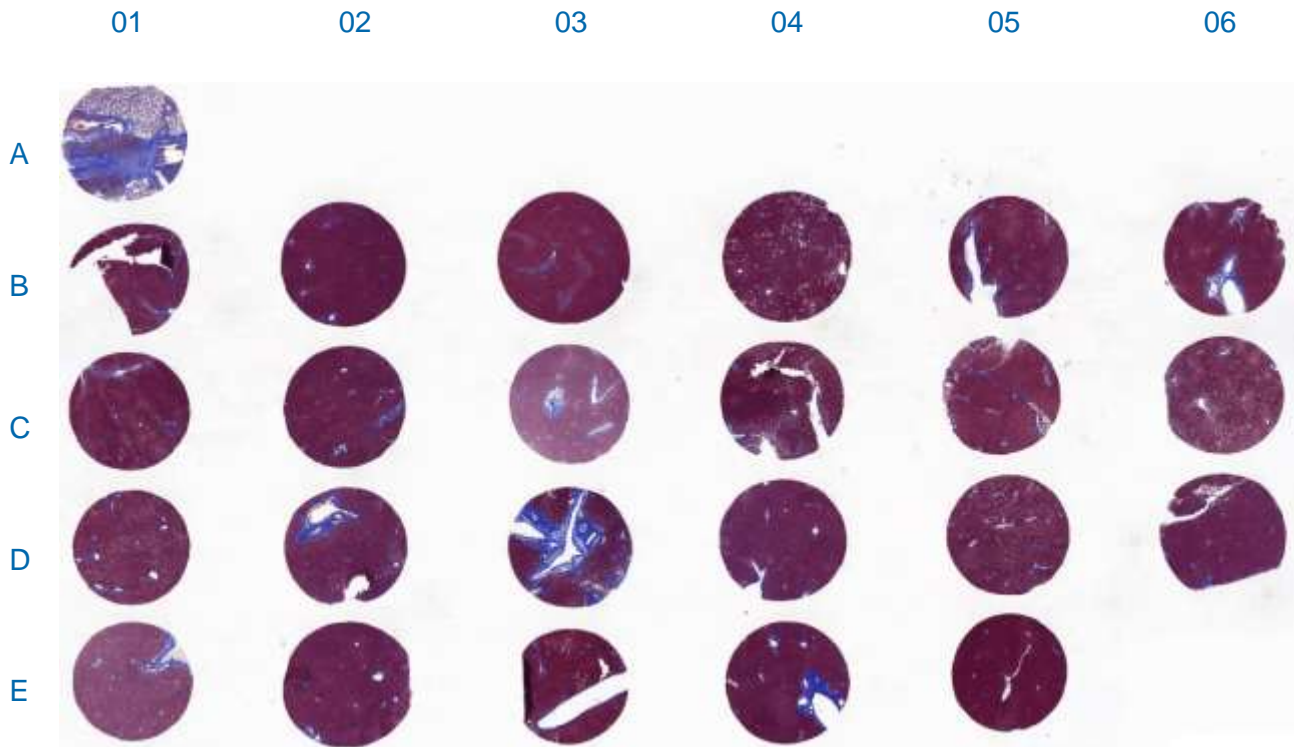


TMA.NORM Lot No. 1910053

Tissue Microarray – Donors with Normal Liver Histology

3 mm diameter cores, 4 µm thick, 23 cases, 1 uncharacterized control



The microarray is provided unstained. The image above is a Masson's trichrome staining of the microarray, which was generated to illustrate histological characteristics of individual samples.



Store at 2-4°C for up to 3 months after date of receipt

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. These data are not to be reproduced, published or distributed without the express written consent of XenoTech.

Datasheet prepared 10 June 2019

Donor Information

Core	Donor	Pathology	Gender	Age (Yrs)	Ethnicity	Macro Fat (%)	BMI	Alcohol Consumption
A01	Control	Uncharacterized	NA	NA	NA	NA	NA	NA
B01	H0510	Normal, Down syndrome	Male	41	African American	0	24.8	None
B02	H0529	Normal	Male	26	Caucasian	0	24.5	None
B03	H0805	Normal	Female	24	African American	0	54.6	None
B04	H0881	Normal, pulmonary fibrosis	Male	48	Caucasian	1	27.3	None
B05	H0898	Normal	Female	58	African American	0	31.3	None
B06	H0919	Normal	Female	39	Caucasian	<1	32.9	Occasional
C01	H1261	Normal	Male	43	Caucasian	0	20.1	Occasional
C02	H1263	Normal	Male	18	Caucasian	0	25.3	None
C03	H1265	Normal	Male	26	Caucasian	0	21.1	Occasional
C04	H1272	Normal	Female	20	Caucasian	0	21.2	Occasional
C05	H1279	Normal	Male	21	Caucasian	1	24.6	None
C06	H1285	Normal, NIDDM	Male	31	Hispanic	3	36.3	Occasional
D01	H1290	Normal	Male	51	Hispanic	0	30.2	Occasional
D02	H1291	Normal	Male	40	Caucasian	<1	29.0	Occasional
D03	H1295	Normal	Male	45	Caucasian	0	29.8	Occasional
D04	H1307	Normal, coronary artery disease	Male	50	Asian	<1	26.4	None
D05	H1314	Normal	Male	43	Caucasian	0	30.0	None
D06	H1318	Normal	Male	63	Caucasian	0	27.4	None
E01	H1321	Normal	Female	49	Caucasian	<1	30.3	None
E02	H1330	Normal, cerebral palsy	Female	56	Asian	0	37.9	None
E03	H1338	Normal, idiopathic lung disease	Female	59	Caucasian	1	32.7	Social
E04	H1362	Normal	Male	56	Caucasian	0	13.7	None
E05	H1400	Normal	Male	41	Caucasian	0	25.7	None

Alcohol Consumption Terms (Listed Highest to Lowest)
Heavy > Moderate > Social > Occasional > None

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.