

H2C19.HA / Lot No. 0710444

Human Liver Microsomes

Male, Individual No. 489

0.5 mL at 20 mg protein / mL

Genotype, specific content and activities ^a

Content / Rate

CYP2C19 Allelic variant

CYP2C19*1/*1

Cytochrome P450 (nmol/mg protein)

0.260

Cytochrome b₅ (nmol/mg protein)

0.344

NADPH-cytochrome c reductase (nmol/mg protein/min)

114 ± 0

Enzyme

Marker substrate reaction (pmol/mg protein/min)

CYP1A2	Phenacetin O-dealkylation	498 ± 42
CYP2A6	Coumarin 7-hydroxylation	601 ± 16
CYP2B6	Bupropion hydroxylation	147 ± 9
CYP2C8	Amodiaquine N-dealkylation	922 ± 63
CYP2C9	Diclofenac 4'-hydroxylation	2670 ± 70
CYP2C19	S-Mephenytoin 4'-hydroxylation	31.9 ± 0.4
CYP2D6	Dextromethorphan O-demethylation	210 ± 2
CYP2E1	Chlorzoxazone 6-hydroxylation	1240 ± 70
CYP3A4/5	Testosterone 6β-hydroxylation	812 ± 16
CYP3A4	Midazolam 1'-hydroxylation	40.2 ± 2.0
CYP4A11	Lauric acid 12-hydroxylation	1520 ± 20

^a Values for enzyme activities are mean ± standard deviation of three or more determinations.

Sample	Gender	Age (yrs)	Race	Cause of Death
H0489	Male	46	Caucasian	Cerebrovascular aneurysm

Serology information

- This donor tested negative for cytomegalovirus
- This donor tested negative for HIV, HTLV, HbsAg, and HCV*
- This donor tested negative for RPR**

* Antibody to Human Immunodeficiency Virus, Antibody to Human T Cell Lymphotropic Virus, Hepatitis B Surface Antigen, Antibody to Hepatitis C Virus, respectively.

** Rapid Plasma Reagin.

Data sheet prepared 2/26/08



Store at -80 °C

For in vitro use only

CAUTION: This liver sample is from a donor 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.

These data were generated by and are the property of XENOTECH, LLC. These data are not to be reproduced, published or distributed without the expressed written consent of XENOTECH, LLC.