

## **CryostaX**

Single Freeze Pooled Cryopreserved Beagle Dog Hepatocytes

## DPCH1000 Lot No. 2210339

Male, Pool of 3

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

74%

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 K8000 OptiThaw Hepatocyte Kit.

Enzyme Activi	ties
---------------	------

7-Ethoxycoumarin O-dealkylation	(pmol/million cells/min)	522 ± 21
7-Hydroxycoumarin glucuronidation	(pmol/million cells/min)	58.5 ± 7.6
7-Hydroxycoumarin sulfonation	(pmol/million cells/min)	40.4 ± 4.5

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

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

Animal Information		
Species:	Dog	
Strain:	Beagle	
Sex:	Male	
Age:	Sexually Mature	
Vendor:	Marshall Bioresources, North Rose, NY	
Animals were housed	in an AAALAC-accredited facility and allowed to acclimate $\geq$ seven days before use.	
Food:	Nutrena ( <i>ad libitum</i> )	
Water:	Automatic watering system, tap water ( <i>ad libitum</i> )	
Light/dark cycle:	Not monitored	
Temperature:	Ranges from 62°F to 82°F	
Humidity:	Not monitored	
Cage:	Indoor/outdoor run cages, plastic coated rod bottom, sanitized at least every 2 weeks	



## 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, Senior Manager. Signature and Date: <u>Stephanic Helmstetter</u> <u>30 November 2022</u>

Data Shee

Rate