

CryostaX

Single Freeze Pooled Cryopreserved Monkey Hepatocytes

PPCH2000 Lot No. 2110025

Cryopreserved Cynomolgus Monkey Hepatocytes

Male, Pool of 3

Assured Minimum Yield: 5.0×10^6 cells per vial
 Viability: 89%

Livers were perfused and subjected to collagenase digestion for the purpose of hepatocyte isolation. Yield and viability are based on experiments performed at Sekisui XenoTech using Sekisui XenoTech's thawing protocol and K8000 OptiThaw CryostaX Kit.

Enzyme Activities		Rate
7-Ethoxycoumarin O-dealkylation	(pmol/million cells/min)	377 ± 8
7-Hydroxycoumarin glucuronidation	(pmol/million cells/min)	98.9 ± 12.5
7-Hydroxycoumarin sulfonation	(pmol/million cells/min)	18.4 ± 0.6

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: Monkey
 Strain: Cynomolgus
 Sex: Male
 Age: 8-9 years
 Vendor: Worldwide Primates, Miami, FL

Animals were housed in an AAALAC-accredited facility.

Imported animals were quarantined for one month prior to shipment into the United States to reduce the risk of importing Ebola virus-infected monkeys. All animals were under veterinary care and were asymptomatic at the time of euthanasia. All of the monkeys tested negative for Simian Retrovirus. None of the animals examined tested positive for any other infectious agents.



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 Sekisui XenoTech. These data are not to be reproduced, published or distributed without the express written consent of Sekisui XenoTech.

Datasheet prepared 15 June 2021