

CryostaX

Single Freeze Pooled Cryopreserved Beagle Dog Hepatocytes

DPCH1500 Lot No. 2210101

Female, Pool of 3

Assured Minimum Yield: 5.0×10^6 cells per vial
 Viability: 75%

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 Activities

Rate

Enzyme Activities	Rate
7-Ethoxycoumarin O-dealkylation (pmol/million cells/min)	388 ± 16
7-Hydroxycoumarin glucuronidation (pmol/million cells/min)	47.3 ± 7.4
7-Hydroxycoumarin sulfonation (pmol/million cells/min)	33.9 ± 3.6

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 (1×10^6 /mL) in suspension were incubated in triplicate at $37 \pm 2^\circ\text{C}$ for 30 minutes in Opti^{INCUBATE} 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: Female
 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 (*ad libitum*)
 Water: Automatic watering system, tap water (*ad libitum*)
 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: Stephanie Helmstetter 17 June 2022