

# H1500.H15C      Lot No. HC5-6

Cryopreserved Human Hepatocytes

## Donor Information

### Demographics

**Gender:** Female  
**Age:** 66 Years  
**Race:** Caucasian  
**Cause of Death:** Anoxia

### Serology

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

**Assured Minimum Yield:**  $6.0 \times 10^6$  per vial

**Viability:** 80.2%

(Yield and viability are based on experiments performed at XenoTech using XenoTech's thawing protocol and K2000 Hepatocyte Isolation Kit.)

## Enzyme

## Marker substrate reaction (pmol/cells per inc/min)

CYP1A2	Phenacetin <i>O</i> -dealkylation	204 ± 7
CYP2A6	Coumarin 7-hydroxylation	14.7 ± 2.0
CYP2B6	Bupropion hydroxylation	18.1 ± 1.5
CYP2C8	Amodiaquine <i>N</i> -dealkylation	100
CYP2C9	Diclofenac 4'-hydroxylation	515 ± 58
CYP2C19	<i>S</i> -Mephenytoin 4'-hydroxylation	9.26
CYP2D6	Dextromethorphan <i>O</i> -demethylation	98.0 ± 14.4
CYP2E1	Chlorzoxazone 6-hydroxylation	379 ± 77
CYP3A4/5	Testosterone 6β-hydroxylation	74.5 ± 3.2
CYP3A4/5	Midazolam 1'-hydroxylation	17.0
CYP4A11	Lauric acid 12-hydroxylation	13.8 ± 1.4

**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.**



Data sheet prepared 3/19/09