

CryostaX

Single Freeze Pooled Cryopreserved Beagle Dog Hepatocytes

DPCH1000 Lot No. 2210338

Male, Pool of 4

Assured Minimum Yield: 5.0 x 10⁶ 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
7-Ethoxycoumarin O-dealkylation 7-Hydroxycoumarin glucuronidation 7-Hydroxycoumarin sulfonation	(pmol/million cells/min) (pmol/million cells/min) (pmol/million cells/min)	517 ± 49 57.1 ± 10.7 45.6 ± 2.4

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 x 10^6 /mL) in suspension were incubated in triplicate at $37 \pm 2^{\circ}$ C for 30 minutes in Opti^{INCUBATE} medium and 7-ethoxycoumarin (500 μ M). 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 > 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.