

XTreme 200

Lot No. 0810405

Human Liver S9

Pool of 200 (100 Male and 100 Female)

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

H2610.S9	0.5 mL at 20 mg/mL
H2620.S9	1.0 mL at 20 mg/mL
H2630.S9	5.0 mL at 20 mg/mL
H2640.S9	50.0 mL at 20 mg/mL

Specific content and activities		Content
Cytochrome P450	(nmol/mg protein)	0.109
Cytochrome b ₅	(nmol/mg protein)	0.077
Marker substrate reaction		
		[S] (μM)
7-Ethoxycoumarin <i>O</i> -dealkylation	(pmol/mg protein/min)	500
4-Methylumbelliferone glucuronidation	(nmol/mg protein/min)	100
CDNB ^a	(nmol/mg protein/min)	1000
		Rate
		211 ± 31
		34.5 ± 7.0
		432 ± 125

^a1-chloro-2,4-dinitrobenzene-glutathione conjugation by Glutathione Transferase.

Values for enzyme activities were determined at a single substrate concentration and are mean ± standard deviation of three or more determinations.

To measure cytochrome P450 (CYP) activity, liver S9 (200 μg/mL) was incubated in triplicate at 37 ± 1 °C for 10 minutes in potassium phosphate buffer (50 mM, pH 7.4), containing MgCl₂ (3.0 mM), EDTA (1.0 mM), NADP (1.0 mM), glucose-6-phosphate (5.0 mM), glucose-6-phosphate dehydrogenase (1 Unit/mL) and marker substrate, at the final concentrations indicated. Metabolite formation was determined by validated LC-MS/MS methods with deuterated metabolites as internal standards.

To measure UDP-glucuronosyltransferase (UGT) activity, liver S9 (100 μg/mL) was incubated in triplicate at 37 ± 1 °C for 5 or 10 minutes in Tris-HCl (100 mM, pH 7.7 at 37 °C), CHAPS (0.5 mM), EDTA (1.0 mM), MgCl₂ (10 mM), D-saccharic acid 1,4-lactone (100 μM), uridine diphosphate-glucuronic acid (8.0 mM) and marker substrate, at the final concentrations indicated.

To measure the rate of glutathione conjugation, liver S9 (20 μg/mL) was incubated at 37 ± 1 °C in potassium phosphate buffer (100 mM, pH 6.5), glutathione (1mM), and CDNB (1mM).

Each donor is equally represented in this pool.



Store at -80 °C

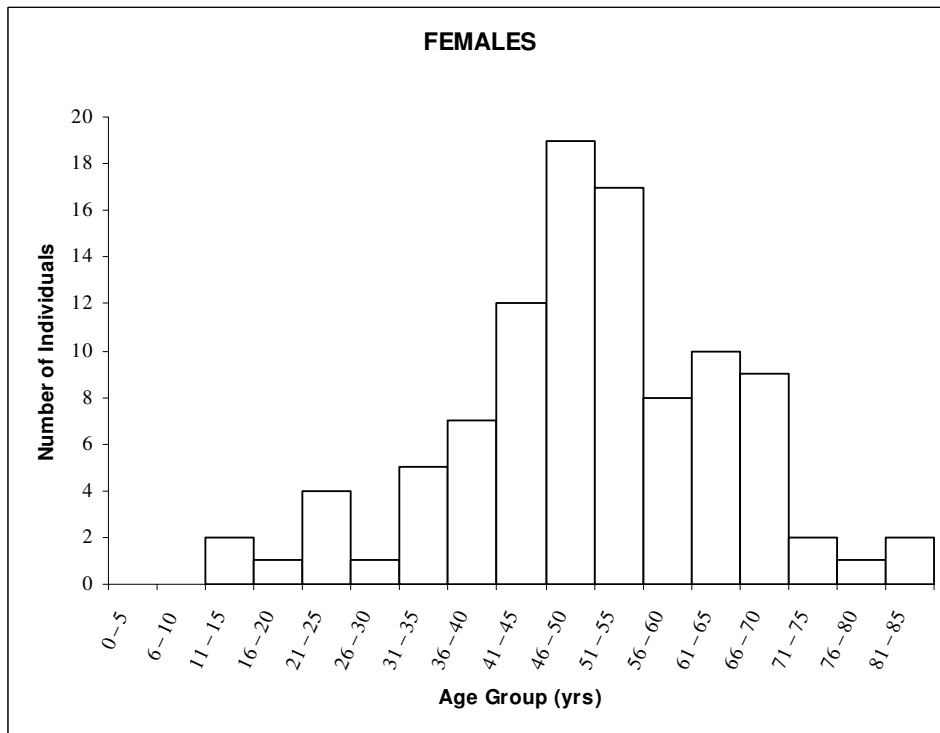
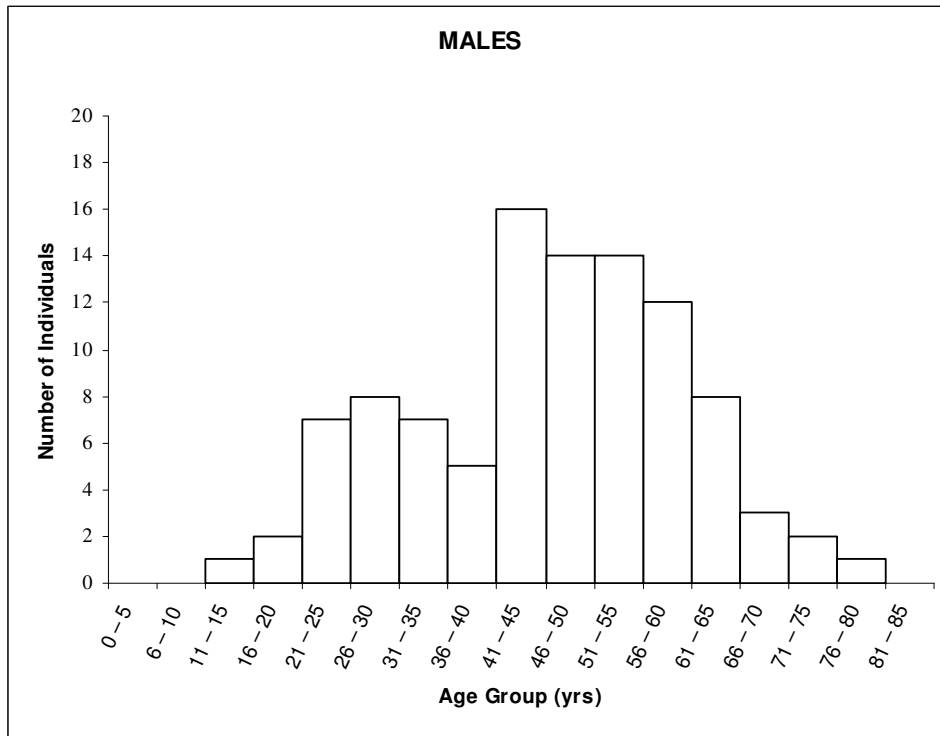
For in vitro use only

CAUTION: This sample should be considered as a potential biohazard and universal precautions should be followed.

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

DATASHEET PREPARED 09-JUNE-09

Age distribution of donors in the pool of 200



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

Human Liver S9

Donor Information for Human Liver S9, Pool of 200

Sample	Gender	Age (Yrs)	Race	Cause of Death
248	M	29	Hispanic	Head trauma
253	M	57	Caucasian	Cerebrovascular accident
255	M	46	Hispanic	Anoxia
309	F	39	Caucasian	Cerebrovascular accident
325	M	54	Caucasian	Head trauma
337	F	57	Caucasian	Cerebrovascular accident
342	F	31	Caucasian	Anoxia
361	M	47	Caucasian	Head trauma
366	M	45	Hispanic	Anoxia
375	M	23	Caucasian	Head trauma
377	M	11	Caucasian	Anoxia
384	M	53	Caucasian	Anoxia
386	M	32	Caucasian	Head trauma
400	F	42	Caucasian	Anoxia
403	F	51	Caucasian	Anoxia
405	M	32	Caucasian	Anoxia
413	M	50	Caucasian	Anoxia
419	M	56	Hispanic	Cerebrovascular accident
427	M	22	Caucasian	Anoxia
444	F	43	Caucasian	Cerebrovascular accident
445	F	44	Asian	Cerebrovascular accident
476	F	46	Caucasian	Cerebrovascular accident
477	F	65	Caucasian	Anoxia
478	F	45	Caucasian	Anoxia
492	F	46	Hispanic	Cerebrovascular accident
506	F	48	Caucasian	Cerebrovascular accident
515	F	59	Caucasian	Cerebrovascular accident
517	F	47	Caucasian	Cerebrovascular accident
530	F	64	Caucasian	Head trauma
531	F	44	Caucasian	Cerebrovascular accident
536	F	34	Caucasian	Anoxia
540	F	54	Caucasian	Head trauma
542	F	53	Caucasian	Cerebrovascular accident
543	F	58	Caucasian	Anoxia
546	F	53	Caucasian	Cerebrovascular accident
550	F	68	Caucasian	Anoxia
554	F	43	Caucasian	Anoxia
556	F	49	Caucasian	Cerebrovascular accident
563	F	52	Caucasian	Cerebrovascular accident
564	F	42	Caucasian	Cerebrovascular accident
565	F	72	Caucasian	Cerebrovascular accident
571	F	45	Caucasian	Anoxia
573	F	47	Caucasian	Cerebrovascular accident
575	M	43	Caucasian	Head trauma
576	F	55	Caucasian	Cerebrovascular accident
577	F	23	Caucasian	Cerebrovascular accident
579	M	46	Caucasian	Cerebrovascular accident
580	M	77	Caucasian	Cerebrovascular accident

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

Human Liver S9

Donor Information for Human Liver S9, Pool of 200

Sample	Gender	Age (Yrs)	Race	Cause of Death
581	M	63	African American	Cerebrovascular accident
582	F	60	Caucasian	Anoxia
583	M	64	Caucasian	Cerebrovascular accident
584	F	44	Caucasian	Head trauma
585	F	44	Caucasian	Anoxia
586	M	57	Caucasian	Anoxia
587	M	44	Caucasian	Cerebrovascular accident
588	M	29	Caucasian	Anoxia
589	F	55	Caucasian	Head trauma
590	F	64	Caucasian	Cerebrovascular accident
591	F	11	Caucasian	Head trauma
592	F	83	Caucasian	Head trauma
593	M	59	Caucasian	Head trauma
594	F	51	Caucasian	Cerebrovascular accident
595	M	37	Caucasian	Head trauma
597	M	41	Caucasian	Anoxia
598	M	55	Caucasian	Anoxia
599	M	51	Caucasian	Cerebrovascular accident
601	M	35	Caucasian	Anoxia
602	F	83	Caucasian	Cerebrovascular accident
603	F	67	Caucasian	Head trauma
605	F	49	Caucasian	Cerebrovascular accident
606	F	76	Caucasian	Cerebrovascular accident
607	M	52	Caucasian	Head trauma
608	F	54	Caucasian	Anoxia
609	F	48	Caucasian	Anoxia
611	F	25	Caucasian	Cerebrovascular accident
612	M	38	Caucasian	Anoxia
613	M	65	Caucasian	Cerebrovascular accident
615	M	55	Caucasian	Cerebrovascular accident
616	M	30	Caucasian	Head trauma
617	F	52	Caucasian	Cerebrovascular accident
618	F	70	Caucasian	Cerebrovascular accident
619	F	45	Caucasian	Cerebrovascular accident
620	M	48	Caucasian	Anoxia
621	M	50	Caucasian	Head trauma
622	M	41	Caucasian	Head trauma
623	F	69	Asian	Cerebrovascular accident
624	M	21	Caucasian	Anoxia
625	F	39	Caucasian	Anoxia
626	M	30	African American	Anoxia
628	F	62	Caucasian	Cerebrovascular accident
629	M	20	Caucasian	Anoxia
630	M	62	Caucasian	Cerebrovascular accident
631	M	24	Caucasian	Head trauma
632	M	57	Caucasian	Anoxia

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

Human Liver S9

Donor Information for Human Liver S9, Pool of 200

Sample	Gender	Age (Yrs)	Race	Cause of Death
633	M	47	African American	Cerebrovascular accident
634	F	63	Caucasian	Cerebrovascular accident
637	F	66	Caucasian	Cerebrovascular accident
659	F	33	Hispanic	Anoxia
686	F	52	Caucasian	Anoxia
708	M	56	Caucasian	Head trauma
715	M	21	Caucasian	Head trauma
717	M	44	African American	Cerebrovascular accident
718	M	29	Caucasian	Head trauma
720	M	49	Caucasian	Head trauma
721	F	52	Caucasian	Cerebrovascular accident
723	F	57	Caucasian	Cerebrovascular accident
726	F	48	Caucasian	Cerebrovascular accident
728	M	39	Caucasian	Cerebrovascular accident
732	M	54	Caucasian	Head trauma
733	M	33	Caucasian	Anoxia
734	M	57	Caucasian	Cerebrovascular accident
735	M	34	Caucasian	Anoxia
736	F	46	Caucasian	Anoxia
737	F	49	African American	Anoxia
738	M	62	Caucasian	Cerebrovascular accident
739	M	66	Caucasian	Cerebrovascular accident
741	F	74	Caucasian	Head trauma
742	M	42	Caucasian	Head trauma
743	M	35	Caucasian	Cerebrovascular accident
745	M	59	Caucasian	Cerebrovascular accident
747	M	43	Caucasian	Anoxia
748	M	64	Caucasian	Cerebrovascular accident
750	F	53	Caucasian	Anoxia
751	M	29	Caucasian	Anoxia
752	M	44	Caucasian	Anoxia
754	F	66	Caucasian	Cerebrovascular accident
755	F	39	Caucasian	Anoxia
756	F	47	Caucasian	Cerebrovascular accident
757	M	54	Caucasian	Cerebrovascular accident
759	F	33	Caucasian	Cerebrovascular accident
760	M	44	Caucasian	Cerebrovascular accident
761	M	70	Caucasian	Cerebrovascular accident
763	F	53	Caucasian	Anoxia
765	F	39	Caucasian	Cerebrovascular accident
766	F	38	Caucasian	Anoxia
767	F	24	Caucasian	Cerebrovascular accident
768	F	63	Caucasian	Cerebrovascular accident
769	M	58	Caucasian	Head trauma
771	F	47	Hispanic	Anoxia
772	M	46	Caucasian	Anoxia

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

Human Liver S9

Donor Information for Human Liver S9, Pool of 200

Sample	Gender	Age (Yrs)	Race	Cause of Death
773	M	46	Caucasian	Head trauma
774	M	55	Caucasian	Cerebrovascular accident
775	M	26	Caucasian	Head trauma
777	M	44	Hispanic	Head trauma
778	M	61	Caucasian	Cerebrovascular accident
779	F	20	Caucasian	Head trauma
780	F	48	African American	Cerebrovascular accident
781	F	15	Caucasian	Cerebrovascular accident
782	M	57	Caucasian	Cerebrovascular accident
783	M	41	Caucasian	Cerebrovascular accident
784	F	63	Caucasian	Head trauma
785	M	47	Caucasian	Head trauma
786	F	27	Caucasian	Head trauma
787	M	58	African American	Cerebrovascular accident
790	M	52	Caucasian	Cerebrovascular accident
794	M	36	Hispanic	Head trauma
795	M	69	Caucasian	Cerebrovascular accident
796	M	29	Caucasian	Anoxia
797	F	60	Caucasian	Cerebrovascular accident
798	M	55	Caucasian	Anoxia
800	F	67	Caucasian	Cerebrovascular accident
802	M	33	Caucasian	Cerebrovascular accident
803	M	23	African American	Head trauma
804	M	57	Caucasian	Cerebrovascular accident
805	F	24	African American	Anoxia
806	M	44	Caucasian	Head trauma
807	F	47	Caucasian	Cerebrovascular accident
809	M	51	Caucasian	Head trauma
810	M	37	Caucasian	Head trauma
811	M	16	Caucasian	Head trauma
812	M	73	Caucasian	Cerebrovascular accident
813	M	45	Caucasian	Anoxia
816	M	55	Hispanic	Head trauma
817	F	31	Caucasian	Anoxia
818	M	48	Hispanic	Cerebrovascular accident
819	F	49	Caucasian	Head trauma
820	F	65	Caucasian	Cerebrovascular accident
821	F	51	Caucasian	Anoxia
822	M	48	Asian	Cerebrovascular accident
823	F	67	Caucasian	Anoxia
824	M	74	Caucasian	Cerebrovascular accident
826	F	55	Caucasian	Cerebrovascular accident
827	M	54	Caucasian	Head trauma
829	M	45	Caucasian	Anoxia
830	M	65	Caucasian	Cerebrovascular accident
831	M	44	Caucasian	Anoxia
832	F	58	Caucasian	Cerebrovascular accident

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

Human Liver S9

Donor Information for Human Liver S9, Pool of 200

Sample	Gender	Age (Yrs)	Race	Cause of Death
833	M	48	Hispanic	Anoxia
834	F	59	Caucasian	Cerebrovascular accident
835	M	23	Hispanic	Head trauma
847	F	55	Caucasian	Cerebrovascular accident
850	F	45	Caucasian	Anoxia
851	F	47	Caucasian	Cerebrovascular accident
858	F	69	Hispanic	Cerebrovascular accident
861	F	61	Caucasian	Head trauma
864	F	46	Caucasian	Cerebrovascular accident
867	F	39	Hispanic	Cerebrovascular accident
869	F	39	Hispanic	Cerebrovascular accident
874	F	65	Caucasian	Cerebrovascular accident
877	F	48	Caucasian	Cerebrovascular accident

Serology information

- Cytomegalovirus: 110 of 200 donors tested positive and 2 donors were not determined.
- RPR*: 198 donors tested negative and 2 donors were 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. These data are not to be reproduced, published or distributed without the express written consent of XenoTech.