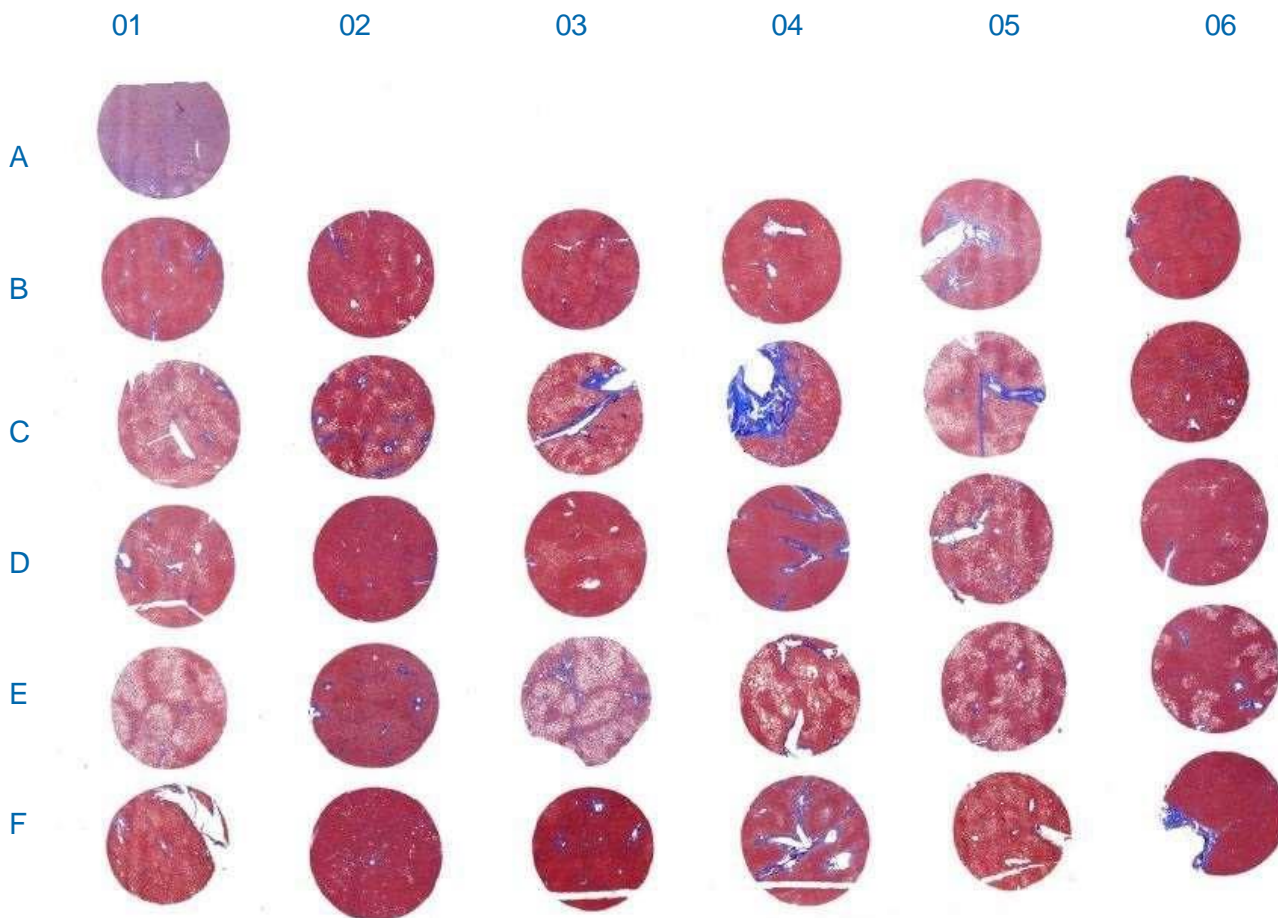


## TMA.NIDDM Lot No. 1910005

Tissue Microarray – Diabetes, Non-insulin dependent, Liver

3 mm diameter cores, 4 µm thick, 25 cases, 5 controls, 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 25 April 2019

## Donor Information

Core	Donor	Pathology	Gender	Age (Yrs)	Ethnicity	Macro Fat (%)	BMI	Diabetes	Alcohol Consumption
A01	Control	Uncharacterized	NA	NA	NA	NA	NA	NA	NA
B01	H0528	Normal, areas of focal necrosis	Male	60	Caucasian	0	27.6	NIDDM	None
B02	H0531	Steatosis, mild	Female	44	Caucasian	20	42.9	NIDDM	None
B03	H0536	Normal	Female	34	Caucasian	1	35.9	NIDDM	Moderate
B04	H0548	Normal	Female	61	Caucasian	0	39.1	NIDDM	Heavy
B05	H0550	Normal, portal inflammation and expansion	Female	68	Caucasian	1	34.2	NIDDM	Occasional
B06	H0572	Normal	Male	68	Caucasian	0	32.2	NIDDM	Moderate
C01	H0574	Steatosis, severe	Male	63	Caucasian	70	34.8	NIDDM	None
C02	H0576	Steatosis, moderate	Female	55	Caucasian	40	48.1	NIDDM	Occasional
C03	H0594	Steatosis, mild	Female	51	Caucasian	20	35.5	NIDDM	None
C04	H0605	Steatohepatitis, bridging fibrosis	Female	49	Caucasian	3	29.7	NIDDM	Occasional
C05	H0608	Steatosis, severe	Female	54	Caucasian	80	40.7	NIDDM	None
C06	H0615	Normal, focal necrosis	Male	55	Caucasian	0	33.5	NIDDM	None
D01	H0630	Normal	Male	62	Caucasian	<5	28.0	NIDDM	Moderate
D02	H0632	Normal	Male	57	Caucasian	1	37.1	NIDDM	Social
D03	H0736	Normal	Female	46	Caucasian	0	39.9	NIDDM	None
D04	H0761	Normal, portal inflammation, infectious monocyctosis	Male	70	Caucasian	0	34.1	NIDDM	None
D05	H0804	Steatosis	Male	57	Caucasian	20	30.5	NIDDM	None
D06	H0833	Normal	Male	48	Hispanic	0	50.0	NIDDM	Social
E01	H0847	Steatohepatitis	Female	55	Caucasian	50-60	36.9	NIDDM	None
E02	H0944	Normal, macro fat in central region	Male	49	Caucasian	1	41.8	NIDDM	Occasional
E03	H0958	Steatohepatitis, occasional ballooned hepatocytes	Male	47	Caucasian	90	40.7	NIDDM	Occasional
E04	H1027	Steatohepatitis, scattered ballooned hepatocytes	Female	63	Caucasian	50	38.9	NIDDM	None
E05	H1081	Steatohepatitis, scattered ballooned hepatocytes	Male	55	Caucasian	80	32.5	NIDDM	Social
E06	H1253	Steatosis	Male	54	Caucasian	5-10	36.1	NIDDM	None
F01	H1285	Normal	Male	31	Hispanic	1	36.3	NIDDM	Occasional
F02	H0881	Control - Normal	Male	48	Caucasian	<1	27.3	No	None
F03	H0898	Control - Normal	Female	58	African American	0	31.3	No	None
F04	H0947	Control - Normal	Female	45	Caucasian	0	35.9	No	None
F05	H1005	Control - Normal	Male	45	Caucasian	3	37.0	No	None
F06	H1008	Control - Normal	Male	36	Hispanic	0	52.6	No	None

Alcohol Consumption Terms (Listed Highest to Lowest)

Heavy &gt; Moderate &gt; Social &gt; Occasional &gt; 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.