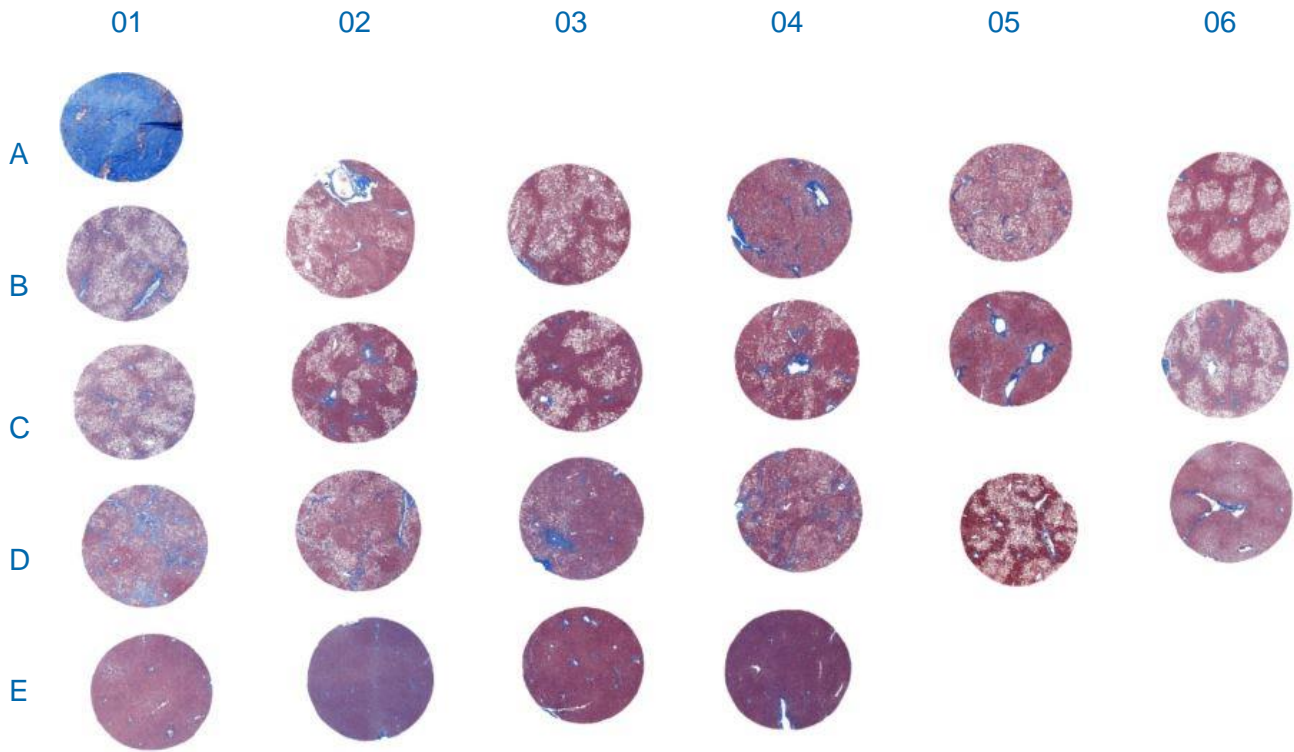


**TMA.NASH Lot No. 1810177**

Tissue Microarray – Liver steatohepatitis with no history of alcohol use  
3 mm diameter cores, 4 µm thick, 12 cases, 10 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 Sekisui XenoTech. These data are not to be reproduced, published or distributed without the express written consent of Sekisui XenoTech.

Datasheet prepared 09 August 2018

## 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	H0448	Steatohepatitis with no history of alcohol use	Female	58	Caucasian	60	32.3	Yes	None
B02	H0499	Steatohepatitis with no history of alcohol use	Female	55	Caucasian	40	41.5	No	None
B03	H0501	Steatohepatitis with no history of alcohol use	Female	58	Caucasian	30	30.2	No	Occasional
B04	H0554	Steatohepatitis with no history of alcohol use	Female	43	Caucasian	25	50.4	No	None
B05	H0656	Steatohepatitis with no history of alcohol use	Female	56	Hispanic	70	34.3	No	None
B06	H0847	Steatohepatitis with no history of alcohol use	Female	55	Caucasian	50 - 60	36.9	Yes	None
C01	H0958	Steatohepatitis with no history of alcohol use	Male	47	Caucasian	90	40.7	Yes	Occasional
C02	H1027	Steatohepatitis with no history of alcohol use	Female	63	Caucasian	50	43.0	Yes	None
C03	H1028	Steatohepatitis with no history of alcohol use	Female	51	African American	70	40.0	Yes	None
C04	H1060	Steatohepatitis with no history of alcohol use	Male	49	Hispanic	50	44.9	Yes	None
C05	H1069	Steatohepatitis with no history of alcohol use	Male	39	Caucasian	25	62.7	No	None
C06	H1097	Steatohepatitis with no history of alcohol use	Male	36	Caucasian	65	37.0	Yes	Occasional
D01	H0515	Control - Steatohepatitis with history of alcohol use	Female	59	Caucasian	80	23.4	No	Heavy
D02	H0530	Control - Steatohepatitis with history of alcohol use	Female	64	Caucasian	40 - 50	17.2	No	Heavy
D03	H0564	Control - Steatohepatitis with history of alcohol use	Female	42	Caucasian	10	25.2	No	Heavy
D04	H0959	Control - Steatohepatitis with history of alcohol use	Male	48	Caucasian	40	32.2	No	Heavy
D05	H1081	Control - Steatohepatitis with history of alcohol use	Male	55	Caucasian	80	32.5	Yes	Social
D06	H1293	Control - Normal	Female	52	Caucasian	0	29.1	No	None
E01	H1307	Control - Normal	Male	50	Asian	0	26.4	No	None
E02	H1310	Control - Normal	Female	27	Caucasian	0	22.0	No	None
E03	H1314	Control - Normal	Male	43	Caucasian	0	30.0	No	None
E04	H1318	Control - Normal	Male	63	Caucasian	0	27.4	No	None

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