

**H0610.RS9 / Lot No. 0510093**

Human Renal S9

Mixed Gender, Pool of 8

1.0 mL at 5 mg protein / mL

Specific activities^a**Rate**

NADPH-Cytochrome <i>c</i> Reductase (nmol/mg protein/min)	7.92 ± 0.37
Lauric Acid 12-hydroxylation (pmol/mg protein/min)	84.2 ± 0.6
Glucuronidation of 4-Methylumbelliferone (nmol/mg protein/min)	18.7 ± 0.1

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

Subcellular fractions were prepared from whole kidney homogenate.

Donor Information for Human Renal S9

Sample	Gender	Age (Yrs)	Race	Cause of Death
1	F	63	Caucasian	Cerebrovascular Stroke
2	M	58	Caucasian	Intracranial Hemorrhage
3	M	60	Caucasian	Head Trauma
4	M	26	Caucasian	Head Trauma
5	M	56	African American	Intracranial Hemorrhage
6	F	63	Caucasian	Intracranial Hemorrhage
7	M	39	Caucasian	Cerebrovascular Stroke
8	F	69	Caucasian	Cerebrovascular Stroke

Serology information

- Five of eight donors tested positive for cytomegalovirus
- All donors tested negative for HIV, HTLV, HbsAg, and HCV*
- One donor tested nonreactive and seven donors tested negative for RPR *

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

**Store at -80°C***For in vitro use only*

CAUTION: These renal 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.

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.

DATASHEET PREPARED 9/7/05