

■ 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

NBP2-30234 - Human Multi-tissue Tissue MicroArray (Cancer)

No.	Age	Sex	Organ	Diagnosis	Associated lesion	Tissue type*
1	70	M	Skin	normal	liposarcoma	2
2	37	F	Breast	normal	breast cancer	3
3	53	F	Spleen	normal	stomach cancer	2
4	67	M	Skeletal muscle	normal	larynx cancer	2
5	60	M	Lung	normal	lung cancer	3
6	42	F	Liver	normal	hepatocellular carcinoma	3
7	41	M	Stomach, body	normal	stomach cancer	3
8	53	M	Colon	normal	colon cancer	3
9	28	F	Kidney	normal	renal cell carcinoma	3
10	70	M	Prostate	normal	bladder cancer	2
11	26	F	Placenta	normal	normal delivery	1
12	0	F	Brain	normal	none	1
13	32	F	Breast	infiltrating duct carcinoma	.	6
14	51	F	Lung	squamous cell carcinoma	.	6
15	56	M	Liver	hepatocellular carcinoma	.	6
16	52	M	Esophagus	squamous cell carcinoma	.	6
17	49	M	Stomach	adenocarcinoma	.	6
18	56	M	Small bowel	malignant stromal tumor	.	6
19	63	F	Rectum	adenocarcinoma	.	6
20	77	M	Kidney	renal cell carcinoma	.	6
21	66	M	Urinary bladder	transitional cell carcinoma	.	6
22	67	F	Uterus	endometrial carcinoma	.	6
23	34	F	Ovary	mucinous cystadenocarcinoma	.	6
24	64	F	Lymph node	metastatic malignant melanoma (from skin)	.	6

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