

H2C19.MA / Lot No. 0710455

Human Liver Microsomes

Female, Individual No. 443

0.5 mL at 20 mg protein / mL

Genotype, specific content and activities ^a

Content / Rate

CYP2C19 Allelic variant

CYP2C19*1/*2

Cytochrome P450 (nmol/mg protein)

0.327

Cytochrome b₅ (nmol/mg protein)

0.229

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

89.8 ± 6.1

Enzyme

Marker substrate reaction (pmol/mg protein/min)

CYP1A2

Phenacetin O-dealkylation

106 ± 17

CYP2A6

Coumarin 7-hydroxylation

132 ± 12

CYP2B6

Bupropion hydroxylation

73.0 ± 5.2

CYP2C8

Amodiaquine N-dealkylation

256 ± 8

CYP2C9

Diclofenac 4'-hydroxylation

1410 ± 70

CYP2C19

S-Mephenytoin 4'-hydroxylation

5.17 ± 0.47

CYP2D6

Dextromethorphan O-demethylation

98.6 ± 5.7

CYP2E1

Chlorzoxazone 6-hydroxylation

941 ± 21

CYP3A4/5

Testosterone 6β-hydroxylation

867 ± 159

CYP3A4

Midazolam 1'-hydroxylation

173 ± 8

CYP4A11

Lauric acid 12-hydroxylation

624 ± 31

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

| Sample | Gender | Age (yrs) | Race | Cause of Death |
|--------|--------|-----------|----------|------------------------|
| H0443 | Female | 64 | Hispanic | Cerebrovascular Stroke |

Serology information

- This donor tested positive 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 2/6/07



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.

16825 West 116th St. | Lenexa KS 66219

913.GET.P450 | fax 913.227.7100 | xenotechllc.com