

■ 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

NBP2-30189 - Human Multi-Organ Tissue MicroArray (Normal)

No.	Age	Sex	Organ	Associated lesion	Tissue type*
1	43	F	Skin	breast cancer	2
2	90	M	Skin	thigh liposarcoma	2
3	49	F	Breast	no abnormal finding	1
4	17	F	Breast	virginal hypertrophy	1
5	62	M	Spleen	pancreas cancer	2
6	42	F	Spleen	pancreas cancer	2
7	81	F	Lymph node	stomach caner	2
8	79	F	Lymph node	stomach caner	2
9	62	F	Skeletal muscle	thyroid cancer	2
10	41	F	Skeletal muscle	breast cancer	2
11	37	F	Lung	metastatic osteosarcoma	2
12	58	F	Lung	lung cancer	3
13	57	M	Heart	rectum cancer	2
14	37	M	Heart	prostatic valve failure	1
15	68	F	Salivary gland	thyroid cancer	2
16	68	M	Salivary gland	alveolar ridge caner	2
17	55	M	Liver	hepatocellular carcinoma	3
18	42	F	Liver	hepatocellular carcinoma	3
19	64	M	Pancreas	ampulla of Vater cancer	3
20	59	F	Pancreas	pancreas cancer	3
21	34	M	Tonsil	chronic tonsillitis	1
22	30	M	Tonsil	chronic tonsillitis	1
23	48	M	Esophagus	esophagus cancer	3
24	71	M	Esophagus	esophagus cancer	3
25	51	M	Stomach, body	stomach caner	3
26	75	M	Stomach, body	stomach caner	3
27	64	M	Small intestine, jejunum	ampulla of Vater cancer	3
28	57	M	Small intestine, jejunum	lung cancer	2
29	83	M	Colon	rectum cancer	3
30	53	M	Colon	colon cancer	3
31	67	M	Rectum	rectum cancer	3
32	74	F	Rectum	rectum cancer	3
33	48	M	Kidney, cortex	renal cell carcinoma	3
34	76	M	Kidney, cortex	renal cell carcinoma	3
35	48	M	Kidney, medulla	renal cell carcinoma	3
36	76	M	Kidney, medulla	renal cell carcinoma	3
37	50	M	Urinary bladder	bladder cancer	3
38	66	M	Urinary bladder	rectum cancer	2
39	32	M	Prostate	bladder cancer	2
40	66	M	Prostate	bladder cancer	2
41	81	M	Testis	no abnormal finding	1
42	65	M	Testis	no abnormal finding	1
43	47	F	Endometrium, proliferative	cervix cancer	3
44	44	F	Endometrium, secretory	cervix cancer	3
45	43	F	Ovary	endometrium cancer	2
46	32	F	Ovary	endometrium cancer	2
47	41	F	Placenta	mature placenta	1
48	30	F	Placenta	mature placenta	1
49	71	M	Adrenal gland	renal cell carcinoma	2
50	50	F	Adrenal gland	renal cell carcinoma	2
51	37	F	Thyroid	thyroid cancer	3
52	60	F	Thyroid	thyroid cancer	3
53	33	F	Thymus	lymphoid hyperplasia	1
54	20	F	Thymus	lymphoid hyperplasia	1
55	19	M	Brain, white matter	pineal germinoma	2
56	32	M	Brain, white matter	renal failure	1
57	19	M	Brain, gray matter	pineal germinoma	2
58	32	M	Brain, gray matter	renal failure	1
59	19	M	Cerebellum	pineal germinoma	2
60	32	M	Cerebellum	renal failure	1