

## ■ 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

## NBP2-30271 - Human Liver Tissue MicroArray (Cancer)

No.	Age	Sex	Organ	Diagnosis	pTNM	Stage	No. of NBP2-30276#	Tissue type*
1	60	F	Liver	hepatocellular carcinoma, moderately differentiated	T2N0M0	II	101	6
2	59	M	Liver	hepatocellular carcinoma, poorly differentiated	T2N0M0	II	102	6
3	45	M	Liver	hepatocellular carcinoma, moderately differentiated	T1N0M0	I	103	6
4	45	M	Skin	metastatic hepatocellular carcinoma	TxNxM1	IVB	.	6
5	54	F	Chest wall	metastatic hepatocellular carcinoma	TxNxM1	IVB	.	6
6	48	F	Bone	metastatic hepatocellular carcinoma	TxNxM1	IVB	.	6
7	66	M	Liver	hepatocellular carcinoma, poorly differentiated	T1N0M0	I	107	6
8	43	M	Bile duct	hepatocellular carcinoma	T4NxM0	IIIC	.	6
9	51	M	Colon	metastatic hepatocellular carcinoma	TxNxM1	IVB	.	6
10	49	M	Bile duct	hepatocellular carcinoma	T4NxM0	IIIC	.	6
11	64	M	Liver	hepatocellular carcinoma, poorly differentiated	T1N0M0	I	.	6
12	59	M	Liver	metastatic adenocarcinoma from colon	TxNxM1	IVB	112	6
13	40	M	Lymph node	metastatic hepatocellular carcinoma	T3aN1M	IVA	.	6
14	56	M	Liver	hepatocellular carcinoma, moderately differentiated	T1N0M0	I	.	6
15	73	M	Liver	hepatocellular carcinoma, moderately differentiated	T1N0M0	I	.	6
16	48	F	Bone and soft tissue	metastatic hepatocellular carcinoma	TxNxM1	IVB	.	6
17	68	M	Liver	cholangiocarcinoma, moderately differentiated	T2N0M0	II	117	6
18	50	M	Femur	metastatic hepatocellular carcinoma	TxNxM1	IVB	.	6
19	45	M	Liver	hepatocellular carcinoma, poorly differentiated	T2N0M0	II	119	6
20	52	M	Liver	metastatic adenocarcinoma from rectum	TxNxM1	IVB	120	6
21	47	M	Liver	metastatic adenocarcinoma from rectum	TxNxM1	IVB	121	6
22	39	F	Liver	hepatocellular carcinoma, moderately differentiated	T3aN0M	IIIA	122	6
23	64	M	Liver	hepatocellular carcinoma, moderately differentiated	T3bN0M	IIIB	.	6
24	60	M	Liver	mixed hepatocellular and cholangiocarcinoma, poorly differentiated	T2N0M0	II	124	6
25	56	M	Liver	hepatocellular carcinoma, well differentiated	T1N0M0	I	125	6
26	55	M	Liver	hepatocellular carcinoma, well differentiated	T2N0M0	II	126	6
27	62	M	Liver	hepatocellular carcinoma, well differentiated	T1N0M0	I	127	6
28	52	M	Liver	hepatocellular carcinoma, moderately differentiated	T2N0M0	II	128	6
29	64	M	Liver	hepatocellular carcinoma, moderately differentiated	T2N0M0	II	129	6
30	66	M	Liver	hepatocellular carcinoma, moderately differentiated	T2N0M0	II	130	6
31	54	F	Liver	hepatocellular carcinoma, moderately differentiated	T2N0M0	II	131	6
32	56	M	Liver	hepatocellular carcinoma, poorly differentiated	T1N0M0	I	132	6
33	57	M	Liver	hepatocellular carcinoma, moderately differentiated	T2N0M0	II	133	6
34	61	M	Liver	hepatocellular carcinoma, moderately differentiated	T2N0M0	II	134	6
35	35	M	Omentum	metastatic hepatocellular carcinoma	TxNxM1	IV	.	6
36	53	M	Liver	hepatocellular carcinoma, moderately differentiated	T2N0M0	II	136	6
37	46	M	Liver	hepatocellular carcinoma, moderately differentiated	T1N0M0	I	.	6
38	65	M	Liver	hepatocellular carcinoma, poorly differentiated	T3aN0M	IIIA	138	6
39	56	F	Liver	hepatocellular carcinoma, poorly differentiated	T4N0M0	IIIC	.	6
40	61	M	Liver	hepatocellular carcinoma, moderately differentiated	T3aN0M	IIIA	140	6
41	52	M	Liver	hepatocellular carcinoma, moderately differentiated	T2N0M0	II	.	6
42	57	M	Liver	hepatocellular carcinoma, moderately differentiated	T3aN0M	IIIA	142	6
43	63	M	Liver	hepatocellular carcinoma, moderately differentiated	T2N0M0	II	143	6
44	54	M	Liver	hepatocellular carcinoma, poorly differentiated	T2N0M0	II	144	6
45	44	M	Liver	hepatocellular carcinoma, poorly differentiated	T2N0M0	II	145	6
46	64	M	Liver	hepatocellular carcinoma, moderately differentiated	T1N0M0	I	.	6
47	54	M	Liver	hepatocellular carcinoma, moderately differentiated	T3bN0M	IIIB	.	6
48	64	M	Liver	hepatocellular carcinoma, moderately differentiated	T3aN0M	IIIA	148	6
49	50	M	Adrenal gland	metastatic hepatocellular carcinoma	TxNxM1	IVB	.	6
50	45	M	Liver	cholangiocarcinoma, moderately differentiated	T2N1M0	IVA	150	6
51	58	F	Liver	cholangiocarcinoma, moderately differentiated	T1N0M0	I	151	6
52	64	M	Liver	cholangiocarcinoma, poorly differentiated	T1N0M0	I	152	6
53	55	M	Liver	hepatocellular carcinoma, well differentiated	T1N0M0	I	.	6
54	49	F	Liver	cholangiocarcinoma, moderately differentiated	T1N1M0	IVA	.	6
55	57	F	Liver	hepatocellular carcinoma, poorly differentiated	T2N0M0	II	155	6
56	67	M	Colon	metastatic hepatocellular carcinoma	TxNxM1	IVB	.	6
57	44	M	Liver	hepatocellular carcinoma, poorly differentiated	T3bN0M	IIIB	.	6
58	62	M	Liver	hepatocellular carcinoma, poorly differentiated	T2N0M0	II	.	6
59	42	F	Liver	hepatocellular carcinoma, moderately differentiated	T3aN0M	IIIA	159	6
60	.	.	Carbon	.	.	.	.	.

TNM and Stage: AJCC Cancer Staging Manual (7th Edition)