

# H1000.H15B Lot No. HC3-5A

Cryopreserved Human Hepatocytes

## Donor Information

### Demographics

**Gender:** Male  
**Age:** 45 Years  
**Race:** Caucasian  
**Cause of Death:** Anoxia

### Serology

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

**Assured Minimum Yield:**  $4.0 \times 10^6$  per vial

**Viability:** 84.0%

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

## Enzyme

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

CYP1A2	Phenacetin <i>O</i> -dealkylation	22.1
CYP2A6	Coumarin 7-hydroxylation	45.5 ± 3.6
CYP2B6	Bupropion hydroxylation	13.3 ± 2.0
CYP2C8	Amodiaquine <i>N</i> -dealkylation	25.6 ± 4.0
CYP2C9	Diclofenac 4'-hydroxylation	279 ± 36
CYP2C19	S-Mephenytoin 4'-hydroxylation	0.717 ± 0.020
CYP2D6	Dextromethorphan <i>O</i> -demethylation	11.9 ± 0.7
CYP2E1	Chlorzoxazone 6-hydroxylation	584 ± 84
CYP3A4/5	Testosterone 6β-hydroxylation	82.0 ± 10.1
CYP3A4/5	Midazolam 1'-hydroxylation	26.6 ± 1.5
CYP4A11	Lauric acid 12-hydroxylation	17.7 ± 1.0

**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 2/17/09