



H1500.H15-3 Lot No. 683

Cryopreserved Human Hepatocytes

Donor Information

Demographics

Gender: Female
Age: 54 Years
Race: Caucasian
Cause of Death: Anoxia

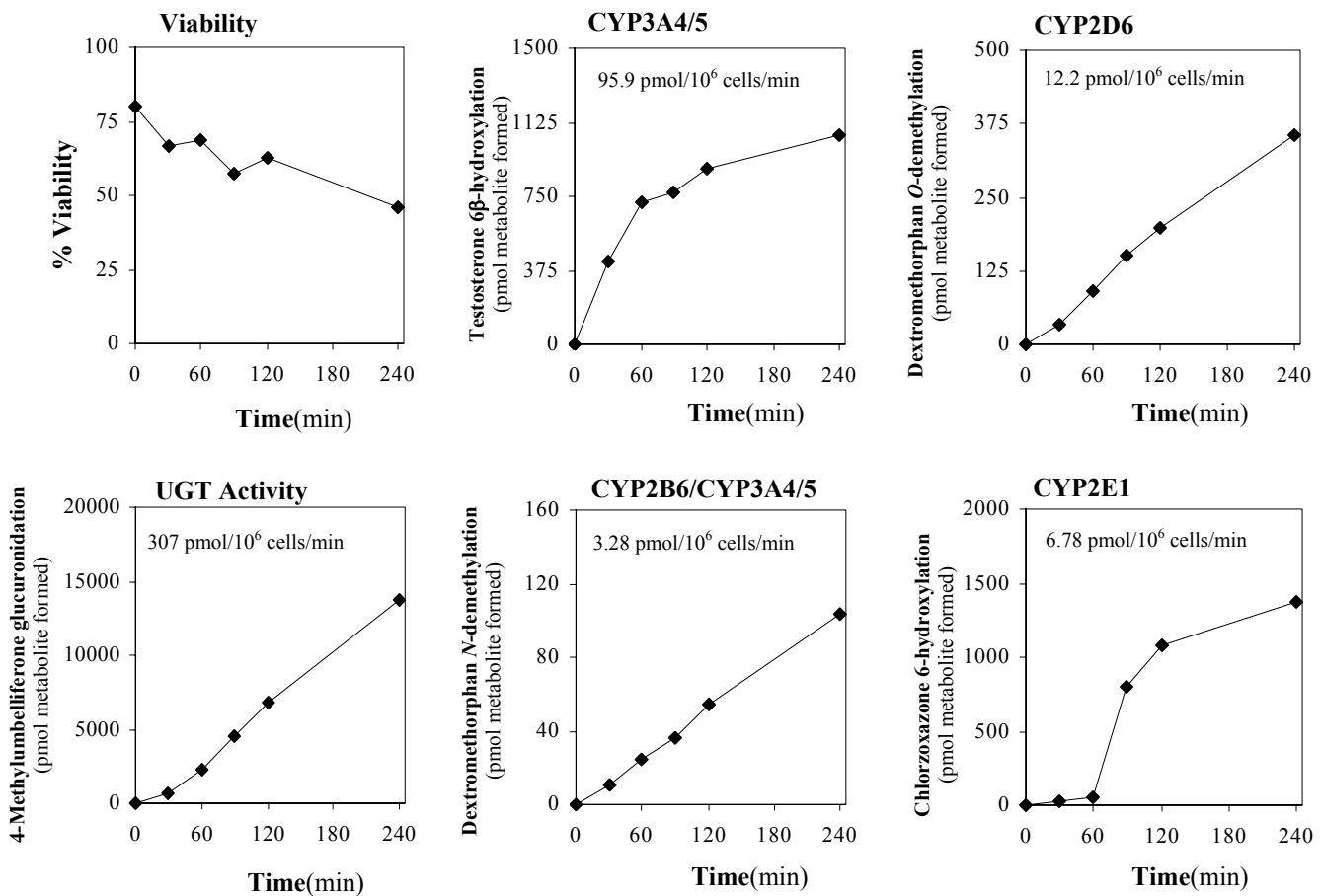
Serology

CMV: (+)
HIV: (-)
HBV: (-)
HCV: (-)

Yield: 3.92 x 10⁶ per vial

(Yield is based on experiments performed at XenoTech using XenoTech's thawing protocol and Hepatocyte Isolation Kit.)

Characterization Data



CAUTION: These hepatocyte samples are from donors who tested negative for HIV and hepatitis. However, we recommend that these samples be considered as potential biohazards and that universal precautions be used when working with human derived products.



Store vials in liquid nitrogen, vapor phase.