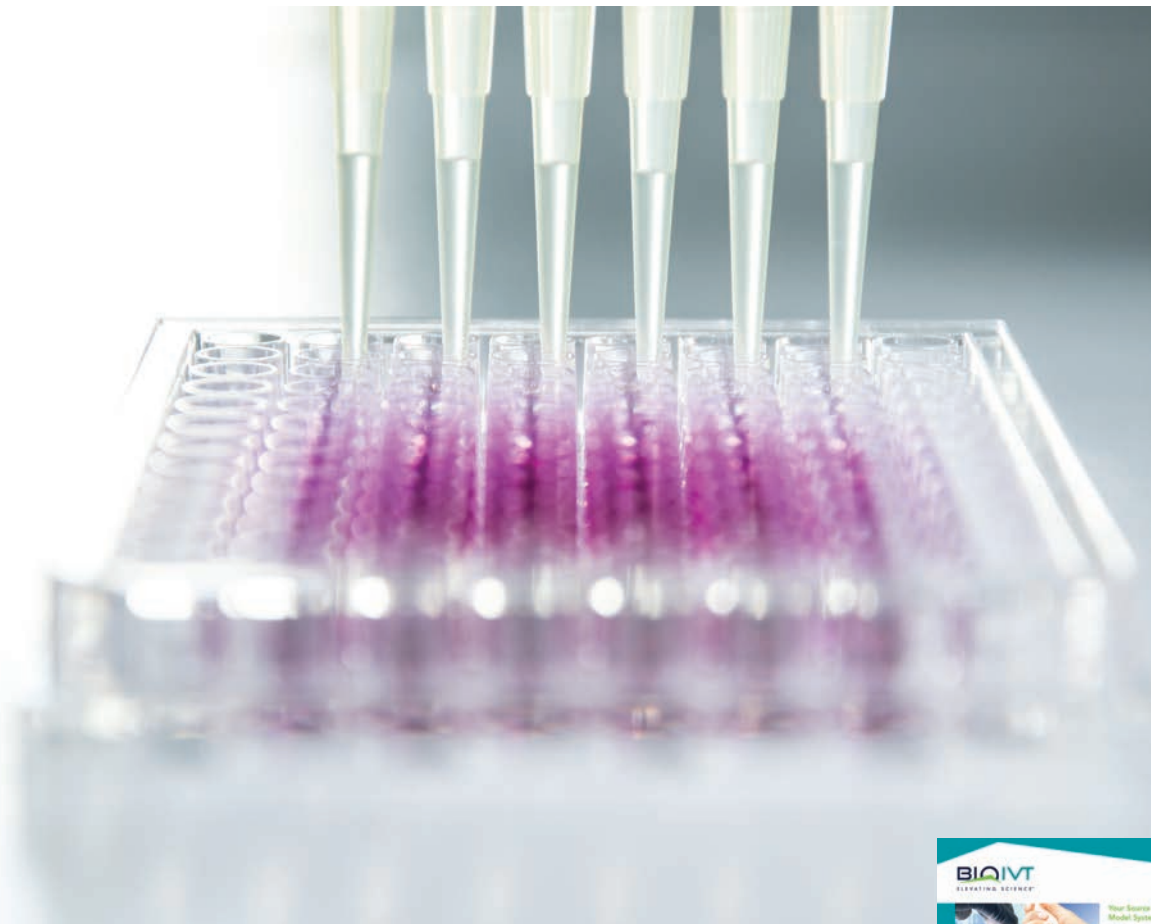
# HepatA!ert™



The HepatAlert is the fastest way to get notified on new lots of hepatocytes :

- TRANSPORTER CERTIFIED™ Hepatocytes
- Cryoplateable Human & Animal Hepatocytes
- Fresh Animal & Human

Sign-up at [www.bioivt.com/hepatalert!](http://www.bioivt.com/hepatalert!)



Visit [BioIVT.com](http://BioIVT.com) to learn more about all the products and services we provide for *in vitro* ADME-Tox research including skin-derived products.



## About BioIVT

BioIVT, formerly BioreclamationIVT, is a leading global provider of high-quality biological specimens and value-added services. We specialize in control and disease state samples including human and animal tissues, cell products, blood and other biofluids. Our unmatched portfolio of clinical specimens directly supports precision medicine research and the effort to improve patient outcomes by coupling comprehensive clinical data with donor samples. Our PHASEZERO® Research Services team works collaboratively with clients to provide target and biomarker validation, phenotypic assays to characterize novel therapeutics, clinical assay development and *in vitro* hepatic modeling solutions. And as the world's premier supplier of ADME-Tox model systems, including hepatocytes and subcellular fractions, BioIVT enables scientists to better understand the pharmacokinetics and drug metabolism of newly discovered compounds and the effects on disease processes. By combining our technical expertise, exceptional customer service, and unparalleled access to biological specimens, BioIVT serves the research community as a trusted partner in ELEVATING SCIENCE.



### For Inquiries:

**North America & Asia Pacific**  
 T +1 516 483 1196 F +1 516 483 4683  
 E [customerservice@bioivt.com](mailto:customerservice@bioivt.com)

**Europe, Middle East & Africa**  
 T +44 (0) 1444 250010 F +44 (0) 1444 250066  
 E [ceurope@bioivt.com](mailto:ceurope@bioivt.com)

[BioIVT.com](http://BioIVT.com)