

TMA.NASH Lot No. 2210347

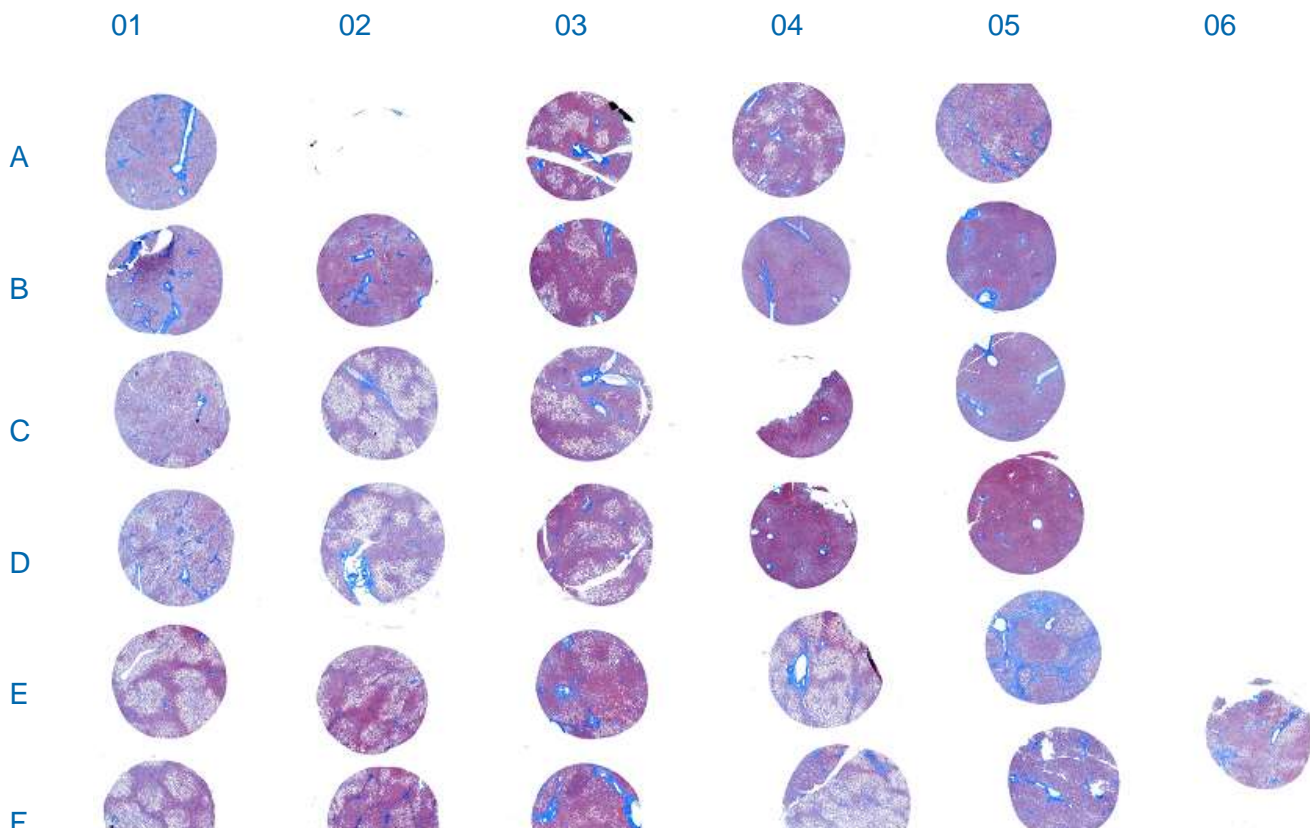
Tissue Microarray – Liver Steatohepatitis with no history of alcohol use

3 mm diameter cores, 4 µm thick

10 cases in duplicate, 11 controls in singlet

Data Sheet

Certificate of Analysis



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 08 December 2022

Donor Information

Core	Donor	Pathology	Gender	Age (Yrs)	Ethnicity	Macro Fat (%)	BMI	Diabetes	Alcohol Consumption
A01	H0554	Steatohepatitis with no history of alcohol use	Female	43	Caucasian	25	50.4	No	None
A02	H1069	Steatohepatitis with no history of alcohol use	Male	39	Caucasian	25	62.7	No	None
A03	H1027	Steatohepatitis with no history of alcohol use	Female	63	Caucasian	50	43.0	Yes	None
A04	H0564	Control - Steatohepatitis with history of alcohol use	Female	42	Caucasian	10	25.2	No	Heavy
A05	H0530	Control - Steatohepatitis with history of alcohol use	Female	64	Caucasian	40 - 50	17.2	No	Heavy
B01	H0554	Steatohepatitis with no history of alcohol use	Female	43	Caucasian	25	50.4	No	None
B02	H1069	Steatohepatitis with no history of alcohol use	Male	39	Caucasian	25	62.7	No	None
B03	H1027	Steatohepatitis with no history of alcohol use	Female	63	Caucasian	50	43.0	Yes	None
B04	H1081	Control - Steatohepatitis with history of alcohol use	Male	55	Caucasian	80	32.5	Yes	Social
B05	H0959	Control - Steatohepatitis with history of alcohol use	Male	48	Caucasian	40	32.2	No	Heavy
C01	H0656	Steatohepatitis with no history of alcohol use	Female	56	Hispanic	70	34.3	No	None
C02	H1097	Steatohepatitis with no history of alcohol use	Male	36	Caucasian	65	37.0	Yes	Occasional
C03	H1028	Steatohepatitis with no history of alcohol use	Female	51	African American	70	40.0	Yes	None
C04	H1307	Control - Normal	Male	50	Asian	0	26.4	No	None
C05	H1336	Control - Normal	Female	60	African American	1-2	24.2	No	Occasional
D01	H0656	Steatohepatitis with no history of alcohol use	Female	56	Hispanic	70	34.3	No	None
D02	H1097	Steatohepatitis with no history of alcohol use	Male	36	Caucasian	65	37.0	Yes	Occasional
D03	H1028	Steatohepatitis with no history of alcohol use	Female	51	African American	70	40.0	Yes	None
D04	H1314	Control - Normal	Male	43	Caucasian	0	30.0	No	None
D05	H1310	Control - Normal	Female	27	Caucasian	0	22.0	No	None
E01	H0847	Steatohepatitis with no history of alcohol use	Female	55	Caucasian	50 - 60	36.9	Yes	None
E02	H0499	Steatohepatitis with no history of alcohol use	Female	55	Caucasian	40	41.5	No	None
E03	H1060	Steatohepatitis with no history of alcohol use	Male	49	Hispanic	50	44.9	Yes	None
E04	H0958	Steatohepatitis with no history of alcohol use	Male	47	Caucasian	90	40.7	Yes	Occasional
E05	H1318	Control - Normal	Male	63	Caucasian	0	27.4	No	None
F01	H0847	Steatohepatitis with no history of alcohol use	Female	55	Caucasian	50 - 60	36.9	Yes	None
F02	H0499	Steatohepatitis with no history of alcohol use	Female	55	Caucasian	40	41.5	No	None
F03	H1060	Steatohepatitis with no history of alcohol use	Male	49	Hispanic	50	44.9	Yes	None
F04	H0958	Steatohepatitis with no history of alcohol use	Male	47	Caucasian	90	40.7	Yes	Occasional
F05	H0515	Control - Steatohepatitis with history of alcohol use	Female	59	Caucasian	80	23.4	No	Heavy
F06	H0530	Control - Steatohepatitis with history of alcohol use	Female	64	Caucasian	40 - 50	17.2	No	Heavy

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.