

■ 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

NBP2-30186 - Primate Multi-Organ Tissue MicroArray (Normal)

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

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