

CryostaX

Single Freeze Pooled Plateable Cryopreserved CD1 Mouse Hepatocytes

MPCH1000+ Lot No. 2410055

Male, Pool of 16

Assured Minimum Yield: 2.0×10^6 cells per vial
Viability: 85%

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)	104 ± 4
7-Hydroxycoumarin glucuronidation	(pmol/million cells/min)	219 ± 15
7-Hydroxycoumarin sulfonation	(pmol/million cells/min)	93.6 ± 13.3

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 (1×10^6 /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: Mouse
Strain: CD1
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*)

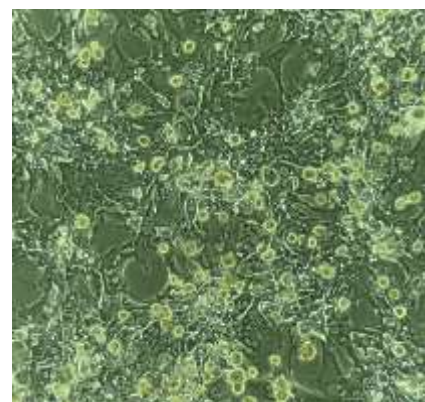
Light/dark cycle: 5:00 am - 5:00 pm, light; 5:00 pm - 5:00 am, dark (12-hour light/dark)

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



MPCH1000+ 2410055 day 3 of culture

	Recommended Seeding Density (million cells/mL)	Recommended Seeding/Feeding Volume Per Well
6 well format	0.6	1.7 mL
12 well format	0.6	650 μL
24 well format	0.6	330 μL
48 well format	Not recommended	
96 well format	Not recommended	



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