

■ 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

NBP2-30227 - Rat Multi-Organ Tissue MicroArray (Normal)

No.	Organ	Remarks
1	Skin, ear lobe	.
2	Spleen	.
3	Skeletal muscle, abdominal wall	.
4	Lung	.
5	Heart	.
6	Salivary gland	.
7	Liver	.
8	Pancreas	.
9	Stomach	.
10	Ileum	.
11	Colon	.
12	Kidney, cortex	.
13	Kidney, medulla	.
14	Prostate	.
15	Seminal vesicle	.
16	Testis	.
17	Epididymis	.
18	Uterus	.
19	Adrenal	.
20	Thymus	.
21	Cerebrum	.
22	Pons	.
23	Cerebellum	.
24	Carbon	.

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

Sprague-Dawley rats, 8 weeks old
 Formalin fixed
 Each specimen is 4.0 mm in diameter