

■ General Information

Applications

- Immunohistochemistry
 - TUNEL for apoptosis
- In situ hybridization (ISH)
 - mRNA
 - miRNA
 - Fluorescent In situ hybridization (FISH)

Storage and stability

- Individual slide is put in an air-tight pack with inert gas.
- If the slides are stored at 4C, they are good for up to one year.

How processed

- Tissues were initially fixed with formalin except for some of the animal tissues
- Then, dehydrated with gradient ethanol; typically 1 hour each progressive steps; 70%, 90%, 95%, 99%, 100% x 3 times.
- Cleared by xylene, three changes for 1 hour each.
- Infiltrated with 60°C paraffin, three changes for 1 hour each
- Sectioned by microtome in 4 µm thickness

Before use

- Dry slides for 1 hour in a oven at 60C.
- Dewax slides in xylene for 4 minutes x 5 times.
- Hydrate slides in 100%, 95% and 75% ethanol for 3 minutes x 2 times each.
- Immerse slides in tap water for 5 minutes.

Slide orientation

- In most of the slides with 59 or 60 cores, the orientation is as below unless indicated otherwise. #60 location is usually filled with carbon for orientation.

Shaded area	1	2	3	4	5	6	7	8	9	10
	11	12	13	14	15	16	17	18	19	20
	21	22	23	24	25	26	27	28	29	30
	30	32	33	34	35	36	37	38	39	40
	41	42	43	44	45	46	47	48	49	50
	51	52	53	54	55	56	57	58	59	60

■ Tissue types*

The "tissue type" column in the data sheet denotes the following categories.

1. normal tissue from a non-cancer patient
2. normal tissue from a cancer patient, but the cancer involves unrelated organ
3. normal tissue adjacent to the cancer
4. benign tumor
5. tumor of borderline malignancy or uncertain malignant potential
6. cancer

NBP2-30221 - Human Liver Tissue MicroArray (Cancer metastasis)

No.	Age	Sex	Organ	Diagnosis	pTNM	Stage	Months*	Tissue type**
1	63	M	Liver	hepatocellular carcinoma, moderately differentiated	T2N0M0	II	.	6
2	64	M	Liver	hepatocellular carcinoma, moderately differentiated	T3N0M0	III	.	6
3	62	M	Liver	hepatocellular carcinoma, well & moderately differentiated	T2N0M0	II	.	6
4	65	M	Liver	small cell neuroendocrine carcinoma	T3N0M0	III	.	6
5	67	F	Liver	hepatocellular carcinoma, moderately differentiated	T2N0M0	II	.	6
6	53	M	Liver	hepatocellular carcinoma, moderately differentiated	T3N0M0	III	.	6
7	71	M	Liver	hepatocellular carcinoma, well differentiated	T1N0M0	I	.	6
8	54	M	Liver	hepatocellular carcinoma, moderately differentiated	T2N0M0	II	.	6
9	69	M	Liver	hepatocellular carcinoma, moderately differentiated	T3N0M0	III	.	6
10	57	M	Liver	hepatocellular carcinoma, moderately differentiated	T1N0M0	I	.	6
11	55	M	Liver	hepatocellular carcinoma, well differentiated	T1N0M0	I	.	6
12	45	M	Liver	hepatocellular carcinoma	T1N0M0	I	.	6
13	55	M	Liver	hepatocellular carcinoma	T1N0M0	I	.	6
14	72	M	Liver	hepatocellular carcinoma, moderately differentiated	T1N0M0	I	.	6
15	66	M	Liver	hepatocellular carcinoma	T1N0M0	I	.	6
16	64	M	Liver	hepatocellular carcinoma	T1N0M0	I	.	6
17	44	M	Liver	hepatocellular carcinoma, moderately differentiated	T2N0M0	II	.	6
18	47	M	Liver	hepatocellular carcinoma, moderately differentiated	T2N0M0	II	.	6
19	41	F	Liver	hepatocellular carcinoma, poorly differentiated	T2N0M0	II	.	6
20	51	M	Liver	hepatocellular carcinoma, poorly differentiated	T2N0M0	II	.	6
21	64	M	Liver	hepatocellular carcinoma, moderately differentiated	T1N0M0	I	.	6
22	54	M	Liver	hepatocellular carcinoma, moderately differentiated	T3bN0M0	IIIB	.	6
23	25	M	Liver	hepatocellular carcinoma	T2N0M0	II	.	6
24	32	M	Liver	hepatocellular carcinoma	T3N0M0	III	.	6
25	58	M	Liver	hepatocellular carcinoma	T1N0M0	I	.	6
26	62	M	Liver	hepatocellular carcinoma	T1N0M0	I	.	6
27	52	M	Liver	hepatocellular carcinoma	T2N0M0	II	.	6
28	52	M	Liver	hepatocellular carcinoma, moderately differentiated	T2N0M0	II	.	6
29	71	M	Liver	hepatocellular carcinoma, moderately differentiated	T4N1M0	IVA	.	6
30	64	M	Liver	hepatocellular carcinoma, moderately differentiated	T3bN0M0	IIIB	.	6
31	52	M	Liver	hepatocellular carcinoma, moderately differentiated	T3bN0M0	IIIB	.	6
32	62	M	Liver	hepatocellular carcinoma	T2N0M0	II	.	6
33	56	M	Liver	hepatocellular carcinoma, well differentiated	T1N0M0	I	.	6
34	77	M	Liver	hepatocellular carcinoma, moderately differentiated	T2N0M0	II	.	6
35	38	M	Liver	hepatocellular carcinoma	T2N0M0	II	.	6
36	75	M	Liver	hepatocellular carcinoma, moderately differentiated	T2N0M1	IVB	.	6
37	43	M	Liver	hepatocellular carcinoma, moderately differentiated	T3aN0M0	IIIA	.	6
38	66	M	Liver	hepatocellular carcinoma, moderately differentiated	T2N0M0	II	.	6
39	60	F	Liver	cholangiocarcinoma	T1N0M0	I	.	6
40	58	M	Liver	cholangiocarcinoma	T2N1M0	IVA	.	6
41	51	M	Colon	metastatic hepatocellular carcinoma	TxNxM1	IVB	.	6
42	53	M	Anterior neck	metastatic hepatocellular carcinoma from #20	TxNxM1	IVB	21	6
43	61	M	Lung	metastatic hepatocellular carcinoma	TxNxM1	IVB	.	6
44	60	M	Adrenal gland	metastatic hepatocellular carcinoma	TxNxM1	IVB	.	6
45	39	M	Lumbar vertebra	metastatic hepatocellular carcinoma from #35	TxNxM1	IVB	22	6
46	73	M	Colon	metastatic hepatocellular carcinoma from #36	T2N0M1	IVB	0	6
47	43	M	Soft tissue	metastatic hepatocellular carcinoma from #37	TXNXM1	IVB	5	6
48	72	M	Lymph node	metastatic hepatocellular carcinoma from #38	T4N1M0	IVA	0	6
49	60	F	Soft tissue	metastatic cholangiocarcinoma from #39	TXNXM1	IVB	16	6
50	57	M	Lymph node	metastatic cholangiocarcinoma from #40	T2N1M0	IVA	0	6
51	63	M	Liver	non-neoplastic liver (adjacent to cancer, match of #1)	.	.	.	3
52	64	M	Liver	non-neoplastic liver (adjacent to cancer, match of #2)	.	.	.	3
53	62	M	Liver	non-neoplastic liver (adjacent to cancer, match of #3)	.	.	.	3
54	65	M	Liver	non-neoplastic liver (adjacent to cancer, match of #4)	.	.	.	3
55	53	M	Liver	non-neoplastic liver (adjacent to cancer, match of #6)	.	.	.	3
56	71	M	Liver	non-neoplastic liver (adjacent to cancer, match of #7)	.	.	.	3
57	69	M	Liver	non-neoplastic liver (adjacent to cancer, match of #9)	.	.	.	3
58	55	M	Liver	non-neoplastic liver (adjacent to cancer, match of #13)	.	.	.	3
59	52	M	Liver	non-neoplastic liver (adjacent to cancer, match of #28)	.	.	.	3
60	.	.	Carbon

Months*: interval between primary and metastatic cancer
 TNM and Stage: AJCC Cancer Staging Manual (7th Edition)