

**CryostaX**

Single Freeze Pooled Cryopreserved Rat Hepatocytes

**RPCH1000 Lot No. 2410059**

Male, Pool of 8

Assured Minimum Yield:  $5.0 \times 10^6$  cells per vial  
Viability: 80%

Livers were perfused and subjected to collagenase digestion for the purpose of hepatocyte isolation. Yield and viability are based on experiments performed at XenoTech using XenoTech's thawing protocol and K8800 Rodent CryostaX OptiThaw Kit.

**Enzyme Activities****Rate**

7-Ethoxycoumarin O-dealkylation	(pmol/million cells/min)	$138 \pm 11$
7-Hydroxycoumarin glucuronidation	(pmol/million cells/min)	$171 \pm 7$
7-Hydroxycoumarin sulfonation	(pmol/million cells/min)	$159 \pm 15$

Values for enzyme activities were determined at a single substrate concentration and are mean  $\pm$  standard deviation of three or more determinations.

To measure metabolic enzyme activities, hepatocytes ( $1 \times 10^6$  /mL) in suspension were incubated in triplicate at  $37 \pm 2^\circ\text{C}$  for 30 minutes in Opti<sup>INCUBATE</sup> medium and 7-ethoxycoumarin (500  $\mu\text{M}$ ). Metabolite formation was determined by validated LC-MS/MS methods with deuterated metabolites as internal standards.

**Animal Information**

Species: Rat  
Strain: International Genetic Standard (IGS), Sprague Dawley  
Sex: Male  
Age: ~ 8-12 weeks  
Vendor: Charles River, Raleigh, NC

Animals were housed in an AAALAC-accredited facility and allowed to acclimate  $\geq$  seven days before use.

Food: Purina 5L79 (*ad libitum*)  
Water: Automatic watering system (*ad libitum*)  
Light/dark cycle: 5:00 am - 5:00 pm, light; 5:00 pm - 5:00 am, dark (12-hour light/dark)  
Temperature:  $70^\circ\text{F} \pm 2^\circ\text{F}$   
Humidity: 30-70 %  
Bedding: Beta Chip (hardwood), NEPCO, Warrensburg, NY  
Cage: Polycarbonate Shoebox Cage, conventional cage

**Store in liquid nitrogen, vapor phase**

CAUTION: This sample should be considered as a potential biohazard and universal precautions should be followed. Intended for *in vitro* use only.

These data were generated by and are the property of XenoTech. These data are not to be reproduced, published or distributed without the express written consent of XenoTech.

This data sheet serves as a Certificate of Analysis and has been approved by **Stephanie Helmstetter, Assistant Director.**  
Signature and Date: Stephanie Helmstetter 13 March 2024