

H2D6.MA / Lot No. 0710413

Human Liver Microsomes

Male, Individual No. 510

0.5 mL at 20 mg protein / mL

Genotype, specific content and activities ^a Content / Rate

CYP2D6 Allelic variant CYP2D6*17/*29

| | |
|--|---------|
| Cytochrome P450 (nmol/mg protein) | 0.277 |
| Cytochrome b ₅ (nmol/mg protein) | 0.424 |
| NADPH-cytochrome c reductase (nmol/mg protein/min) | 188 ± 3 |

| Enzyme | Marker substrate reaction (pmol/mg protein/min) | |
|----------|---|-------------|
| CYP1A2 | Phenacetin O-dealkylation | 226 ± 20 |
| CYP2A6 | Coumarin 7-hydroxylation | 14.6 ± 1.2 |
| CYP2B6 | Bupropion hydroxylation | 1120 ± 10 |
| CYP2C8 | Amodiaquine N-dealkylation | 1400 ± 50 |
| CYP2C9 | Diclofenac 4'-hydroxylation | 1230 ± 50 |
| CYP2C19 | S-Mephenytoin 4'-hydroxylation | 6.08 ± 0.09 |
| CYP2D6 | Dextromethorphan O-demethylation | 98.9 ± 3.8 |
| CYP2E1 | Chlorzoxazone 6-hydroxylation | 1180 ± 30 |
| CYP3A4/5 | Testosterone 6β-hydroxylation | 2470 ± 40 |
| CYP3A4 | Midazolam 1'-hydroxylation | 224 ± 22 |
| CYP4A11 | Lauric acid 12-hydroxylation | 1470 ± 30 |

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

| Sample | Gender | Age (yrs) | Race | Cause of Death |
|--------|--------|-----------|------------------|----------------|
| H0510 | Male | 41 | African American | Anoxia |

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 2/12/08



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.