

H0610.S9 0.5 mL at 20 mg/mL
H0620.S9 1.0 mL at 20 mg/mL
H0630.S9 5.0 mL at 20 mg/mL
H0640.S9 50.0 mL at 20 mg/mL

Lot No. 0810471

Pool of Human Liver S9

Mixed Gender, Pool of 50

Suspension medium: 50 mM Tris·HCl,
150 mM KCl,
2 mM EDTA

Specific content and activities	Content / Rate
Cytochrome P450 (nmol/mg protein)	0.118
Cytochrome b ₅ (nmol/mg protein)	0.066
7-Ethoxycoumarin <i>O</i> -dealkylation (pmol/mg protein/min) ^a	186 ± 1
Glucuronidation of 4-methylumbelliferone (nmol/mg protein/min)	15.4 ± 0.5
CDNB (nmol/mg protein/min) ^b	542 ± 8

Please see attached donor information sheets for more details regarding the individual donors in this pool.

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

^a Assay performed at 500 μM substrate concentration.

^b 1-chloro-2,4-dinitrobenzene-glutathione conjugation by Glutathione Transferase



Store at -80°C

CAUTION: This sample should be considered as a potential biohazard and universal precautions should be followed. Intended for in vitro use only.

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

DATASHEET PREPARED 21-OCTOBER-09

Human Liver S9

Donor Information for Human Liver S9, Pool of 50

Sample	Gender	Age (Yrs)	Race	Cause of Death
152	F	67	Caucasian	Cerebrovascular Accident
153	F	63	Caucasian	Cerebrovascular Accident
192	M	37	Caucasian	Head Trauma
227	F	55	Hispanic	Cerebrovascular Accident
229	M	62	Caucasian	Cerebrovascular Accident
264	M	44	Hispanic	Anoxia
279	F	49	Caucasian	Cerebrovascular Accident
285	F	49	Caucasian	Anoxia
291	F	18	Caucasian	Anoxia
310	F	52	Caucasian	Cerebrovascular Accident
311	M	21	Hispanic	Head Trauma
313	F	49	Caucasian	Cerebrovascular Accident
320	F	69	Caucasian	Cerebrovascular Accident
321	M	58	Caucasian	Cerebrovascular Accident
343	M	26	Caucasian	Cerebrovascular Accident
344	M	58	Caucasian	Anoxia
347	F	49	Caucasian	Cerebrovascular Accident
348	F	59	Caucasian	Anoxia
349	M	56	Caucasian	Cerebrovascular Accident
350	M	33	Hispanic	Head Trauma
353	F	63	Caucasian	Head Trauma

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

Human Liver S9

Donor Information for Human Liver S9, Pool of 50

Sample	Gender	Age (Yrs)	Race	Cause of Death
356	F	48	Caucasian	Cerebrovascular Accident
357	M	29	Caucasian	Head Trauma
364	F	42	Caucasian	Head Trauma
365	M	46	Caucasian	Cerebrovascular Accident
379	M	50	Caucasian	Cerebrovascular Accident
380	M	47	Caucasian	Head Trauma
381	F	58	Caucasian	Cerebrovascular Accident
382	F	19	Caucasian	Head Trauma
387	M	60	Caucasian	Cerebrovascular Accident
388	F	63	Caucasian	Cerebrovascular Accident
389	F	42	Caucasian	Cerebrovascular Accident
391	M	76	Caucasian	Cerebrovascular Accident
393	F	30	Caucasian	Anoxia
394	M	53	Asian	Cerebrovascular Accident
397	M	60	Caucasian	Anoxia
398	F	55	Caucasian	Cerebrovascular Accident
399	M	48	Caucasian	Cerebrovascular Accident
401	F	45	Caucasian	Cerebrovascular Accident
410	F	11	Caucasian	Head Trauma
420	M	42	Caucasian	Anoxia

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

Human Liver S9

Donor Information for Human Liver S9, Pool of 50

Sample	Gender	Age (Yrs)	Race	Cause of Death
422	M	69	Caucasian	Cerebrovascular Accident
428	F	57	Caucasian	Head Trauma
432	M	60	African American	Cerebrovascular Accident
433	M	47	Caucasian	Head Trauma
438	M	56	Caucasian	Head Trauma
441	M	63	Caucasian	Cerebrovascular Accident
442	M	49	Caucasian	Head Trauma
445	F	44	Asian	Cerebrovascular Accident
482	F	67	Caucasian	Anoxia

Serology information

- Cytomegalovirus: 29 of 50 donors tested positive.
- RPR*: 49 donors tested negative and 1 donor was not determined.
- HIV, HTLV, HbsAg, and HCV**: All donors tested negative.

* Rapid Plasma Reagin

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

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