

**H1000.H15-3      Lot No. HC3-5**

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:**  $3.0 \times 10^6$  per vial**Viability:** 85.0%

(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	$21.7 \pm 0.6$
CYP2A6	Coumarin 7-hydroxylation	$43.7 \pm 0.3$
CYP2B6	Bupropion hydroxylation	$9.43 \pm 1.84$
CYP2C8	Amodiaquine <i>N</i> -dealkylation	$16.5 \pm 2.6$
CYP2C9	Diclofenac 4'-hydroxylation	$200 \pm 42$
CYP2C19	S-Mephenytoin 4'-hydroxylation	$0.883 \pm 0.081$
CYP2D6	Dextromethorphan <i>O</i> -demethylation	$12.8 \pm 1.3$
CYP2E1	Chlorzoxazone 6-hydroxylation	$397 \pm 31$
CYP3A4/5	Testosterone 6 $\beta$ -hydroxylation	$77.9 \pm 6.6$
CYP3A4/5	Midazolam 1'-hydroxylation	$20.8 \pm 0.8$
CYP4A11	Lauric acid 12-hydroxylation	$37.3 \pm 4.8$

**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