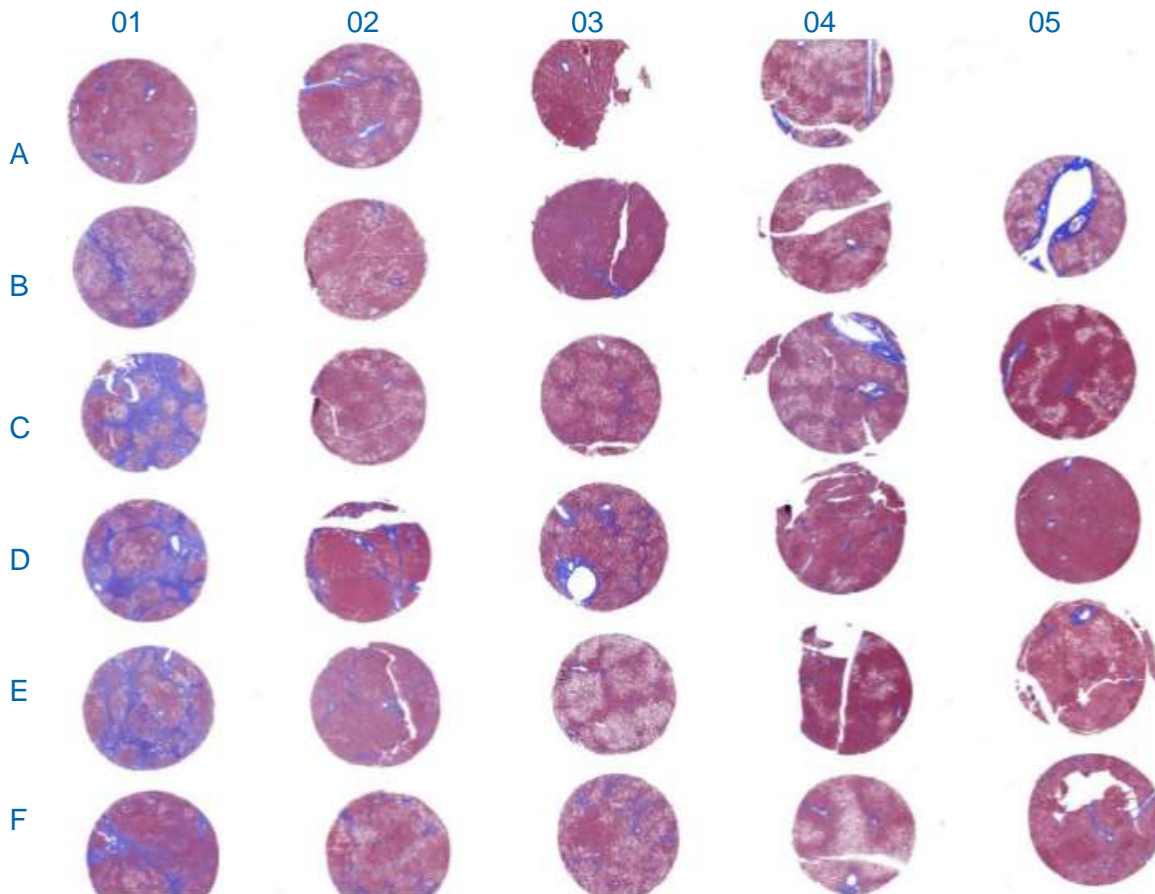


## TMA.FIB Lot No. 2210004

Tissue Microarray – Donors with Liver Fibrosis

3 mm diameter cores, 4 µm thick, 25 cases, 4 controls



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.

This data sheet serves as a Certificate of Analysis and has been approved by **Stephanie Helmstetter, Senior Manager.**  
Signature and Date: Stephanie Helmstetter 25 March 2022

## Donor Information

Core	Donor	Pathology/Fibrosis*	Gender, Age (Yrs), Ethnicity	Macro Fat (%)	BMI	Diabetes	Alcohol Consumption
A01	H0527	Steatohepatitis, perisinusoidal fibrosis, bridging fibrosis (3)	F, 36, C	50	17.6	No	Heavy
A02	H0708	Steatosis, mild, perisinusoidal (zone 1 -3), portal/periportal (2)	M, 56, C	10	48.2	No	Heavy
A03	H0588	Steatosis, mild, diffused portal/periportal (1)	M, 29, C	30	33.8	No	Moderate
A04	H0599	Steatosis, severe, very mild portal/periportal (1)	M, 51, C	80	37.3	No	Heavy
B01	H0515	Steatohepatitis, ballooned hepatocytes, bridging fibrosis (3)	F, 59, C	80	23.4	No	Heavy
B02	H0603	Steatosis, mild, perisinusoidal (zone 1 -3), portal/periportal (2)	F, 67, C	20	50.8	No	None
B03	H0522	Steatohepatitis suspected, diffused portal/periportal (1)	F, 56, C	3	34.2	Yes, type unknown	None
B04	H0896	Steatosis, moderate, portal/periportal (1)	M, 37, C	50	38.6	No	None
B05	H0838	Steatosis, severe, portal/periportal (1)	M, 51, C	50	41.6	No	None
C01	H0598	Cirrhosis (4)	M, 55, C	40	25.1	No	Heavy
C02	H0507	Steatosis, moderate, perisinusoidal (zone 1 -3), portal/periportal (2)	M, 84, C	50	28.8	No	Occasional
C03	H0342	Steatosis, mild, focal perisinusoidal (zone 1 -3), portal/periportal (2)	F, 31, C	25	42.0	No	Moderate
C04	H0518	Steatosis, moderate, focal perisinusoidal (zone 1 -3), portal/periportal (2)	M, 38, C	50	37.2	No	Social
C05	H0778	Steatosis, mild, near normal – Control	M, 61, C	30	33.5	No	None
D01	H0589	Cirrhosis (4)	F, 55, C	5	20.4	No	Heavy
D02	H0605	Steatohepatitis, bridging fibrosis (3)	F, 49, C	3	30.2	NIDDM	Occasional
D03	H0959	Steatohepatitis, ballooned hepatocytes, bridging fibrosis (3)	M, 48, C	40	32.2	No	Heavy
D04	H1069	NASH, scattered ballooned hepatocytes, focal perisinusoidal (zone 1 -3), portal/periportal (2)	M, 39, C	25	62.7	IDDM	None
D05	H0606	Normal – Control	F, 76, C	0	21.0	IDDM	None
E01	H0563	Cirrhosis (4)	F, 52, C	20	24.2	No	Heavy
E02	H0564	Steatohepatitis, bridging fibrosis (3)	F, 42, C	10	25.2	No	Heavy
E03	H0958	NASH, occasional ballooned hepatocytes, focal perisinusoidal (zone 1 -3), portal/periportal (2)	M, 47, C	90	40.7	NIDDM	Occasional
E04	H0774	Steatosis, mild, mild, diffused periportal and pericentral (1)	M, 55, C	10	30.5	No	None
E05	H0594	Steatosis, mild, near normal – Control	F, 51, C	20	35.5	NIDDM	None
F01	H0525	Cirrhosis (4)	F, 56, C	5	23.1	No	Heavy
F02	H0530	Steatohepatitis, ballooned hepatocytes, bridging fibrosis (3)	F, 64, C	50	17.2	No	Heavy
F03	H0878	Steatosis, mild, focal perisinusoidal (zone 1 -3), portal/periportal (2)	M, 43, C	40	35.4	No	None
F04	H0608	Steatosis, severe, diffused portal/periportal (1)	F, 54, C	80	40.7	NIDDM	None
F05	H0541	Steatosis, moderate, near normal – Control	M, 40, C	40	22.1	No	Heavy

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

\*Fibrosis stage was assigned according to Brunt EM, et al., American Journal of Gastroenterology 9, 2467-2474, 1999

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.