

■ 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

NBP2-30232 - Human Multi-tissue Tissue MicroArray (Normal)

No.	Age	Sex	Organ	Associated lesion	Tissue type*
1	60	F	Skin	breast cancer	2
2	43	F	Skin	breast cancer	2
3	49	F	Subcutis	fatty abdomen	1
4	53	F	Breast	breast cancer	3
5	43	F	Breast	breast cancer	3
6	63	M	Spleen	stomach cancer	2
7	53	F	Spleen	stomach cancer	2
8	49	M	Lymph node	stomach cancer	2
9	46	M	Lymph node	stomach cancer	2
10	61	M	Skeletal muscle	angiosarcoma	2
11	25	M	Nasal mucosa	chronic sinusitis	1
12	55	M	Lung	metastatic cancer of lung (from stomach)	2
13	42	F	Lung	lung cancer	3
14	69	F	Bronchus	lung cancer	3
15	58	M	Heart	no abnormal finding	1
16	50	M	Salivary gland	oropharyngeal cancer	2
17	56	M	Liver	hepatocellular carcinoma	3
18	53	F	Liver	stomach cancer	2
19	42	F	Liver	hepatocellular carcinoma	3
20	64	M	Gallbladder	rectal cancer	2
21	63	M	Pancreas	stomach cancer	2
22	62	M	Pancreas	pancreas islet cell tumor	3
23	33	F	Tonsil	chronic tonsillitis	1
24	54	M	Esophagus	esophageal cancer	3
25	74	M	Esophagus	esophageal cancer	3
26	42	M	Stomach, body	stomach cancer	3
27	49	F	Stomach, body	stomach cancer	3
28	40	M	Stomach, antrum	stomach cancer	3
29	36	M	Stomach, smooth muscle	stomach cancer	3
30	57	F	Duodenum	ampulla of Vater cancer	3
31	77	M	Small bowel	pseudomyxoma peritonei	2
32	56	M	Small bowel	colonic diverticulosis	1
33	72	F	Appendix	metastatic cancer of ovary (from stomach)	2
34	72	F	Colon	rectal cancer	3
35	71	M	Colon	colon cancer	3
36	49	F	Rectum	rectal cancer	3
37	34	M	Kidney cortex	renal cell carcinoma	3
38	44	F	Kidney cortex	renal cell carcinoma	3
39	34	M	Kidney medulla	renal cell carcinoma	3
40	53	M	Urinary bladder	invasive bladder carcinoma	3
41	47	M	Prostate	bladder cancer	2
42	59	M	Prostate	bladder cancer	2
43	59	M	Seminal vesicle	bladder cancer	2
44	72	M	Testis	prostate cancer	2
45	42	F	Endometrium, proliferative	benign ovarian neoplasm	1
46	41	F	Endometrium, secretory	ovarian cancer	2
47	53	F	Myometrium	adenomyosis	1
48	42	F	Uterine cervix	leiomyoma	1
49	53	F	Salpinx	cervix cancer	2
50	39	F	Ovary	ovary cancer	3
51	30	F	Placenta	mature placenta	1
52	uk	uk	Placenta	mid-trimester placenta	1
53	30	F	Umbilical cord	mature placenta	1
54	59	M	Adrenal gland	renal cell carcinoma	2
55	39	M	Thyroid	thyroid cancer	3
56	20	F	Thymus	lymphoid hyperplasia	1
57	0	F	Brain, white matter	no abnormal finding	1
58	0	F	Brain, gray matter	no abnormal finding	1
59	0	F	Cerebellum	no abnormal finding	1
60	.	.	Carbon	.	.