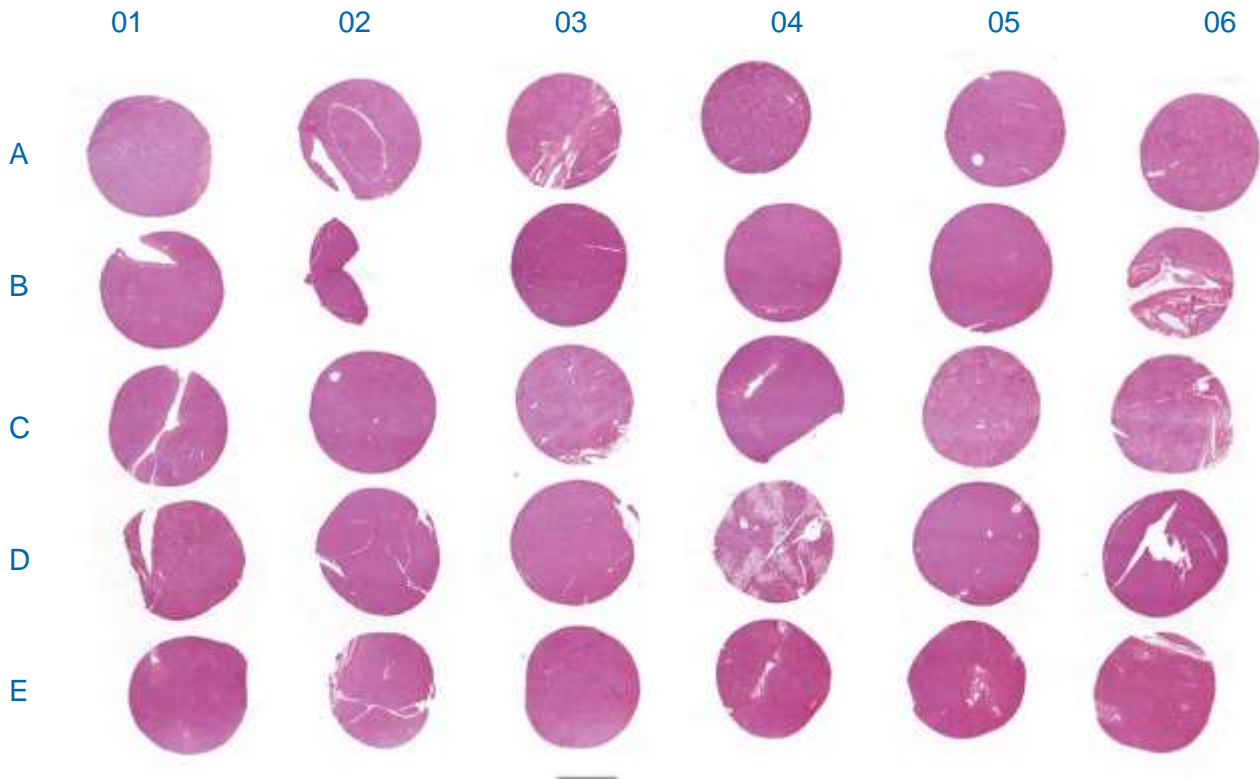


TMA.PED Lot No. 1810266

Tissue Microarray – Pediatric

3 mm diameter cores, 4 µm thick, 25 cases, 5 controls



The microarray is provided unstained. The image above is a hematoxylin and eosin staining of the microarray, which was generated to illustrate histological characteristics of individual samples. This is a representative image.



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 31 January 2019

Donor Information

Core	Donor	Pathology	Gender	Age	Ethnicity	BMI	Alcohol Consumption
A01	H1393	Very near normal	Male	4 months	C	18.8	None
A02	H0395	Hepatocyte degeneration 15-20%	Male	5.5 months	C	17.3	None
A03	H1383	Normal	Male	12 months	C	19.7	None
A04	H0322	Diffused feathery degeneration	Male	19 months	H	14.9	None
A05	H1351	Patchy diffused hepatic cell damage (ischemia)	Female	20 months	H	23.5	None
A06	H1334	Patchy cholestasis, feathery degeneration	Female	21 months	AA	22.8	None
B01	H1397	Feathery and ballooning degeneration	Female	21 months	H	18.9	None
B02	H0551	Near normal	Male	2 years	C	15.5	None
B03	H0852	Diffused hepatic cell death (ischemia)	Male	2 years	H	16.1	None
B04	H0872	Massive liquefaction necrosis	Male	2 years	C	19.3	None
B05	H0346	Necrotic and apoptotic cells ~5%	Male	3 years	C	15.0	None
B06	H0776	Massive necrosis, rapid ischemic event	Female	4 years	AA	18.0	None
C01	H1096	Near normal	Female	6 years	C	13.9	None
C02	H1301	Ischemic changes ~20%, steatosis ~7-10%	Female	7 years	C	14.5	None
C03	H1092	Ischemic damage, single cell necrosis	Female	9 years	A	18.8	None
C04	H0485	Diffused patchy feathery degeneration	Male	10 years	C	20.5	None
C05	H0996	NAFLD (potentially lowest grade of NASH)	Female	10 years	C	32.4	None
C06	H0377	NAFLD, steatosis ~50%	Male	11 years	C	32.5	None
D01	H0591	Feathering degeneration	Female	11 years	C	19.8	None
D02	H0072	Near normal (no inflammation)	Male	11 years	C	19.2	None
D03	H0703	Near normal, scattered microvesicular steatosis	Male	13 years	AA	12.1	None
D04	H0326	NAFLD, steatosis ~70% in zone 3	Female	14 years	C	32.2	None
D05	H0781	Focal mild lobular infiltration	Female	15 years	C	24.2	None
D06	H0707	Near normal	Male	17 years	C	25.4	None
E01	H1263	Near normal	Male	18 years	C	25.3	None
E02	H1279	Near normal - Control	Male	21 years	C	24.6	None
E03	H1357	Microvesicular steatosis ~20% - Control	Male	22 years	H	22.0	Occasional
E04	H1238	Normal - Control	Female	22 years	C	21.0	Occasional
E05	H1265	Normal - Control	Male	26 years	C	21.1	Occasional
E06	H1325	Normal - Control	Female	27 years	C	20.3	None

There is no history of diabetes for any of these donors.

A: Asian
 AA: African-American
 C: Caucasian
 H: Hispanic

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

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.