

■ 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

NBP2-30187 - Primate Multi-Organ Tissue MicroArray (Normal)

No.	Sex	Organ	Remark
1	M	Skin, abdominal	
2	M	Spleen	
3	M	Lymph node, cervical	
4	M	Bone, skull	
5	M	Bone, rib	
6	M	Bone, sternum	
7	M	Cartilage, ear lobe	
8	M	Cartilage, rib	
9	M	Cartilage, sternum	
10	M	Skeletal muscle, thigh	
11	M	Skeletal muscle, diaphragm	
12	M	Larynx, mucosa	
13	M	Trachea	
14	M	Lung	
15	M	Pericardium	
16	M	Artrium	
17	M	Right ventricle	
18	M	Left ventricle	
19	M	Aorta	
20	M	Salivary gland, parotid	
21	M	Salivary gland, submandibular	
22	M	Liver	
23	M	Gallbladder	
24	M	Pancreas	
25	M	Esophagus	
26	M	Stomach, antral	
27	M	Stomach, fundic	
28	M	Omentum	
29	M	Duodenum	
30	M	Jejunum	
31	M	Ileum	
32	M	Cecum	
33	M	Ascending colon	
34	M	Transverse colon	
35	M	Descending colon	
36	M	Rectum	
37	M	Anus	
38	M	Kidney, cortex	
39	M	Kidney, medulla	
40	M	Ureter	
41	M	Urinary bladder	
42	M	Prostate	
43	M	Seminal vesicle	
44	M	Testis	
45	M	Epidydimis	
46	M	Vas deferens	
47	M	Adrenal	
48	M	Thyroid	
49	M	Thymus	
50	M	Meninx	
51	M	Frontal lobe	
52	M	Parietal lobe	
53	M	Occipital lobe	
54	M	Temporal lobe	
55	M	Thalamus	
56	M	Brain stem	
57	M	Pons	
58	M	Cerebellum	
59	M	Spinal cord	
60	M	Eyeball	

Rhesus macaque (*Macaca mulatta*)
 Perfusion fixation by 4% formaldehyde