



*XenoTech offers a wide selection of media to supplement hepatocyte and subcellular products.*

## XenoTech's Newly Optimized Hepatocyte Media!

- **Improved Attachment Efficiency**
- **Extended Expiration Dates**
- **Improved Confluency / Morphology**
- **Convenient Storage Conditions**

### *OptiThaw* Hepatocyte Kit (K8000)

### Opti**THAW**

Sekisui XenoTech's Opti**Thaw** hepatocyte kit contains a nutrient-rich media used in thawing and isolating cryopreserved cells. The Opti**Thaw** kit makes it easy to achieve maximum cell yield and viability for each preparation. The kit includes cell culture media optimized for thawing cryopreserved hepatocytes plus our Opti**Count** solution for determining cell yield and viability counts.

### *OptiThaw* Mouse Hepatocyte Kit (K8100)

### Opti**THAW**

XenoTech's Opti**Thaw** mouse kit is similar to the K8000 kit, but it is specifically formulated to give the optimal yield and percent viability when used with mouse hepatocytes.

### *OptiPlate* Hepatocyte Media (K8200)

### Opti**PLATE**

Sekisui XenoTech's Opti**Plate** media is used when plating attachable and inducible cryopreserved human, dog, monkey, mouse, rat, etc. hepatocytes. It is necessary to re-suspend the cell pellet in Opti**Plate** hepatocyte media before seeding the cells onto culture dishes. Plateable cells resuspended in this media attach to most collagen-coated surfaces.

### *OptiCulture* Hepatocyte Media (K8300)

### Opti**CULTURE**

XenoTech's Opti**Culture** hepatocyte media is recommended for maintaining attachment of cryopreserved human, Beagle dog, Cynomolgus monkey, CD1 mouse, Sprague-Dawley rat cultured hepatocytes, as well as fresh plated hepatocytes. This media is recommended for feeding and dosing of cultured hepatocytes.

### *OptiIncubate* Hepatocyte Media (K8400)

### Opti**INCUBATE**

Sekisui XenoTech's Opti**Incubate** hepatocyte media is recommended for both fresh and cryopreserved hepatocyte suspension and plated incubations.

### *OptiMatrix* Hepatocyte Overlay (K8600-K8650)

### Opti**MATRIX**

XenoTech's Opti**Matrix** increases confluency, preserves viability, structure and appearance, assists network formation, and improves excretory function when plating hepatocytes.

### *OptiThaw* Kupffer Cell Media (K8700)

### Opti**THAW**

Sekisui XenoTech's Opti**Thaw** Kupffer Cell media is a nutrient-rich media used to both thaw and culture Kupffer cells.

## NADPH Regenerating System

### *RapidStart*<sup>™</sup>: NADPH-Regenerating System

Say goodbye to ordering separate components for an NADPH regenerating system. XenoTech's RapidStart is a one-step regenerating system. Simply add water and the system is ready to go!

RapidStart is packaged in a single 7mL, durability-tested polycarbonate vial. Use the concentration chart included with the product data sheet to determine the amount of water to add and vortex, that's it. RapidStart has been activated!

The RapidStart<sup>™</sup> system is the most flexible pre-packaged NADPH-regenerating system on the market, allowing the user to easily manipulate the system concentration by varying the amount of water added. Once activated, RapidStart maintains a suitable concentration of NADPH for up to 12 hours at room temperature or 48 hours at 2-4°C, allowing for multi-day use of the regenerating system.



*Add. Agitate. Activate!*

RapidStart is available in two different volumes:

K5000	RapidStart (1.5mL) single vial	
K5000-10	RapidStart (1.5mL) 10-vial box	<b>Save 10%</b>
K5100	RapidStart (0.3mL) single vial	
K5100-5	RapidStart (0.3mL) 5-vial box	<b>Save 15%</b>

*For more information, ask your account representative for the RapidStart flyer*

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