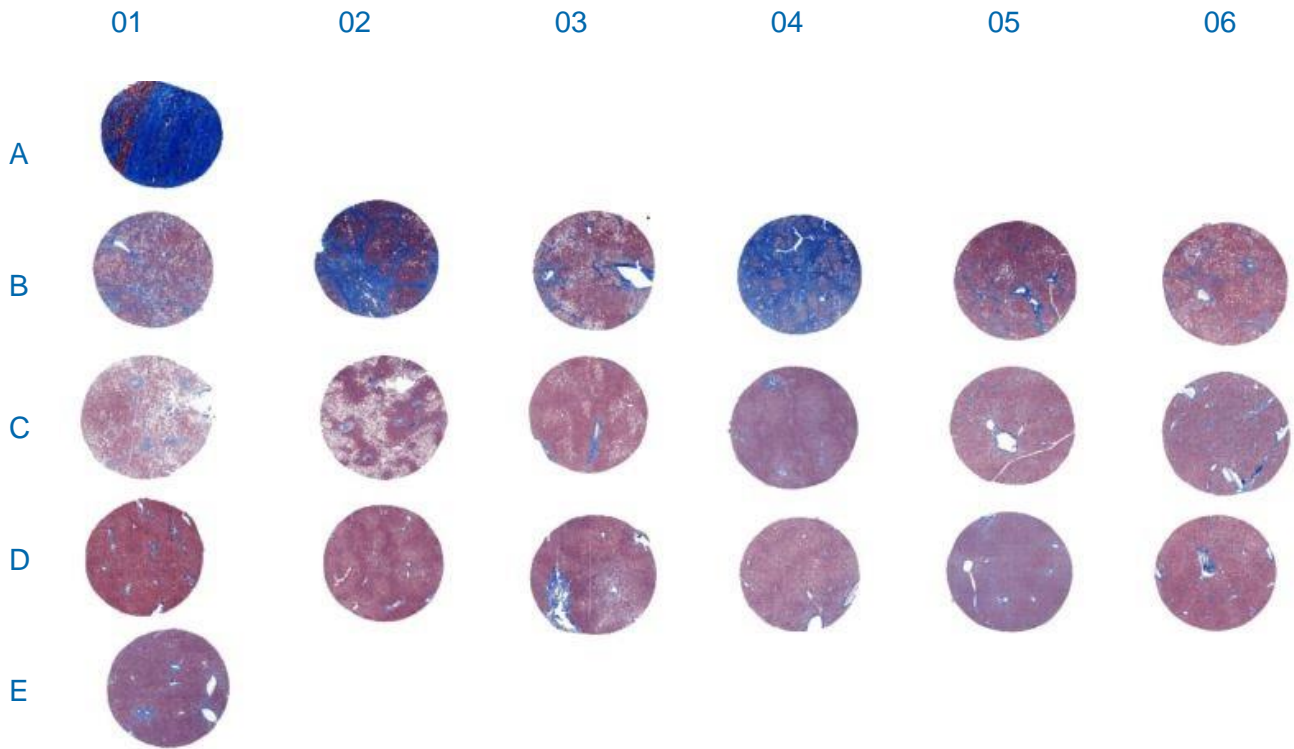


**TMA.ASH Lot No. 1810053**

Tissue Microarray – Liver steatohepatitis with a history of alcohol use  
3 mm diameter cores, 4 µm thick, 9 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 March 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	H0515	Steatohepatitis with a history of alcohol use	Female	59	Caucasian	80	23.4	No	Heavy
B02	H0525	Steatohepatitis with a history of alcohol use	Female	57	Caucasian	5	23.1	No	Heavy
B03	H0530	Steatohepatitis with a history of alcohol use	Female	64	Caucasian	40 - 50	17.2	No	Heavy
B04	H0563	Steatohepatitis with a history of alcohol use	Female	52	Caucasian	20	24.2	No	Heavy
B05	H0564	Steatohepatitis with a history of alcohol use	Female	42	Caucasian	10	25.2	No	Heavy
B06	H0959	Steatohepatitis with a history of alcohol use	Male	48	Caucasian	40	32.2	No	Heavy
C01	H1063	Steatohepatitis with a history of alcohol use	Male	43	Hispanic	80	19.5	No	Heavy
C02	H1081	Steatohepatitis with a history of alcohol use	Male	55	Caucasian	80	32.5	Yes	Social
C03	H1259	Steatohepatitis with a history of alcohol use	Male	64	Caucasian	20	33.2	Yes	Moderate
C04	H0402	Control - Normal with history of alcohol use	Male	32	Caucasian	0	25.7	No	Heavy
C05	H0514	Control - Normal with history of alcohol use	Male	38	African American	0	24.4	No	Heavy
C06	H1277	Control - Normal with history of alcohol use	Female	51	Caucasian	0	22.0	No	Heavy
D01	H1286	Control - Normal with history of alcohol use	Female	56	Caucasian	0	27.2	No	Heavy
D02	H1289	Control - Normal with history of alcohol use	Male	52	Caucasian	0	24.5	No	Moderate
D03	H1293	Control - Normal	Female	52	Caucasian	0	29.1	No	No
D04	H1307	Control - Normal	Male	50	Asian	0	26.4	No	No
D05	H1310	Control - Normal	Female	27	Caucasian	0	22.0	No	No
D06	H1314	Control - Normal	Male	43	Caucasian	0	30.0	No	No
E01	H1318	Control - Normal	Male	63	Caucasian	0	27.4	No	No

Alcohol Consumption Terms (Listed Highest to Lowest)

Heavy > Moderate > Social > Occasional > None