

## ■ 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

## NBP2-30170 - Human Multi-Organ Tissue MicroArray (Normal)

No.	Age	Sex	Organ	Remarks	Tissue type*
1	M	57	Skin	.	1
2	F	38	Skin	.	1
3	M	57	Subcutis	.	1
4	F	38	Subcutis	.	1
5	M	18	Breast	.	1
6	F	35	Breast	.	1
7	M	59	Spleen	.	1
8	M	28	Spleen	.	1
9	M	46	Lymph node	.	1
10	M	60	Lymph node	.	1
11	M	47	Skeletal muscle	.	1
12	F	29	Skeletal muscle	.	1
13	F	40	Lung	.	1
14	F	49	Lung	.	1
15	M	71	Heart	.	1
16	F	31	Heart	.	1
17	F	18	Aorta	.	1
18	F	31	Aorta	.	1
19	F	28	Salivary gland	.	1
20	F	48	Salivary gland	.	1
21	F	29	Liver	.	1
22	M	51	Liver	.	1
23	F	44	Gallbladder	.	1
24	M	58	Gallbladder	.	1
25	M	54	Pancreas	.	1
26	F	31	Pancreas	.	1
27	M	29	Tonsil	.	1
28	M	66	Tonsil	.	1
29	F	18	Esophagus	.	1
30	M	37	Esophagus	.	1
31	M	67	Stomach	.	1
32	M	47	Stomach	.	1
33	M	47	Small intestine	.	1
34	F	67	Small intestine	.	1
35	F	30	Colon	.	1
36	M	37	Colon	.	1
37	F	65	Kidney, cortex	.	1
38	F	47	Kidney, cortex	.	1
39	F	65	Kidney, medulla	.	1
40	F	47	Kidney, medulla	.	1
41	F	40	Uterus	.	1
42	F	62	Uterus	.	1
43	M	85	Prostate	.	1
44	M	58	Prostate	.	1
45	F	29	Placenta	.	1
46	F	25	Placenta	.	1
47	F	32	Umbilical cord	.	1
48	F	25	Umbilical cord	.	1
49	F	18	Adrenal	.	1
50	F	31	Adrenal	.	1
51	F	39	Thyroid	.	1
52	M	51	Thyroid	.	1
53	F	56	Thymus	.	1
54	M	48	Thymus	.	1
55	M	37	Gray matter, cerebrum	.	1
56	F	61	Gray matter, cerebrum	.	1
57	M	37	White matter, cerebrum	.	1
58	F	61	White matter, cerebrum	.	1
59	F	61	Cerebellum	.	1
60	M	51	Cerebellum	.	1