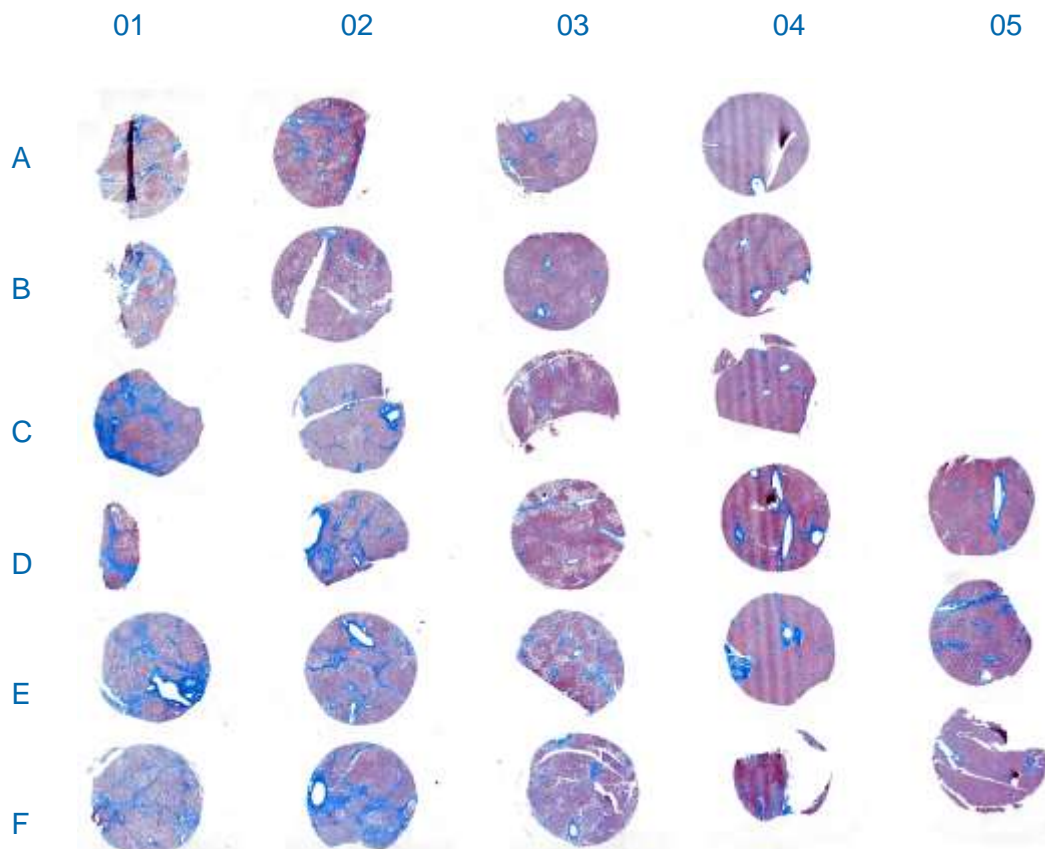


TMA.ASH Lot No. 2210241

Tissue Microarray – Liver steatohepatitis with a history of alcohol use

3 mm diameter cores, 4 µm thick, 9 cases, 8 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.

This data sheet serves as a Certificate of Analysis and has been approved by Stephanie Helmstetter, Senior Manager.

Signature and Date: Stephanie Helmstetter 16 November 2022

Donor Information

Core	Donor	Pathology	Gender	Age (Yrs)	Ethnicity	Macro Fat (%)	BMI	Diabetes	Alcohol Consumption
A01	H0530	Steatohepatitis with a history of alcohol use	Female	64	Caucasian	40 - 50	17.2	No	Heavy
A02	H0959	Steatohepatitis with a history of alcohol use	Male	48	Caucasian	40	32.2	No	Heavy
A03	H1259	Steatohepatitis with a history of alcohol use	Male	64	Caucasian	20	33.2	Yes	Moderate
A04	H1307	Control - Normal	Male	50	Asian	0	26.4	No	No
B01	H0530	Steatohepatitis with a history of alcohol use	Female	64	Caucasian	40 - 50	17.2	No	Heavy
B02	H0959	Steatohepatitis with a history of alcohol use	Male	48	Caucasian	40	32.2	No	Heavy
B03	H1259	Steatohepatitis with a history of alcohol use	Male	64	Caucasian	20	33.2	Yes	Moderate
B04	H1289	Control - Normal with history of alcohol use	Male	52	Caucasian	0	24.5	No	Moderate
C01	H0525	Steatohepatitis with a history of alcohol use	Female	57	Caucasian	5	23.1	No	Heavy
C02	H0564	Steatohepatitis with a history of alcohol use	Female	42	Caucasian	10	25.2	No	Heavy
C03	H1081	Steatohepatitis with a history of alcohol use	Male	55	Caucasian	80	32.5	Yes	Social
C04	H1286	Control - Normal with history of alcohol use	Female	56	Caucasian	0	27.2	No	Heavy
D01	H0525	Steatohepatitis with a history of alcohol use	Female	57	Caucasian	5	23.1	No	Heavy
D02	H0564	Steatohepatitis with a history of alcohol use	Female	42	Caucasian	10	25.2	No	Heavy
D03	H1081	Steatohepatitis with a history of alcohol use	Male	55	Caucasian	80	32.5	Yes	Social
D04	H1277	Control - Normal with history of alcohol use	Female	51	Caucasian	0	22.0	No	Heavy
D05	H1318	Control - Normal	Male	63	Caucasian	0	27.4	No	No
E01	H0515	Steatohepatitis with a history of alcohol use	Female	59	Caucasian	80	23.4	No	Heavy
E02	H0563	Steatohepatitis with a history of alcohol use	Female	52	Caucasian	20	24.2	No	Heavy
E03	H1063	Steatohepatitis with a history of alcohol use	Male	43	Hispanic	80	19.5	No	Heavy
E04	H0514	Control - Normal with history of alcohol use	Male	38	African American	0	24.4	No	Heavy
E05	H1314	Control - Normal	Male	43	Caucasian	0	30.0	No	No
F01	H0515	Steatohepatitis with a history of alcohol use	Female	59	Caucasian	80	23.4	No	Heavy
F02	H0563	Steatohepatitis with a history of alcohol use	Female	52	Caucasian	20	24.2	No	Heavy
F03	H1063	Steatohepatitis with a history of alcohol use	Male	43	Hispanic	80	19.5	No	Heavy
F04	H0402	Control - Normal with history of alcohol use	Male	32	Caucasian	0	25.7	No	Heavy
F05	H1310	Control - Normal	Female	27	Caucasian	0	22.0	No	No

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.