

## ■ 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

### NBP2-30228 - Rat Multi-Organ Tissue MicroArray (Normal)

No.	Sex	Organ	Abnormal finding
1	F	Skin	.
2	M	Skin	.
3	F	Spleen	.
4	M	Spleen	.
5	F	Lymph node	.
6	F	Skeletal muscle	.
7	M	Skeletal muscle	.
8	F	Trachea	.
9	M	Trachea	.
10	F	Lung	.
11	M	Lung	.
12	M	Lung	.
13	F	Heart	.
14	M	Heart	.
15	M	Heart	.
16	F	Salivary gland (mucinous)	.
17	M	Salivary gland (mucinous)	.
18	F	Salivary gland (mucinous)	.
19	M	Salivary gland (mucinous)	.
20	F	Liver	.
21	M	Liver	.
22	M	Liver	.
23	M	Pancreas	.
24	F	Esophagus	.
25	M	Esophagus	.
26	F	Stomach	.
27	M	Stomach	.
28	F	Small intestine	.
29	M	Small intestine	.
30	F	Small intestine	.
31	M	Small intestine	.
32	M	Cecum	.
33	F	Colon	.
34	F	Colon	.
35	M	Colon	.
36	F	Diaphragm	.
37	M	Diaphragm	.
38	F	Kidney (cortex)	.
39	M	Kidney (cortex)	.
40	F	Kidney (medulla)	.
41	M	Kidney (medulla)	.
42	F	Bladder	.
43	M	Bladder	.
44	M	Prostate	.
45	M	Seminal vesicle	.
46	M	Testis	.
47	M	Testis	.
48	F	Uterus	.
49	F	Ovary	.
50	F	Adrenal gland	.
51	M	Adrenal gland	.
52	F	Thymus	.
53	M	Thymus	.
54	F	Brain	.
55	M	Brain	.
56	M	Brain	.
57	F	Cerebellum	.
58	M	Cerebellum	.
59	M	Eye ball	.
60	.	Carbon	.